

thermo scientific



Thermo

FreeStyle

User Guide

Software Version 1.7

XCALI-98191 Revision A • January 2020

ThermoFisher
SCIENTIFIC

© 2020 Thermo Fisher Scientific Inc. All rights reserved.

Foundation, FreeStyle, and mzVault are trademarks, and Orbitrap, Thermo Scientific, TSQ Endura, TSQ Quantiva, mzCloud, and Xcalibur are registered trademarks of Thermo Fisher Scientific Inc. in the United States.

The following are registered trademarks in the United States and other countries: Adobe, Flash, and Reader are registered trademarks of Adobe Systems Incorporated. Microsoft, Windows, PowerPoint, Paint, Outlook, and Excel are registered trademarks of Microsoft Corporation. NIST is a registered trademark of the National Institute of Standards and Technology. Chemical Abstracts Service is a registered trademark of the American Chemical Society in the United States.

All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

Thermo Fisher Scientific Inc. provides this document to its customers with a product purchase to use in the product operation. This document is copyright protected and any reproduction of the whole or any part of this document is strictly prohibited, except with the written authorization of Thermo Fisher Scientific Inc.

The contents of this document are subject to change without notice. All technical information in this document is for reference purposes only. System configurations and specifications in this document supersede all previous information received by the purchaser.

This document is not part of any sales contract between Thermo Fisher Scientific Inc. and a purchaser. This document shall in no way govern or modify any Terms and Conditions of Sale, which Terms and Conditions of Sale shall govern all conflicting information between the two documents.

Release history: Revision A, January 2020

Software version: (Thermo) FreeStyle 1.7 and later

For Research Use Only. Not for use in diagnostic procedures.

Contents

	Preface	ix
	Related Documentation	x
	System Requirements	x
	Special Notices	xi
	Contacting Us	xii
Chapter 1	Introduction	1
	Mass Spectra	1
	Base Peak	2
	Neutral Losses	3
	Effect of Ionization Modes	4
	Adduct Formation	5
	Effect of Isotopes	6
	Isotope Patterns in High-Resolution Data	7
	Analysis Modes for the Mass Spectrometer	7
	Full Scan	7
	Selected Ion Monitoring (SIM)	8
	MS/MS	8
	Qualitative Analysis Tools	9
	New Features and Enhancements	9
Chapter 2	Using the FreeStyle Window	11
	Startup Window	12
	File Menu	12
	Communicator Bar	14
	Opening Raw Data Files or Sequence Files	15
	Factory Default Layout	17
	Hierarchy of the Chromatogram and Spectrum Views	19
	Creating, Positioning, Previewing, Refreshing, and Closing Workspaces	20
	Arranging Views	23
	Selecting the Columns to Display in a View or Dialog Box with Tabular Data	25
	Saving and Applying Layout Templates	26
	Restoring the Default Settings	27
	Using the Pointer to Update the Timebase or Rescale a Graph	27

Adding Text, Graphic, and Structure Annotations to a Graphical View	28
Copying, Exporting, and Printing Graphical Images and Tabular Data.	34
Copying an Image of a Graphical View to the Clipboard.	34
Copying Tabular Data to the Clipboard	34
Exporting or Printing the Contents of a View or Workspace	35
Setting FreeStyle as the Default Data Visualization Application	36
Selecting How You Open the Floating Toolboxes	37
Toolbars	38
Workspace Options Toolbar	38
Workspace Processing Toolbar	42
Display Options Toolbar.	43
Zoom Options Toolbar.	43
Text and Graphic Annotation Toolbar	44
Views.	46
Info Bar Pages	47
Rearranging the Info Bar Pages	48
Hiding the Info Bar Pages	48
Chapter 3 Reviewing Chromatographic Data.	49
Adding Chromatogram Views	50
Closing Chromatogram Views.	50
Adding Chromatogram Traces	51
Adding Chromatogram Traces Manually.	52
Adding Chromatogram Traces with the Auto Filter Feature	53
Deleting Chromatogram Traces	54
Defining a Chromatogram Trace from the Chromatogram Ranges View	56
Manually Defining Chromatogram Ranges	56
Defining Chromatogram Ranges by Using a Spreadsheet File	59
Displaying Mass Defect Filtered Chromatogram Traces	60
Applying Mass Defect filter in a Chromatogram	62
Linking a Spectrum View to an MDF Chromatogram Trace.	65
Setting up Instrument Status Traces	66
Displaying an EIC Trace by Using the Spectrum Toolbox	67
Using the Scan Filters Page to Display a Filtered Chromatogram	71
Setting Up the Display Options for a Chromatogram Trace	72
Formatting Chromatogram Plots.	72
Labeling Chromatogram Peaks or Local Maxima.	73
Filling the Chromatogram Peaks	75
Normalizing Chromatogram Traces	76
Changing the Y-Axis Title and Scale of the Chromatogram View	81
Changing the Zoom Level of a Chromatogram	82
Automatically Detecting and Integrating Chromatographic Peaks.	84
Selecting the Manual Noise Region for the Genesis and ICIS Algorithms	87

Manually Adding, Undoing, and Deleting Chromatographic Peaks	88
Manually Adding Chromatographic Peaks	88
Undoing the Last Manually Added Peak	89
Deleting Chromatographic Peaks	90
Importing Components from a Processing Method	90
Exporting the Scans for a Filtered Chromatogram to a New RAW File	91
Working with Sequences	92
Creating a New Sequence	93
Comparing Chromatogram Traces in a Sequence	93
Chromatogram-Specific Toolbars	94
Chromatogram – Display Options Toolbar	95
Chromatogram – Workspace Processing Toolbar	99
Sequence Toolbar	100
Chromatogram-Specific Views	101
Chromatogram View	102
Chromatogram Ranges View	104
Peaks List View	110
Chromatogram Toolbox	114
Chromatogram-Specific Pages in the Info Bar	116
Detector Type Page	117
Trace Type Page	118
Scan Filters Page	119
Peak Detection Page	121
Genesis Peak Detection Page	122
ICIS Peak Detection Page	127
Avalon Peak Detection Page	130
PPD Peak Detection Page	133
Sequence File Page	136
Chapter 4 Reviewing Spectral Data	139
Displaying the Scan for a Time Point in a Chromatogram	140
Adding Spectrum Views to the Workspace	140
Selecting Spectra from the Spectrum Ranges Dialog Box	141
Creating a MultiSpectrum View and Changing Its Spectrum Plots	142
Formatting a MultiSpectrum View	144
Linking and Releasing Spectra to and from a Chromatogram	145
Linkage States for a Spectrum Plot	145
Releasing a Linked Spectrum from the Chromatogram	146
Linking a Spectrum View to a Chromatogram	146
Averaging Spectra	146
Creating an Average Spectrum from the Chromatogram Toolbox	147
Creating an Average Spectrum from the Spectrum Ranges Dialog Box	148

Subtracting Background Spectra	149
Using the Chromatogram Toolbox to Subtract Background Spectra	149
Using the Spectrum Ranges Dialog Box to Subtract Background Spectra	151
Exporting a Selected Spectrum to a New RAW File	151
Using Precursor Markers to Open Data-Dependent Scans	152
Using the Precursor Flag Marker	152
Using the Nearby Precursors Marker	155
Selecting Spectra from an MSn Tree	158
Setting Up the Display Options for a Spectrum View	163
Modifying the Scan Header for the Spectrum View	164
Applying Local or Global Normalization in a MultiSpectrum View	165
Labeling Spectrum Peaks	166
Labeling Mass Spectrum Peaks with Chemical Formulas	167
Changing the Y-Axis Scale of a Spectrum View	170
Changing the Zoom Level in a Spectrum View or MultiSpectrum View	171
Spectra-Specific Toolbars	172
Spectrum – Workspace Processing Toolbar	173
Spectrum – Display Options Toolbar	174
MultiSpectrum – Display Options Toolbar	178
Spectra-Specific Views	180
Spectrum View	180
MultiSpectrum View	183
Spectrum Ranges Dialog Box	183
Spectrum Toolbox	185
MSn Browser Page	186
Chapter 5 Reviewing Map Data	191
Working with the Map View	191
Adding Map Views to the Workspace	192
Linking and Releasing Map to and from a Chromatogram	193
Linking Status of a Map View	193
Releasing a Linked Map View from the Chromatogram	194
Linking a Map View to a Chromatogram	195
Changing the Zoom level of a Map View	195
Map View - Specific Toolbars	196
Map View - Workspace Options Toolbar	196
Map View - Display Options Toolbar	197
Map-Specific Views	199
Map Ranges Dialog Box	199
Setting Map Ranges	200
Defining the Retention Time and Mass Ranges	201
Saving a Map View	201

	Setting the Map Display Options	202
	Setting the Map Format Options.	202
	Setting the Map Axis Options	204
	Setting the Map Color Options.	204
	Setting the Band Width.	206
Chapter 6	Analyzing MS Data using Data Analytics View	207
	Adding Data Analytics View to the Workspace	207
	Plotting the Data as a Histogram.	208
	Plotting the Data in a Trend Chart	209
	Data Analytics View - Specific Toolbars	209
	Data Analytics View - Workspace Options Toolbar.	209
	Data Analytics View - Display Options Toolbar	210
	Saving a Data Analytics Plot	212
	Data Analytics Ranges	212
	Filtering Data Analytics Views by Trace Types	213
	Creating a Data Analytics Plot.	217
Chapter 7	Determining the Elemental Composition of Ions.	219
	Overview of an Elemental Composition Analysis.	219
	Starting an Elemental Composition Analysis	221
	Starting an Elemental Composition Analysis from the Toolbar	221
	Starting an Elemental Composition Analysis from the Toolbox.	222
	Using the Fragments Matching Algorithm for Confirmation.	224
	Reviewing the Elemental Composition Results	225
	Reviewing the Best-Matching Formulas.	225
	Reviewing the MS/MS Coverage Score for Matching Fragments.	226
	Reviewing MS/MS Annotation Results	227
	Modifying the Elements in the Elements in Use Table	227
	Adding Elements to the Elements in Use Table.	227
	Removing Elements from the Elements In Use Table	228
	Using Custom Periodic Table for Analysis.	229
	Elemental Composition Results View	231
	Elemental Composition Page.	234
Chapter 8	Searching Mass Spectrum Libraries.	241
	Performing a Local NIST or mzVault Library Search	241
	Setting Up the Default Library Search Parameters.	241
	Selecting the Query Spectrum	243
	Starting a Library Search	243
	Reviewing the Results of a Local NIST or mzVault Library Search	244
	Reviewing the Results of a NIST Library Search	244
	NIST Search Results View.	245
	Reviewing the Results of an mzVault Search	246
	mzVault Search Results View.	248

	Modifying the Settings for a Local NIST or mzVault Library Search	249
	Modifying a NIST Search from the NIST Search Page	250
	Modifying an mzVault Search from the mzVault Search Page	255
	Searching the Online mzCloud Mass Spectral Database	262
	Exporting a Mass Spectrum to the NIST MS Search Application	263
	Managing Libraries	264
	Spectrum Workspace Processing Toolbar – Library Search Buttons	266
Chapter 9	Simulating Isotope Distributions	269
	Simulating the Isotope Distribution for a Chemical Formula or Peptide	269
	Displaying Centroid and Profile Spectra in the Same Window	275
	Isotope Simulation Page	277
Chapter 10	Working with CV Plots for FAIMS Data	281
	CV Traces Dialog Box	281
	CV Plot View	282
Chapter 11	Peptide Fragment Annotation	287
	Overview of Peptide Fragments	287
	Fragmentation Methods	288
	Annotating Peptide Fragments	288
	Peptide Fragments Info Bar	289
	Peptide Results View	291
	Filtering the Fragmented Ions	293
Chapter 12	Viewing Experiment and Instrument Information	295
	Workspace Options Toolbar – Report Area	295
	Spectrum List – Display Options Toolbar	297
	Scan Header View	297
	Spectrum List View	298
	Status Log View	300
	Instrument Method View	301
	Tune Method View	303
	Sample Information View	304
	File Header View	304
	Editing the File Header Values	305
	Error Log View	307
Chapter 13	Deconvolving and Deisotoping Spectra with the Xtract Algorithm	309
	Understanding the Xtract Algorithm	309
	Applying the Xtract Algorithm to an Isotopically Resolved Spectrum	310
	Xtract Page	311
	Deconvolved Spectrum View	317
	Xtract Results View	317

Appendix A FreeStyle Default Settings	323
Modifying, Saving, and Resetting the Default Configuration Options	323
Default Mass Precision Page	325
Default Peak Detection Page	327
Default Recently Used Files Page	327
Default Raw Data Files Directory Page	328
Default Elemental Composition Page	329
Factory Default Template Page	330
Default Library Search Page	332
Default Workspace Options Page	333
Default Xtract Options Page	336
Appendix B Scan Filters and Scan Headers	341
Scan Filter Parameters	341
Scan Headers and Scan Header Abbreviations	344
Appendix C One- and Three-Letter Abbreviations for Amino Acid Residues	347
Appendix D Common Isotopes	351
Index	349

Preface

This guide describes how to use the Thermo FreeStyle™ application to view and analyze raw mass spectrometry data.

For contact information and information about related documentation and system requirements, see these topics.

Contents

- [Related Documentation](#)
- [System Requirements](#)
- [Special Notices](#)
- [Contacting Us](#)

For a list of new features, see [New Features and Enhancements](#). To learn how to use the FreeStyle application, begin with the demonstration animations in Help.

❖ To suggest changes to documentation or to Help


Complete a brief survey about this document by clicking the button below. Thank you in advance for your help.



❖ To open the FreeStyle application

- From the Microsoft™ Windows™ taskbar, choose **Start > All Programs > Thermo FreeStyle > FreeStyle**.

—or—

- Click the **FreeStyle** icon,  , on the desktop.

Related Documentation

The FreeStyle application includes complete documentation. In addition to this guide, you can also access the application Help.

❖ **To view the product manual**

From the taskbar, choose **Start > All Programs > Thermo FreeStyle > Manual > FreeStyle User Guide**.

❖ **To open the Help system from the FreeStyle window**

Choose **File > Help**.

❖ **To view context-sensitive Help**

If information is available for a specific view, page, dialog box, or toolbar, click **Help** or press the F1 key.

System Requirements

Your system must meet the following system requirements.

System	Requirements
Computer	<ul style="list-style-type: none"> • 3.6 GHz quad core processor with a minimum of 8 GB RAM (16 GB recommended) • CD/R-ROM or DVD drive • 1 TB hard drive • Video card and monitor capable of 1920 × 1080 resolution
Software	<ul style="list-style-type: none"> • Adobe™ Flash™ Player (to view the demonstration animations in Help) • Adobe Reader™ 10.1 or later • Microsoft .NET Framework 4.7.2 • Microsoft Office 2013 (for exported data) • Microsoft Windows 7 SP1 (64-bit), Windows 10 (64-bit), or Windows 10 Enterprise 2016 LTSB • Thermo Xcalibur™ 4.3 and later (for the NIST Library Browser)

Special Notices

Make sure you follow the special notices presented in this guide. Special notices appear in boxes; those concerning safety or possible system damage also have corresponding caution symbols.




This guide uses the following types of special notices.

IMPORTANT Highlights information necessary to prevent damage to software, loss of data, or invalid test results; or might contain information that is critical for optimal performance of the system.

Note Highlights information of general interest.

Tip Highlights helpful information that can make a task easier.

Contacting Us

Contact	Email	Telephone	QR Code
U.S. Technical Support	us.techsupport.analyze@thermofisher.com	(U.S.) 1 (800) 532-4752	
U.S. Customer Service and Sales	us.customer-support.analyze@thermofisher.com	(U.S.) 1 (800) 532-4752	
Global support	<ul style="list-style-type: none"> ❖ To find global contact information or customize your request <ol style="list-style-type: none"> 1. Go to thermofisher.com. 2. Click Contact Us, select the country, and then select the type of support you need. 3. At the prompt, type the product name. 4. Use the phone number or complete the online form. ❖ To find product support, knowledge bases, and resources <p>Go to thermofisher.com/us/en/home/technical-resources.</p> ❖ To find product information <p>Go to thermofisher.com/us/en/home/brands/thermo-scientific.</p> 		

Note To provide feedback for this document, go to surveymonkey.com/s/PQM6P62 or send an email message to Technical Publications (techpubs-lcms@thermofisher.com).

Introduction

Use the FreeStyle application to visualize and qualitatively analyze mass spectrometry data. A qualitative analysis focuses on identifying unknown compounds and confirming the presence of target (expected or known) compounds.

With the FreeStyle application, you can display chromatograms and spectra, detect and integrate chromatographic peaks, search mass spectral libraries, simulate mass spectra, subtract background spectra, apply scan filters, annotate plots with text and graphics, create and save layouts, view the status of various instrument parameters during data acquisition, and create a 2D or 3D representation of an analysis displaying the acquired mass/wavelength scans. You can also export spectral data to the NIST™ application or [mzCloud.org](https://mzcloud.org), calculate the elemental composition of a component from its exact mass, set the parameters for defining the charge state for peptide fragment matching, and perform Xtract deconvolution.

For a general understanding of mass spectrometry data and to get started with the application, see these topics.

Contents

- [Mass Spectra](#)
- [Analysis Modes for the Mass Spectrometer](#)
- [Qualitative Analysis Tools](#)
- [New Features and Enhancements](#)

Mass Spectra

There are many different types of mass spectrometry (MS) detectors, but the basic principles are the same in all cases: the MS ionizes the sample, separates the ions according to their mass¹, and moves the separated ions toward a detector where they are counted. The data system compiles a spectrum showing the mass distribution of the ions produced from the sample—a snapshot of ion intensities plotted against their mass-to-charge (m/z) ratios.

¹ In the majority of cases $z=1$ and the x axis becomes equivalent to mass, m .

Ionization initially produces molecular ions, but complex secondary processes can cause the molecular ions to fragment. Together with molecular ions, these fragment ions make up the mass spectrum. For individual chemical substances, a mass spectrum can be a characteristic molecular fingerprint.

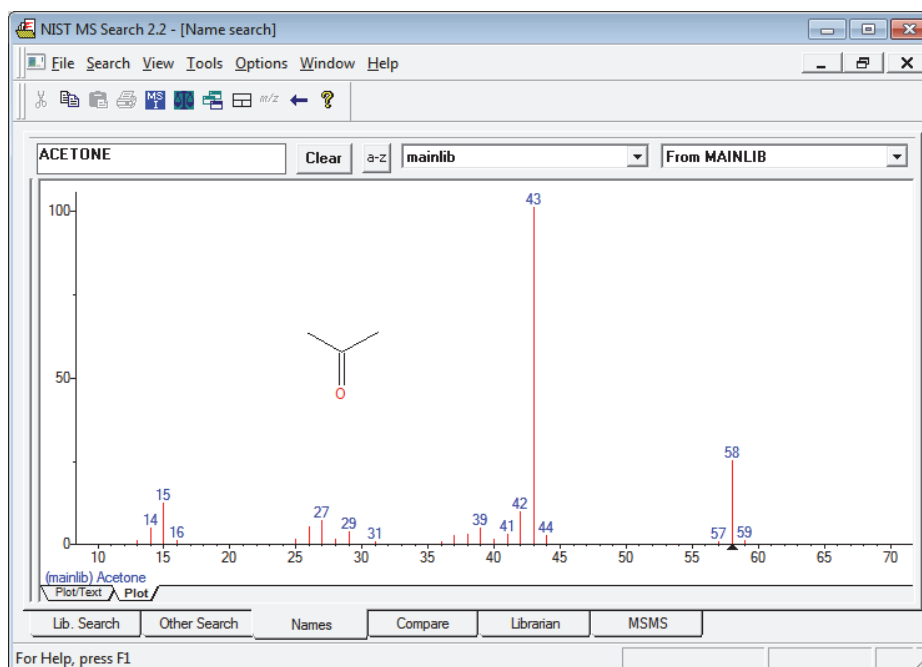
Mass spectra have these common features:

- [Base Peak](#)
- [Neutral Losses](#)
- [Effect of Ionization Modes](#)
- [Adduct Formation](#)
- [Effect of Isotopes](#)
- [Isotope Patterns in High-Resolution Data](#)

Base Peak

To plot the MS detector's response, the most abundant ion, called the base peak, is given an arbitrary abundance or intensity of 100. The application reports other peaks as a percentage of the size of the base peak. After this normalization, the data system can compare spectra directly.

[Figure 1](#) is an example of a NIST library spectrum showing the fragmentation of acetone C_3H_6O (molecular weight = 58 Da). The mass-to-charge labels appear above the spectrum peaks for the most abundant ions. In this example, the molecular ion (58 Da) is not the most abundant ion. The most abundant ion is the acetyl ion CH_3CO (molecular weight = 43 Da).

Figure 1. 70 eV electron ionization (EI) mass spectrum of acetone

Neutral Losses

You can use fragmentation patterns, similar to the pattern displayed in [Figure 1](#) for acetone, to determine the molecular structure of a compound. For example, the neutral loss of 15 Da from the molecular ion of acetone indicates the presence of a methyl group in the original molecule. A subsequent loss of 28 Da corresponds to the loss of CO. [Table 1](#) lists commonly observed neutral losses, measured by the molecular weight of the compound. Assign such losses to help deduce the structure of an unknown compound. A full structural analysis generally relies on the presence of a molecular ion and the measurement of the molecular weight of the compound.

Table 1. Common neutral losses (Sheet 1 of 2)

Loss	Fragment
15	CH ₃
18	H ₂ O
19	F
28	CO
29	C ₂ H ₅ or CHO
35	Cl
46	NO ₂

Table 1. Common neutral losses (Sheet 2 of 2)

Loss	Fragment
59	C ₃ H ₇ O, COOCH ₃ or CH ₂ COOH
77	C ₆ H ₅

In some cases, fragmentation is extensive, leaving little or no trace of a molecular ion. With no molecular ion, determining either the molecular weight or the structure is difficult.

Effect of Ionization Modes

The ionization mode affects the spectrum characteristics of a compound. The ionization modes for LC/MS (liquid chromatograph/mass spectrometer) instruments are different from those used with GC/MS (gas chromatograph/mass spectrometer) instruments:

- [Ionization Modes for LC/MS Instruments](#)
- [Ionization Modes for GC/MS Instruments](#)

Ionization Modes for LC/MS Instruments

LC/MS instruments use a variety of ionization techniques, collectively called atmospheric pressure ionization (API). Detectors of this type can detect positive or negative ions.

API techniques offer soft ionization, usually with little or no fragmentation. An API spectrum typically contains peaks for only the protonated or deprotonated molecule. Compounds with basic sites (such as amines) can form protonated molecules [M+H]⁺. In the positive ion detection mode (polarity +), these ions produce a spectrum peak at the *m/z* value M+1 (where M represents the molecular weight of the neutral compound).

Compounds with acidic sites (sulphonic acids, for example) can form deprotonated molecules [M-H]⁻. In the negative ion detection mode (polarity -), these ions produce a spectrum peak at the *m/z* value M-1.

Ionization Modes for GC/MS Instruments

GC/MS instruments offer two techniques: electron ionization (EI) and chemical ionization (CI).

EI is commonly used because it is simple and reproducible. The fragmentation pattern is effectively determined by the energy of the impacting electrons alone (electron energy, measured in eV). Different types of mass spectrometers that use EI can produce virtually identical spectra as long as they have same electron energy.

This reproducibility has led to an extensive library compilation for 70 eV EI spectra. With the FreeStyle application, you can access the NIST/EPA/NIH Mass Spectral Library with over 108 000 reference EI spectra. You can use library data to select confirmatory ions for your target compounds.

Note You can purchase the NIST Mass Spectral Search application from the National Institute of Science and Technology. Thermo Fisher Scientific provides local versions of the NIST application and its libraries with the Xcalibur data system.

Chemical ionization (CI) offers a softer method of forming ions. In CI, a controlled flow of a reagent gas, commonly ammonia, methane, or isobutane, is introduced into the area where ionization occurs (the ion source). Energetic electrons that pass through the source ionize the reagent gas, as in EI. These ions can then collide with neutral molecules, causing hydrogen transfer. This process is repeated when the reagent gas ions collide with analyte molecules.

CI usually produces protonated molecules, generally at a mass one unit greater than the molecular mass of the compound. Significantly less fragmentation occurs than in comparable EI spectra. Depending on your choice of reagent gas, adduct ions can form. For example, when you use ammonia as the reagent gas, $M+NH_4$ is a typical adduct ion.

Under certain conditions, CI produces negative molecular ions formed by electron capture. The sensitivity of negative ion CI for certain classes of compounds (those containing double bonds, sulfur, phosphorus, chlorine, or bromine) can be orders of magnitude greater than positive CI or EI modes for those compounds.

For more information about the ionization modes available on your instrument, read the hardware manual and the instrument manual on how to get started.

Adduct Formation

If ionization takes place in the presence of contaminants or additives, such as ammonium or sodium ions, some compounds are susceptible to adduct formation. These spectra show other ions in addition to, or instead of, the molecular ion ([Figure 2](#)).

Note The FreeStyle application can automatically add elemental composition and m/z annotations to the mass spectrum peaks (see [Labeling Spectrum Peaks](#)). To add custom annotations, such as those shown in [Figure 2](#), you can use the application's text annotation tools.

Figure 2. Mass spectrum showing sodium and acetonitrile adducts

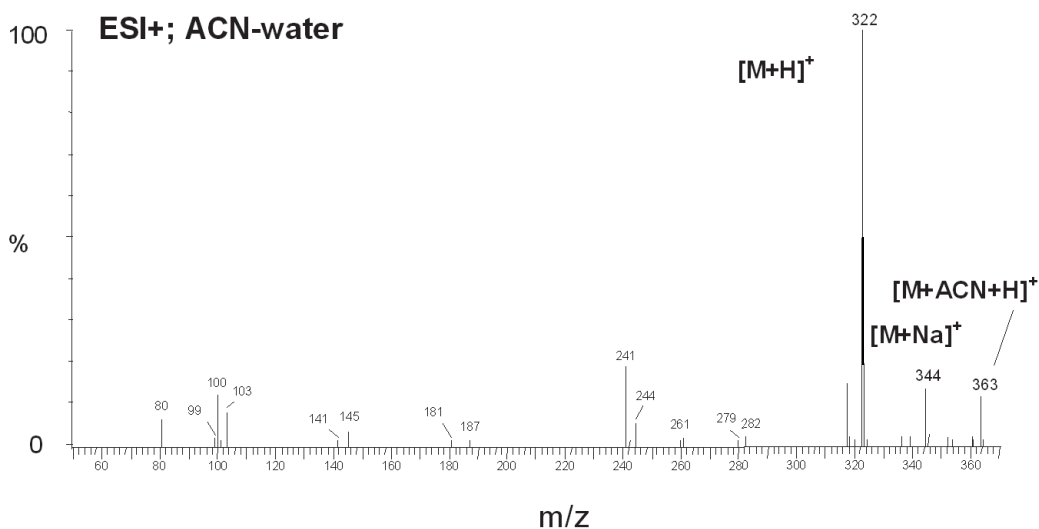


Table 2 lists common adducts for the positive and negative ESI modes.

Table 2. Common adduct ions

Cationized adducts (positive mode)		Anionized adducts (negative mode)	
$[M+\text{NH}_4]^+$	M+18	$[M+\text{OAc}]^-$	M+18
$[M+\text{Na}]^+$	M+23	$[M+\text{Na}]^-$	M+21
$[M+\text{CH}_3\text{OH}+\text{H}]^+$	M+33	$[M+\text{Cl}]^-$	M+35
$[M+\text{K}]^+$	M+39	$[M+\text{K}]^-$	M+37
$[M+\text{CH}_3\text{CN}+\text{H}]^+$	M+42	$[M+\text{HCOO}]^-$	M+59

Take care when determining molecular weights to account for possible adduct ions.

Effect of Isotopes

In some cases, the effect of less abundant isotopes might cause you to use an average molecular weight rather than one based on the most abundant isotopes. When the molecular structure of the target compound contains large numbers of certain elements, the less abundant isotopes become significant. This situation might result in a shift in the mass peaks from their expected m/z values.

For example, the most abundant isotope of chlorine is Cl^{35} . However, Cl^{37} occurs with a natural abundance of 24.47 percent. If a compound contains four chlorine atoms, its molecular ion is two mass units greater than that expected from a calculation based solely on Cl^{35} . Using chlorine's average atomic weight (35.453), you can correctly identify the molecular ion. Also, you observe a distribution of molecular ions across eight mass units from molecules containing between zero and four Cl^{37} atoms.

Isotope Patterns in High-Resolution Data

The mass spectra acquired with a high-resolution, accurate-mass (HRAM) mass spectrometer include isotope clusters for analytes with elements that have more than one stable isotope. The mass difference between the isotopic peaks is proportional to the mass difference of the isotopes, and the relative intensity of the isotopic peaks is proportional to the natural abundance of the isotopes.

To confirm the identity of an unknown analyte, compare the theoretical isotope pattern for its proposed chemical formula to the experimental mass spectrum. Use the [Isotope Simulation Page](#) in the Info Bar to predict the isotope pattern for any chemical formula or peptide sequence.

Analysis Modes for the Mass Spectrometer

A Thermo Scientific mass spectrometer has these analysis modes:

- [Full Scan](#)
- [Selected Ion Monitoring \(SIM\)](#)
- [MS/MS](#)

Full Scan

In full-scan operation, the MS detector scans repetitively over a wide mass range throughout the analysis and sends the data to the data system computer.

With the FreeStyle application, you can display the chromatograms (measured intensity versus analysis time) for full-scan MS data in these ways (plot types):

- As a total ion current (TIC) chromatogram. A TIC chromatogram represents the summed intensities of all the ions in the scanned mass range (mass spectrum) plotted against the chromatographic retention time. Each peak in the TIC represents one or more eluting compounds, which can be identified from the mass spectra recorded across the peak.
- As a mass chromatogram for a range of masses within the scan range. Mass chromatograms show the ion intensities of selected mass-to-charge ratios (m/z). The application extracts these mass spectra from each stored scan and plots them against the analysis time. Use this technique to increase selectivity by displaying an m/z value that is characteristic of the compound of interest but not present in other sample components.
- As a base peak chromatogram. Base peak chromatograms show the ion intensities of the most intense ions for each time point in the chromatogram.

Note The FreeStyle application uses the accurate mass and isotope pattern information in the full-scan MS1 data to calculate the elemental composition of unknown compounds. It then uses the accurate mass data for the fragment ions in the data-dependent MS2 scans to confirm the best matching formulas.

Selected Ion Monitoring (SIM)

In the selected ion monitoring (SIM) mode, the MS detector monitors a limited number of m/z values that are characteristic of a targeted compound or compounds. During an analytical run, the mass analyzer repeatedly switches between the selected m/z values and monitors each m/z value for a programmed dwell time before averaging the measured ion intensities and moving on to the next value.

SIM generates mass chromatograms only of the monitored m/z values, not complete mass spectra as in the full-scan mode. Without a complete mass spectrum, you cannot perform a library search to identify an unknown.

SIM is ideally suited to trace analysis and offers reduced file sizes compared to full-scan operation because SIM records only the information of interest.

MS/MS

Depending on your instrument, you might also be able to do additional stages of mass analysis called MS/MS.

In an MS/MS experiment, you select specific ions for further fragmentation while discarding all other masses. The selected ion is called a precursor (parent) ion and its fragment ions are called product ions. An ion trap mass spectrometer (a mass spectrometer with an ion trap mass analyzer) can perform additional stages of MS (called MSⁿ), up to MS¹⁰.

The MS detector can monitor the product ions in either the full-scan mode or the SIM mode. When you set up the MS detector to monitor a specific product ion of a specific precursor (parent) ion, the scan type is called selective reaction monitoring (SRM).

You can create your own libraries of full-scan MS/MS data to use for matching.

You can display the chromatograms for full-scan MS/MS data in these plot types: TIC, mass range, base peak, neutral fragment, or mass defect filter. With the neutral fragment plot type, you must specify the neutral fragment.

Note The query spectrum for a library search against your local mzVault™ database file or the online mzCloud™ mass spectral library must be an MS2 spectrum from a data-dependent full-scan MS/MS experiment.

Qualitative Analysis Tools

The FreeStyle application includes the following qualitative analysis tools:

- Automated library searches of local mzVault and NIST mass spectral databases and the online mzCloud database. See [Chapter 8, “Searching Mass Spectrum Libraries.”](#)
- Automated peak detection algorithms for the evaluation of chromatographic data (see [Automatically Detecting and Integrating Chromatographic Peaks](#)).
- Elemental composition tool that determines the chemical formulas of ions by using their exact mass and isotope pattern. Includes confirmation of the best match formula by comparing the m/z values of the ions in the data-dependent fragmentation scans of the precursor ion to the possible set of fragment ions based on their m/z values. See [Chapter 7, “Determining the Elemental Composition of Ions.”](#)
- Isotope simulation tool that displays a simulated mass spectrum from the following input: chemical formula, adduct species, and charge distribution. You can compare the simulated spectrum to an experimental spectrum. See [Chapter 9, “Simulating Isotope Distributions.”](#)
- Xtract tool for the deisotoping and deconvolution of mass spectra. See [Chapter 13, “Deconvolving and Deisotoping Spectra with the Xtract Algorithm.”](#)
- Spectrum averaging for noise reduction and mass accuracy (see [Averaging Spectra](#)).
- Background subtraction for the removal of spectral peaks—for example, matrix-related peaks—that are not related to the target components (see [Subtracting Background Spectra](#)).

New Features and Enhancements

The following features are new in FreeStyle 1.7:

- Data Analytics feature that displays options to analyze the MS trending information from mass spectrometers. The analysis is available as a Histogram or Trend plot in a separate window.
- A Mass Defect Filter (MDF) feature is introduced to create a mass defect filtered chromatogram for one or more user-defined mass and mass defect ranges.
- Modifies the Elemental Composition Info page with the addition of a custom periodic table that enables adding custom isotope abundance to the elements for elemental composition calculation and isotope simulations.
- Includes an option to compare the theoretical isotope pattern for the proposed chemical formula by displaying the matched and missed isotopes in the Spectrum view.

- A new customizable File Header view that displays information from the acquisition sequence, the autosampler, and the mass spectrometer.

In addition, these enhancements let you do the following:

- Includes an option to save text and structure annotations in the Spectrum View.
- Includes Ranges as a right-click option in the Map View.
- Modifies the Map Ranges dialog box with the addition of Mass range and Retention Time value fields.
- Includes a new peak identification group section in the Peak Detection Info bar page to identify the nearest, highest, or all peaks.
- Includes an option to import component parameters from the processing method to the current or copy of the selected trace.

The following features are introduced in FreeStyle 1.6:

- Map View feature that displays a map of the currently selected raw file shown in the active Chromatogram plot. The Map view is created in a separate window with retention time (RT) on the x axis and m/z on the z axis.
- Enables the Peptide Fragments button under the Protein Analysis group in the ribbon menu when the spectrum is active. Clicking the Peptide Fragments button displays a new tab in the Info Bar titled Peptide Fragments that lets you to set the parameters for defining the charge state for peptide fragment matching, mass tolerance for annotating the fragments in the spectra, and so on.
- Modifies the Annotation feature with a new option to annotate a spectral peak with a structure, provided as a .mol file.
- Enables the Refresh button under the Create group in ribbon menu when a raw data file in the acquisition queue is opened in the workspace. Clicking the Refresh button refreshes the data of the raw file.

In addition, these enhancements let you do the following:

- Edits the text label and move to any specific position in the view.
- Includes the following new filter options in mzVault search:
 - Collision Energy
 - Compound Class
 - Curation Type
 - Fragmentation Type
 - Peptide Sequence
 - Precursor Mass Range

- Retention Time
- Retention Time Range
- Scan Filter
- Review the results of an MS2 scan in the MSMS Annotation result window.

Using the FreeStyle Window

The FreeStyle window includes multiple toolbars, views, and Info Bar pages. These topics describe the FreeStyle window, the default Workspace layout, the hierarchy of the interactive views, and some common tasks.

Contents

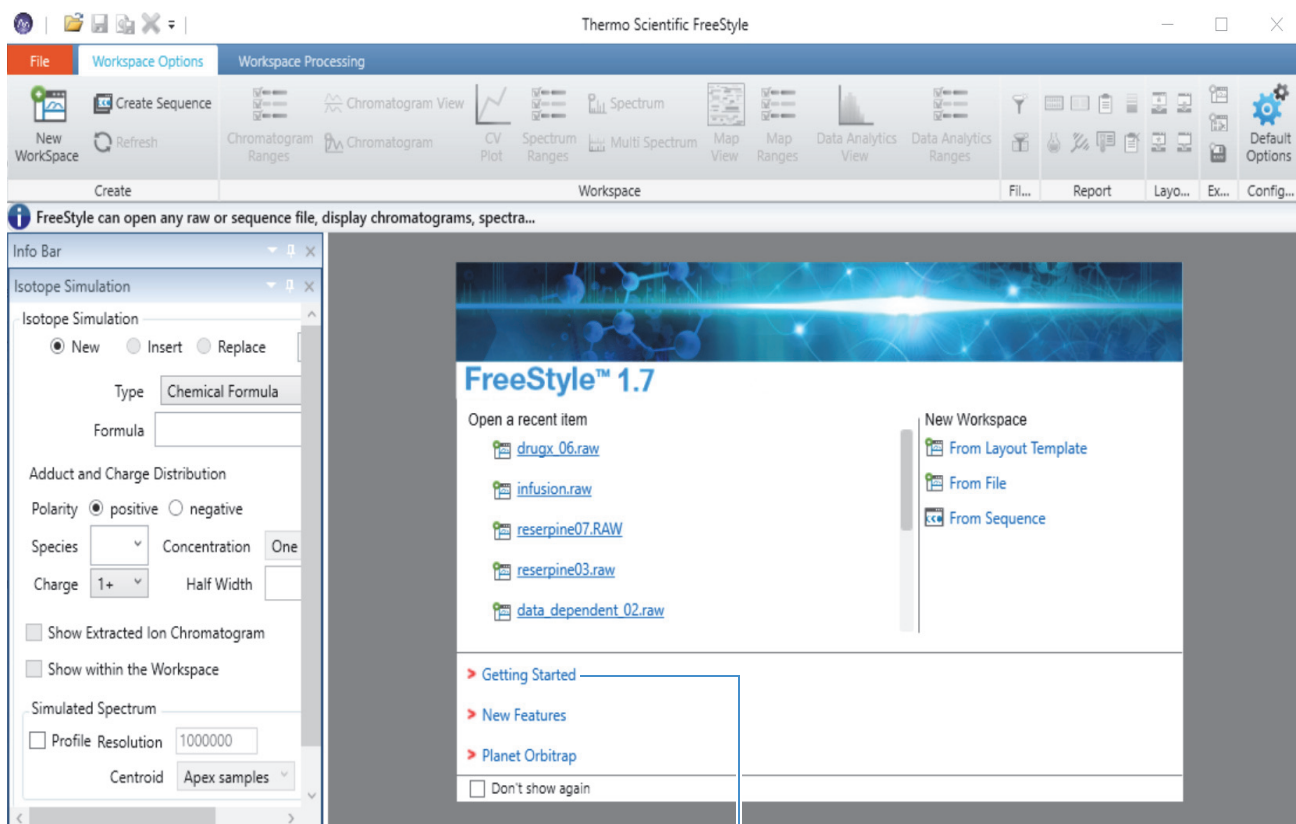
- [Startup Window](#)
- [File Menu](#)
- [Communicator Bar](#)
- [Opening Raw Data Files or Sequence Files](#)
- [Factory Default Layout](#)
- [Hierarchy of the Chromatogram and Spectrum Views](#)
- [Creating, Positioning, Previewing, Refreshing, and Closing Workspaces](#)
- [Arranging Views](#)
- [Selecting the Columns to Display in a View or Dialog Box with Tabular Data](#)
- [Saving and Applying Layout Templates](#)
- [Restoring the Default Settings](#)
- [Using the Pointer to Update the Timebase or Rescale a Graph](#)
- [Adding Text, Graphic, and Structure Annotations to a Graphical View](#)
- [Setting FreeStyle as the Default Data Visualization Application](#)
- [Selecting How You Open the Floating Toolboxes](#)
- [Toolbars](#)
- [Views](#)
- [Info Bar Pages](#)

Startup Window

The FreeStyle window opens to the Isotope Simulation page in the Info Bar to the left and the getting started hyperlinks to the right (Figure 3).

You can use the Isotope Simulation page without opening a raw data file (see [Isotope Simulation Page](#)).

Figure 3. Startup window



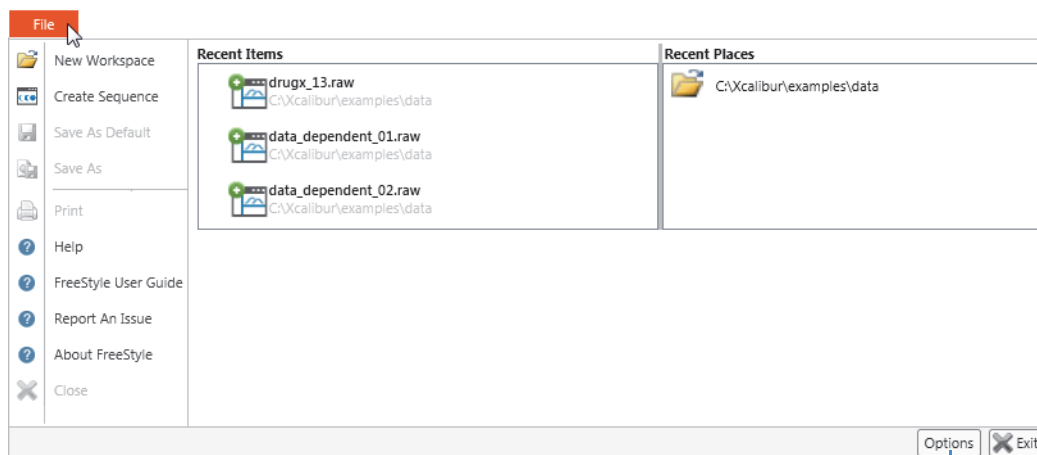
Opens the Help to a list of demonstration animations

File Menu

Use the File menu (see [Figure 4](#)) of the FreeStyle window to open raw data files or sequence files, create a sequence list, save the active Workspace as a template, view the list of recently opened raw data files and folders, and access the default settings.

For more information about Workspaces, see [Creating, Positioning, Previewing, Refreshing, and Closing Workspaces](#). For more information about templates, see [Saving and Applying Layout Templates](#). For more information about default settings, see [Appendix A, “FreeStyle Default Settings.”](#)

Figure 4. File menu



Click to open the Default Options Configuration dialog box

[Table 3](#) describes the File menu commands. For information about using each command, see the related topics.



Table 3. File menu commands (Sheet 1 of 2)

Item	Description
Commands	
New Workspace	Opens a dialog box where you can choose existing raw data files (RAW) or sequence files (SLD) to open and create new Workspaces. See Creating, Positioning, Previewing, Refreshing, and Closing Workspaces .
Create Sequence	Opens a dialog box where you can choose a set of raw data files to open and create a sequence in a new Workspace. See Working with Sequences .
Save as Default	Automatically saves the current layout of the active Workspace to the default template file. See Saving and Applying Layout Templates .
Save As	Opens a dialog box where you can save the current layout to a template file with a different name. In the File Name box, type the new name and click Save . See Saving and Applying Layout Templates .
Print	Prints a report showing chromatogram and spectrum information for the currently selected Workspace. See To save text and structure annotations in the Spectrum view .

Table 3. File menu commands (Sheet 2 of 2)

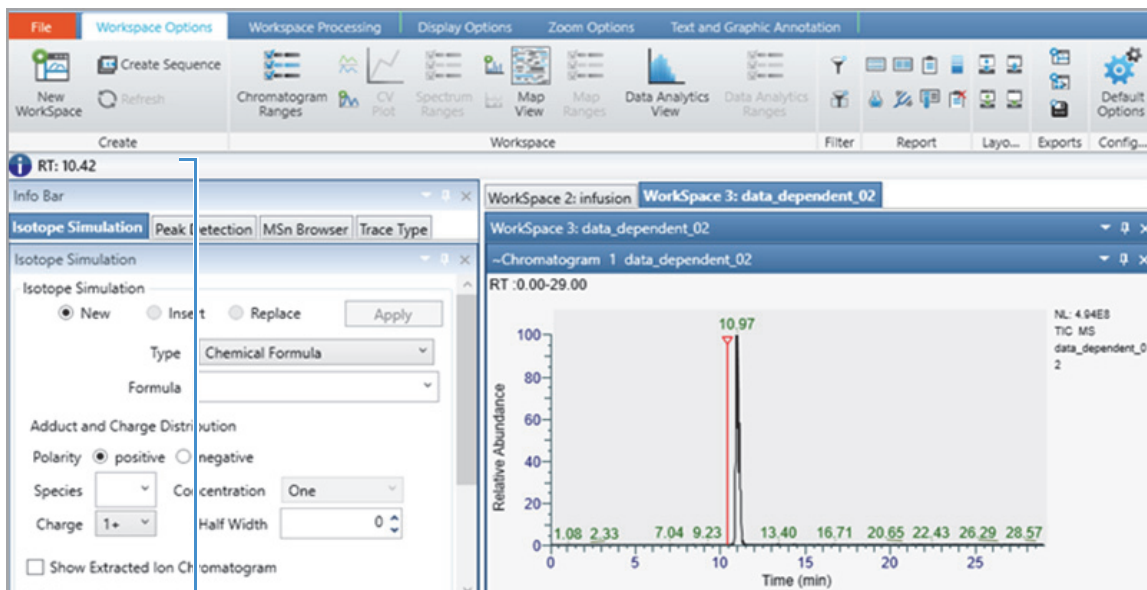
Item	Description
Help	Opens the FreeStyle Help window.
FreeStyle User Guide	Opens the FreeStyle User Guide as a PDF file.
Report an Issue	Opens a mail dialog box in the Microsoft Outlook™ application with a default email address. You can enter a detailed description of the issue, along with the steps to reproduce it, and send this information to Thermo Fisher Scientific.
About FreeStyle	Displays the FreeStyle version information and the release and copyright dates, and shows the version information of other Thermo Scientific applications and instruments installed on your system.
Close	Closes the File menu.
Button	
Options	Displays the current configuration of the default values. See Appendix A, “FreeStyle Default Settings.”
Exit	Closes the FreeStyle application.

Communicator Bar

The communicator bar, which is located immediately below the toolbar, provides information about the selected retention time in the Chromatogram view, general information about the current task, or warning messages. This symbol, , precedes general information, and this symbol, , precedes warnings.

[Figure 5](#) shows the location of the communicator bar.

Figure 5. Communicator bar with informational message



Communicator bar with information about the selected retention time

Opening Raw Data Files or Sequence Files

In the FreeStyle application, you can open two file types—raw data files (RAW) or sequence files (SLD). A raw data file contains the data from a single acquisition run but a sequence file contains a list of associated raw data files from multiple runs. When you open either of these file types, the application creates a new Workspace view.

Follow these procedures:

- To open a raw data file or a sequence file
- To open a raw data file to a custom layout
- To open a file from the startup window

❖ To open a raw data file or a sequence file

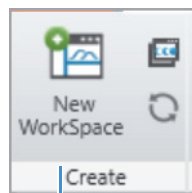
1. Do one of the following:
 - In the menu bar, choose **File > New Workspace**.

—or—

2 Using the FreeStyle Window

Opening Raw Data Files or Sequence Files

- In the Create area of the Workspace Options toolbar, click **New Workspace**.



New Workspace toolbar button

2. In the Open Raw File dialog box, browse to and select a raw data file (RAW) or a sequence file (SLD), and then click **Open**.
 - For a raw data file, a new Workspace appears to the left of the Info Bar and has two stacked views. If the file includes mass spectrometry data, the Chromatogram view at the top displays the TIC trace, and the Spectrum view at the bottom displays the first scan.
 - For a sequence file, the Sequence File page appears in the Info Bar with a list of raw data files, and the Workspace displays data from the first raw data file in the list.

For information about viewing a trace other than the TIC trace, see [Defining a Chromatogram Trace from the Chromatogram Ranges View](#).

❖ To open a raw data file to a custom layout

1. In the startup window, under New Workspace, click **From Layout Template**.

Tip The From Layout Template command is only available in the startup window. To apply a custom layout to an active workspace, from the Workspace Options toolbar, choose **Apply > Named Layout Template**.

2. In the Browse Templates dialog box, select a template (XML) and click **Open**.
3. In the Open Raw File dialog box, select a raw data file and click **Open**.

❖ To open a file from the startup window

Do the following:

- To open a recent file, under Open a Recent Item, click the hyperlink to the file.
- To open a raw data file or a sequence file, under New Workspace, click **From File**.
- To only open a sequence file, under New Workspace, click **From Sequence**.

Factory Default Layout

Figure 6 shows the default FreeStyle window with two raw data files. Each raw data file appears as a separate tabbed Workspace, with only one tabbed Workspace view displayed at a time.

The default layout displays the tabbed Isotope Simulation, Peak Detection, MSn Browser, and Detector Type pages in the Info Bar to the left and the stacked Chromatogram and Spectrum views in the workspace area to the right.

In the title bar for a Chromatogram view or a Spectrum view, the tilde symbol (~) to the left of the view's name indicates that actions in an interactive view or the toolbar affect this view.

Note To open a new Workspace view, see the previous topic [“Opening Raw Data Files or Sequence Files.”](#)

Figure 6. FreeStyle default window



Hierarchy of the Chromatogram and Spectrum Views

A Workspace can contain multiple Chromatogram, Spectrum, and MultiSpectrum views, one each of the report views, and one Chromatogram Ranges view. You can select only one view in the Workspace at a time. Selecting a view turns its title bar a darker shade and adds a tilde to the left of the view's name.

There are three hierarchical states for the Chromatogram, Spectrum, and MultiSpectrum views: selected, active, and inactive (Table 4).

Table 4. Hierarchical states for the Chromatogram, Spectrum, and MultiSpectrum views

State	Indicated by	Description
Selected (and active)	Darker title bar with a tilde (~) to the left of the view's name and a green bar with the selected view's name—Chromatogram, Spectrum, or MultiSpectrum—above the toolbar	<p>Determines the view-specific toolbars and the view-specific tools in the Workspace Processing toolbar.</p> <p>The actions in the view-specific toolbars and the tools in the Workspace Processing toolbar affect the selected view.</p>
Active	Lightly shaded title bar with a tilde (~) to the left of the view's name. The last selected view of each view type is active.	<p>Actions in the selected interactive view affect the active view:</p> <ul style="list-style-type: none"> The actions performed in a linked and selected Spectrum or MultiSpectrum view or the Chromatogram Ranges view affect the active chromatogram plot. The actions performed in the selected chromatogram plot affect the active and linked spectrum. <p>Clicking Chromatogram in the Workspace Options toolbar adds a chromatogram trace to the active or selected Chromatogram view.</p> <p>Clicking Multi Spectrum in the Workspace toolbar changes the active or selected Spectrum view to a Multi Spectrum view.</p> <p>Clicking Chromatogram View or Spectrum in the Workspace Options toolbar adds a copy of the active Chromatogram view or the active Spectrum view, respectively, to the Workspace.</p>
Inactive	Lightly shaded title bar without a tilde	Unaffected by the toolbar actions or actions in other views.

Creating, Positioning, Previewing, Refreshing, and Closing Workspaces

The FreeStyle application creates a new Workspace view whenever you open a raw data file or a sequence of raw data files. The Workspace view displays the processed data (see [Factory Default Layout](#)).

Follow these procedures:

- [To create multiple workspaces](#)
- [To reposition the Workspace view](#)
- [To display a Workspace when multiple Workspaces are open](#)
- [To refresh the Chromatogram and Map views for a Workspace](#)
- [To close a Workspace view](#)

❖ To create multiple workspaces

1. Choose **File > New Workspace**.
2. Do one of the following:
 - Select multiple raw data files and click **Open**.

Tip Use the SHIFT key (for consecutive files) or the CTRL key (for nonconsecutive files).

–or–

- Drag the files into the FreeStyle window.

Note If you drag files into the Chromatogram Ranges view, this action adds new traces to the view instead of adding new workspaces.

A separate Workspace view appears for each raw data file. In the default factory layout, additional workspaces appear as horizontal tabs above the Chromatogram view.

❖ To reposition the Workspace view


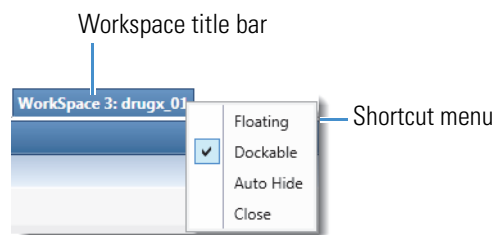
1. Right-click the workspace title bar or click its **Window Position** icon, , to open the shortcut menu.

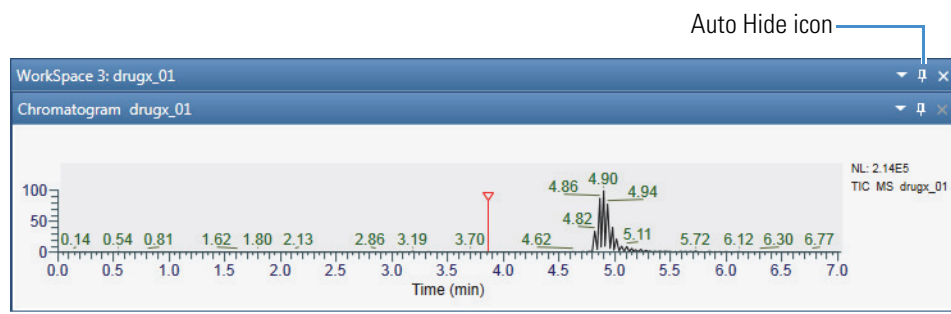
Figure 7. Workspace shortcut menu

2. Do one of the following:

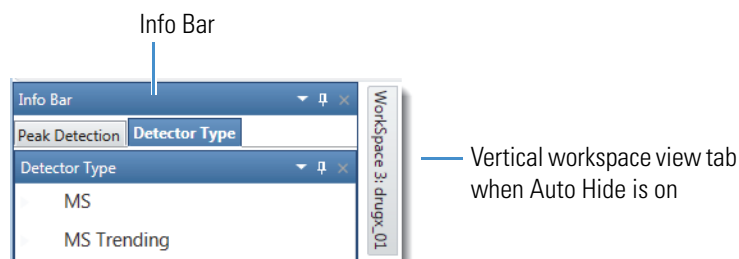
- Choose **Floating** to detach the Workspace view into a floating window.
- Choose **Dockable** to return the Workspace view back to its default position in the FreeStyle window.



—or—

- Choose **Auto Hide** or click the **Auto Hide** (vertical pin) icon, , in the title bar (Figure 8) to temporarily hide the Workspace view.

Figure 8. Workspace vertical Auto Hide icon

The Workspace view collapses as a vertical tab to the right of the Info Bar (Figure 9).

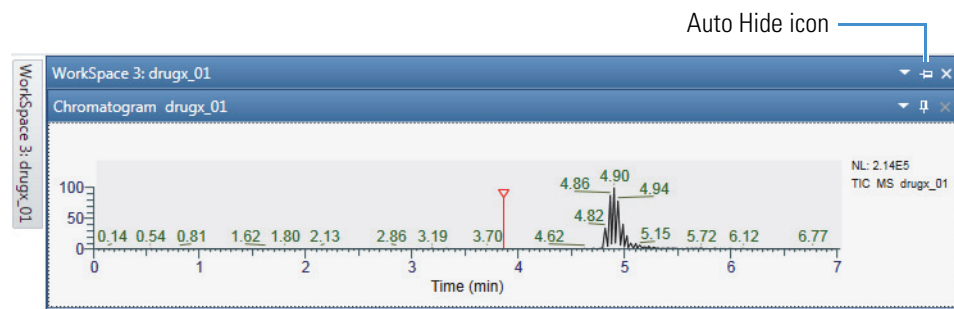
Figure 9. Workspace tab

Pointing to the tab displays the Workspace view again. When Auto Hide is on, the Auto Hide icon is horizontal,  (Figure 10). To turn off the Auto Hide function, choose **Auto Hide** in the shortcut menu again to clear the check mark, or click the Auto Hide icon to change it back to vertical, .

2 Using the FreeStyle Window

Creating, Positioning, Previewing, Refreshing, and Closing Workspaces

Figure 10. Workspace horizontal Auto Hide icon



❖ To display a Workspace when multiple Workspaces are open

- Click the target workspace tab.

–or–

- Click the workspace preview button for the target Workspace.

❖ To refresh the Chromatogram and Map views for a Workspace

1. Open a raw data file in the acquisition queue.
2. In the Workspace Options toolbar, in the Create area, click **Refresh**.
3. (Optional) Press the F5 key on your computer.

The application refreshes the data of the raw data file.

As the acquisition is in progress, the application:

- Updates associated Chromatogram traces.
- Updates Map view, Scan Filters, and MSn Browser that have an associated Time range.
- Automatically applies peak detection to update the retention time of views.
- Updates associated peak list table.

Note When you apply Refresh, the application refreshes all the views of data that is currently being acquired. For files that are not in the acquisition mode in workspace, the Refresh button is inactive.

When acquisition is completed, the Refresh button is inactive and the F5 key does not produce any effect.

❖ To close a Workspace view

- Right-click the title bar and choose **Close**.

–or–

- Click the **Close** icon, .

Arranging Views






You can move or reposition the views to change the layout of the workspace, and then save the modified layout to a template. For information about templates, see [Saving and Applying Layout Templates](#).

❖ To arrange views with the mouse

1. Drag the title bar of the view that you want to move to a second view until the view arranger tool appears.



2. Do one of the tasks in the following table.

Task	Procedure
Move the first view above the second view.	Drag the title bar to the up icon,  .
Move the first view below the second view.	Drag the title bar to the down icon,  .
Move the first view to the left of the second view.	Drag the title bar to the left icon,  .
Move the first view to the right of the second view.	Drag the title bar to the right icon,  .
Make both views tabbed.	Drag the title bar to the tabs icon,  The application displays the first view and creates a tab for the second view.

[Figure 11](#) shows the Spectrum view being dragged to the right of a Chromatogram view, and [Figure 12](#) shows the end result.

Figure 11. Spectrum view being dragged to the right of a Chromatogram view

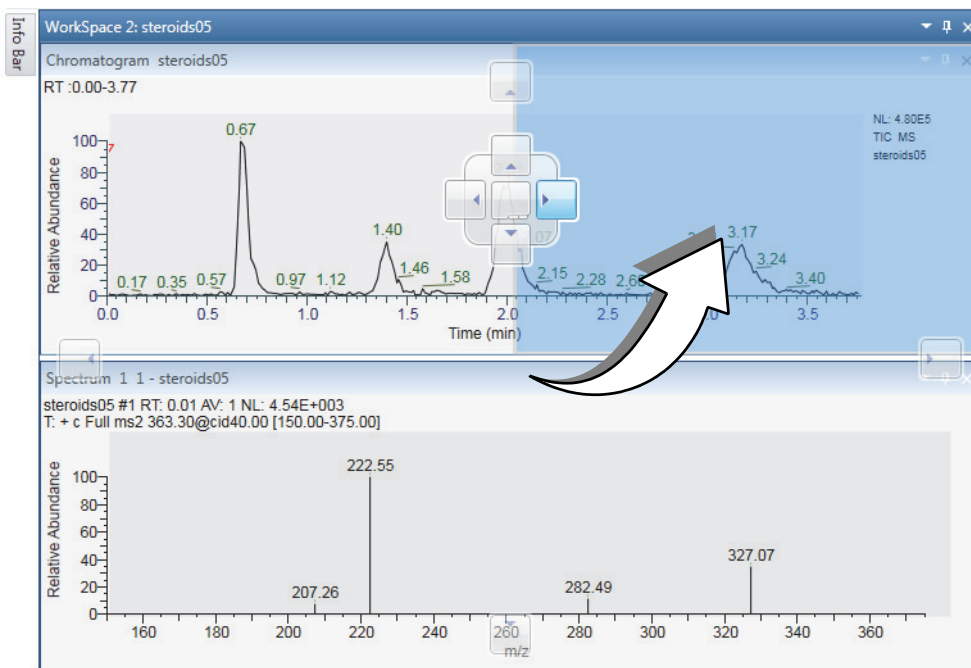
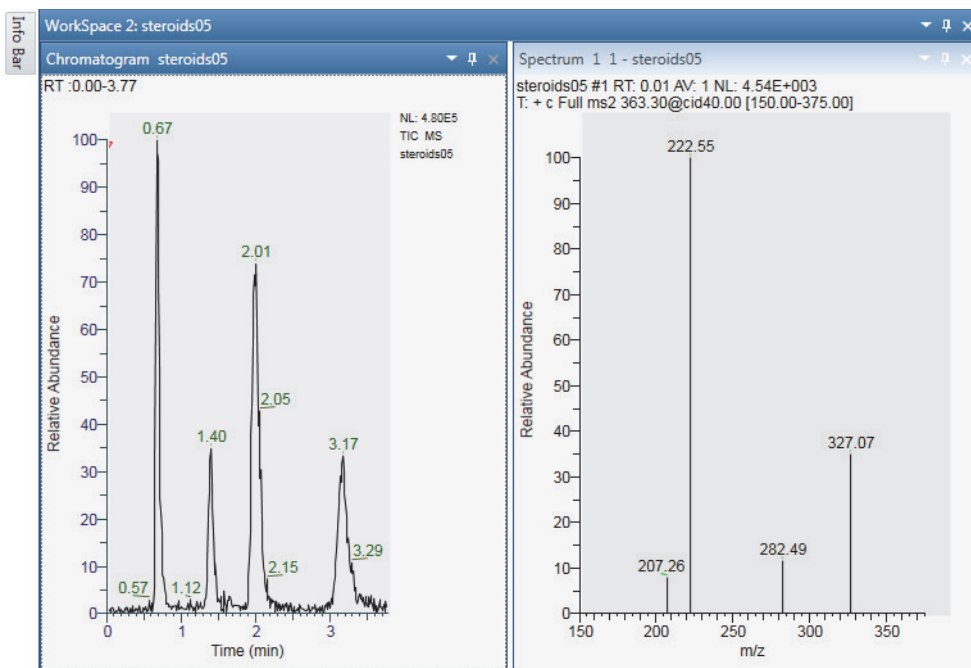


Figure 12. Result of repositioning a Spectrum view to the right of a Chromatogram view



Selecting the Columns to Display in a View or Dialog Box with Tabular Data

To minimize the display of infrequently used information, some of the table columns in a view or dialog box are hidden by default. You can change which columns to display or hide in a specific table by accessing the table's Field Chooser dialog box.

❖ To open a Field Chooser dialog box and change the column selections


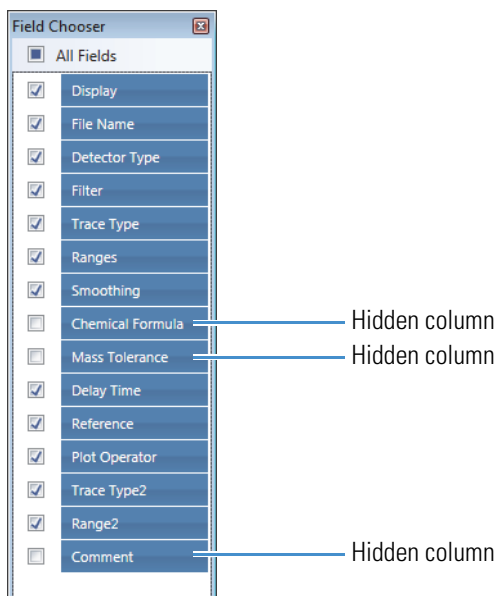
1. Click the **Field Chooser** icon, , to the left of the table heading row.
2. Do the following:
 - To display a column, select its associated check box.
 - To hide a column, clear its associated check box.

Figure 13 shows the Field Chooser dialog box for the Chromatogram Ranges view. In this view, the following columns are hidden by default—Chemical Formula, Mass Tolerance, and Comment.

Figure 13. Field Chooser dialog box for the Chromatogram Ranges view



You can display or hide table columns in these views and dialog boxes:

- [Chromatogram Ranges View](#)
- [Spectrum List View](#)
- [Status Log View](#)
- [Sample Information View](#)
- [Peaks List View](#)

- [Elemental Composition Results View](#)
- [Spectrum Ranges Dialog Box](#)
- Fill Down dialog box (see [Using the Fill Down Feature](#))

Saving and Applying Layout Templates

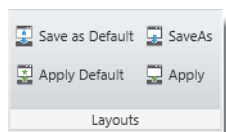
A FreeStyle template contains the settings that define the workspace layout. The template specifies what views are in a Workspace, how the views are arranged, and what labeling the views contain. The template also specifies which pages the Info Bar initially displays. For future use, you can save the template as a new XML file, or save it as the default template.

The FreeStyle application stores all templates in the following folder:

drive:\Users\your name\AppData\Local\Thermo Scientific\FreeStyle\Templates

❖ To save the template as the default template

1. Click the **Workspace Options** toolbar tab (see [Figure 19](#)).
2. Click **Save As Default** in the Layouts area.



Note You can also choose **File > Save As Default**.

The application saves the layout as the default template file, DefaultTemplate.xml.

❖ To apply the default template

1. Click the **Workspace Options** toolbar tab.
2. Click **Apply Default** in the Layouts area.

❖ To name the template and save it as an XML file

1. Click the **Workspace Options** toolbar tab.
2. Click **Save As** in the Layouts area.
3. Type a name for the template and click **Save**.

IMPORTANT Do not navigate to a different folder. The application reads the template files only from the default template folder.

❖ **To apply a saved template**

1. Click the **Workspace Options** toolbar tab.
2. Click **Apply** in the Layouts area to open the DataBrowser Templates dialog box.
3. Select a template and click **OK**.

Restoring the Default Settings

The FreeStyle application includes a dialog box with a set of pages containing initial default settings upon installation. You can modify these settings as needed or restore them to the factory default values.

For information about these settings, see [Appendix A, “FreeStyle Default Settings.”](#)

❖ **To open the Default Options Configuration dialog box**

- In the Workspace Options toolbar, click **Default Options**.

–or–

- a. Choose **File** to open the [File Menu](#).
- b. In the lower right corner, click **Options**.

❖ **To restore the settings of a particular page to the factory default settings**

1. Open the Default Options Configuration dialog box.
2. In the left pane, select the appropriate page of default settings.
3. In the upper right corner, click **Revert to Factory Default Values**.

❖ **To restore all settings to the factory default settings**

1. Open the Default Options Configuration dialog box.
2. In the bottom left corner, click **Revert All to Factory Default Values**.

Using the Pointer to Update the Timebase or Rescale a Graph

Within a Chromatogram View, Spectrum view, or MultiSpectrum view, use the pointer in three ways:

- To select a retention time or scan number, click the linked chromatogram trace at that retention time.
- To select a range, drag a line parallel to any axis.
- To select an area, drag in any diagonal direction.

Pointer actions scale the view according to the dimensions of the dragged line or area (see Table 5).

Table 5. Pointer effects

Pointer action	Effect
Drag parallel to the x axis	Rescales the graph to the selected x -axis range. The y -axis range might rescale depending on the selected normalization display options.
Drag parallel to the y axis	Rescales the graph to the selected y -axis range, with the same x -axis range.
Drag diagonally over the x and y axes	Rescales the graph to the selected x - and y -axes ranges.

In addition, when you select a retention time in a Chromatogram view, the application automatically synchronizes to display data at that selected RT for the linked Spectrum view, or the active spectrum in the linked MultiSpectrum view, the Scan Header view, and the Status Log view in the Workspace.

Adding Text, Graphic, and Structure Annotations to a Graphical View


To add custom text, graphics, and structures to the graphical views, use the Text and Graphic Annotation toolbar.

Note Except for the Structure button, the toolset for Text and Graphic Annotations toolbar are same for the Chromatogram and Spectrum views. To add labels that the application generates from the data, use the Labels area of the Display Options toolbar.





After you select the view and open the Text and Graphic Annotation toolbar, follow these procedures:

- To place a drawing object behind the plot
- To draw a line
- To draw a box or a filled box
- To add a text label
- To edit a text label
- To remove one or more labels
- To save text and structure annotations in the Spectrum view
- To import text and structure annotations
- To annotate a spectral peak with a structure





❖ To place a drawing object behind the plot

Click the **Behind Graph** icon, , before you use the pointer to draw the object.


❖ To draw a line

1. To select the line color, click the **Line** icon, , and then select a color from the graphic.
2. Do one of the following:
 - To draw a horizontal line, click the **Horizontal Line** icon, , and then drag the pointer horizontally across the view.
 - To draw a vertical line, click the **Vertical Line** icon, , and then drag the pointer vertically across the view.
 - To draw a diagonal line, click the **Diagonal Line** icon, , and then drag the pointer diagonally across the view.

❖ To draw a box or a filled box

1. To select the border color, click the **Line** icon, , and then select a color from the graphic.
2. Click the **Box** icon, , or the **Filled Box** icon, .
3. For a Filled Box, click the **Fill** icon, , and then select a color from the graphic.
4. Click the view where you want to place the top left corner of the box, and then drag the pointer across the view to size the box. You can continue to increase or decrease the box dimensions until you release the mouse button.

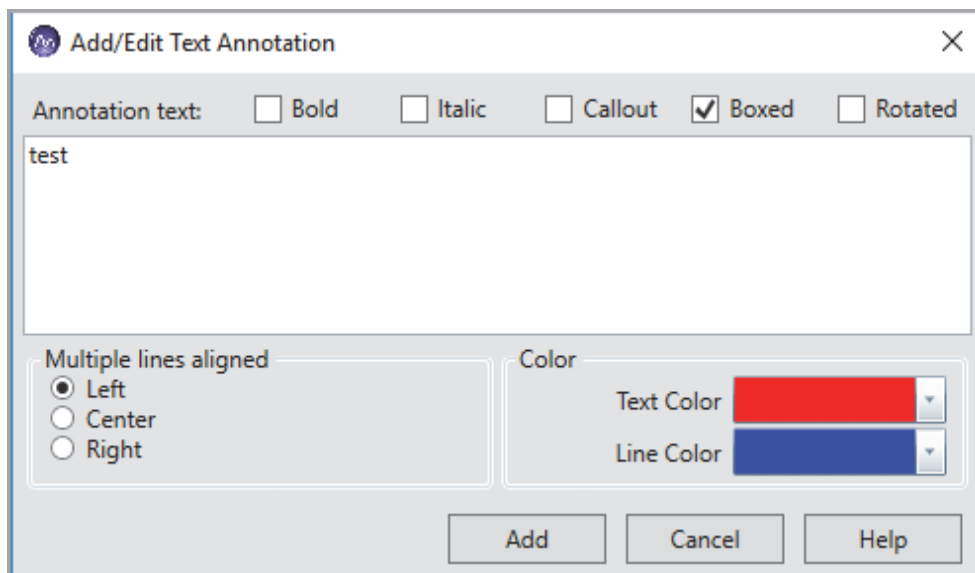
❖ To add a text label

1. Click the **Add Text** icon () to add a text label.
The Add/Edit Text Annotation dialog box opens.

2 Using the FreeStyle Window

Adding Text, Graphic, and Structure Annotations to a Graphical View

Figure 14. Add/Edit Text Annotation



2. Type the text in the text box.
You can enter several lines of text by pressing **ENTER** after each line.
3. To format the text, do the following:
 - To display the text in bold and italic, select **Bold** and **Italic**.
 - To display the text in box or rotate the text, select **Boxed** or **Rotated**.
 - To set up a text label with a callout line, select **Callout**.
 - To align the text, select Left, Center, or Right. By default, if there are multiple lines, the text is left aligned. You can change the alignment to Center or Right.
 - To select the text and line color, under **Color**, click the dropdown button. The default line color is blue and text color is red.
4. To add the text in the Spectrum or Chromatogram view, click the location where you want to add the text.
The text appears in the selected location.
5. To move the text, point to the text, and when the red box appears, drag the text to the new location.

Note When you move the text with a callout line, the callout anchor remains at the original location and only the text moves.

❖ To edit a text label

1. Double-click the text label that you want to edit.
The Add/Edit Text Annotation dialog box appears with the text format options.

2. Select the options to edit, and click **Edit**.

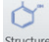
❖ **To remove one or more labels**

Do any of the following:

- To clear selected text labels, click **Selected Text** in the toolbar. Then, drag the pointer across the text that you want to remove.
- To clear a single label, double-click the text label. Then, remove the text from the Add/Edit Text Annotation dialog box, and click **Edit**.
- To clear selected graphic labels, click **Selected Graphics** in the toolbar. Then, drag the pointer across the graphics that you want to remove.
- To clear all the labels, click **All** in the toolbar.

❖ **To annotate a spectral peak with a structure**

1. Select the Spectrum or Multi Spectrum view.
2. To add a structure annotation, do one of the following:

- In the Structure area, click **Structure**  .

The Structure Annotation dialog box appears.

-or-

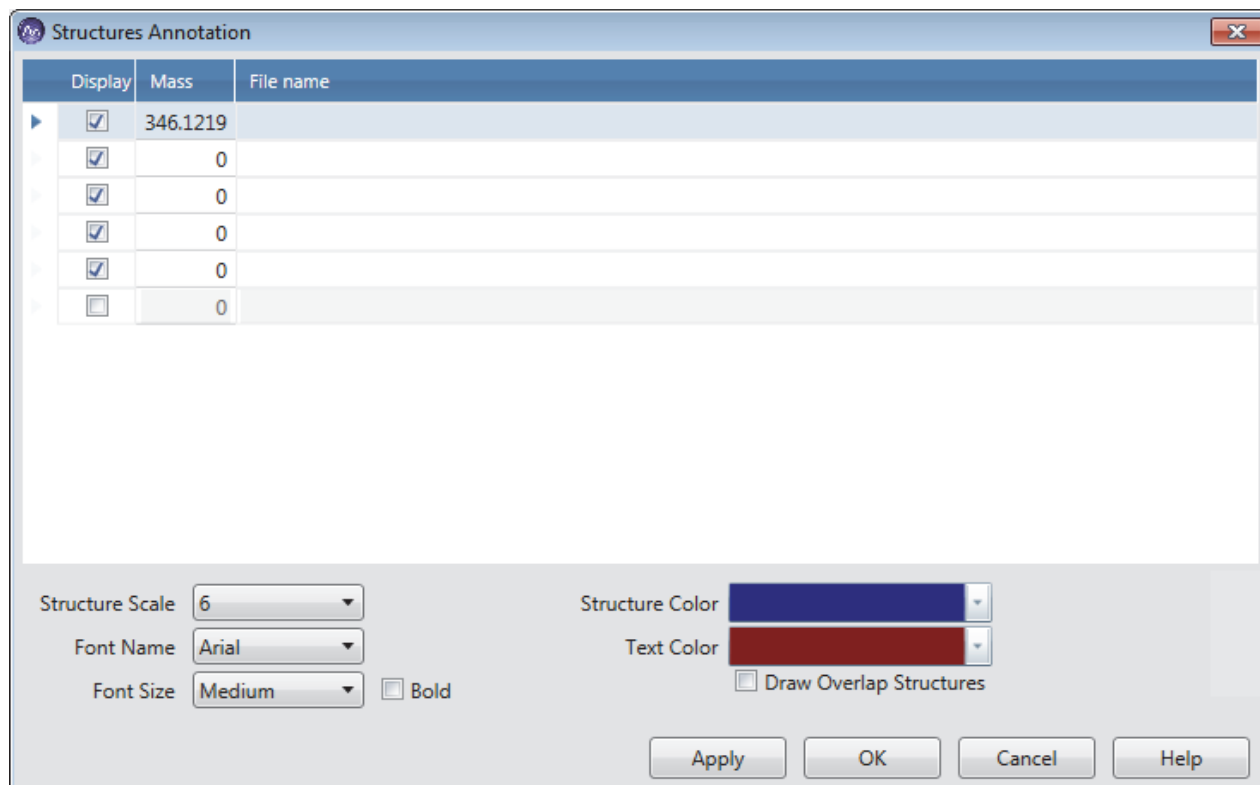
- Right-click the view, select **Show Toolbox**, and click **Structure**.

The Structure Annotation dialog box appears. If you click an m/z label and open the toolbox, the toolbox displays the mass you have selected. The Structure Annotation dialog box also displays the m/z value in the Mass box.


2 Using the FreeStyle Window

Adding Text, Graphic, and Structure Annotations to a Graphical View

Figure 15. Structure Annotation



Note By default, the Structure Annotation dialog box displays five rows. The application automatically adds additional rows to add more annotations. The structure list added is specific for each spectrum plot and you cannot copy the structures to a new plot.

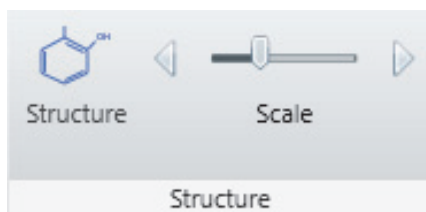
3. To enter or edit the m/z value, click the **Mass** box.
4. In the File name box, click  and browse to the location of the mol file saved in your system,
5. Select the mol file and click **Open** to import the structure.
6. Specify the appearance of the two-dimensional molecular structure as follows:
 - From the Structure Scale dropdown list, select the relative size.
 - From the Font Name dropdown list, select the font for the structure's atoms.
 - From the Font Size dropdown list, select the relative size of the text for the structure's atoms.
 - To bold the text for the structure's atoms, select the **Bold** check box.
 - To select different colors for the structure's bonds and atoms, open respective color palettes and select the colors

- To display the overlapped structures when the structure is bigger than the workspace and does not appear, select the **Draw Overlap Structures** check box. By default, the check box is not selected.
- Click **Apply**, and then click **OK**.

The annotated structure is displayed at the selected mass point in the spectra.

- To change relative size of the spectrum directly from the Text and Graphic Annotation toolbar, click and move the **Scale** slider.

Figure 16. Structure Scale Slider



- To change a structure's position, drag and drop the structure within the view.

The structure changes only the position but remains linked to the original mass value until you edit the mass.

- To clear the structure annotations, right-click the spectrum and choose **Clear Structure Annotations**.

Choosing **Clear Structure Annotations** removes the annotations for the Spectrum view and clears the list from the Structures Annotation dialog box.

Note The application does not import the annotations to a copy of the spectrum within the workspace, but when you copy the spectrum to the Clipboard and paste it into other applications (such as PowerPoint, Paint, and so on), the application also copies the annotations to the new file.

❖ To save text and structure annotations in the Spectrum view

- Select the Spectrum or Multi Spectrum view.
- Click the **Save Annotations button**.

The Save File dialog box opens.

- Select the file to save the annotation, and click **Save**.

The XML file saves the following details:

- Text annotation(s)
- Structure annotation(s)
- Boxed label

2 Using the FreeStyle Window

Copying, Exporting, and Printing Graphical Images and Tabular Data

- Bold and Italic styles
 - Callout link
 - Rotated format
 - Text color
 - Line color
 - Alignment
- ❖ **To import text and structure annotations**
1. In the Save area, click **Apply Annotations**.
 2. Select the annotation XML file you want to import to the workspace.
 3. Click **Open**.

The application imports the text and structure annotation to the spectrum view.

Copying, Exporting, and Printing Graphical Images and Tabular Data

You can copy graphical views and tabular data to the Clipboard, send the contents of a Workspace or view to a printer, or export the contents of a Workspace or view to a CSV or an EMF file.

Follow these topics:

- [Copying an Image of a Graphical View to the Clipboard](#)
- [Copying Tabular Data to the Clipboard](#)
- [Exporting or Printing the Contents of a View or Workspace](#)

Copying an Image of a Graphical View to the Clipboard

You can copy an image of the following views or floating window to paste into other Microsoft Office applications: Chromatogram view, Spectrum view, MultiSpectrum view, and Isotope Simulation window.

- ❖ **To copy an image of a graphical view or floating window to the Clipboard**

Right-click the view or window and choose **Copy To Clipboard**.

Copying Tabular Data to the Clipboard

You can copy the contents of the eight report views (see [Reports](#)) to the Clipboard.

❖ **To copy all the contents of a report view to the Clipboard**

1. Click the view.
2. Press CTRL+C.

Exporting or Printing the Contents of a View or Workspace

To print, copy, or export the contents of a view or workspace, use the Selection As button or Workspace As button, respectively, in the Exports area of the Workspace Options toolbar.

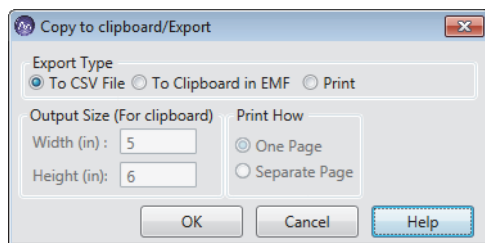
Clicking either button opens the Copy to Clipboard/Export dialog box where you select the export type, the output size for a Clipboard image, and the number of pages to print.

❖ **To export, copy, or print the contents of a view or workspace**

1. Do one of the following:
 - For the contents of a view, click the view, and then in the Workspace Options toolbar, click **Selection As**.
 - For the contents of a workspace, in the Workspace Options toolbar, click **Workspace As**.

The Copy to Clipboard/Export dialog box opens (Figure 17).

Figure 17. Copy to Clipboard/Export dialog box



2. In the Export Type area, select one of these options:
 - To send an enhanced metafile image to the Clipboard, select the **To Clipboard in EMF** option. Then, go to [step 3](#).
 - To send the printing information to the selected printer, select the **Print** option. Then, go to [step 4](#).
 - To export the data in tabular format to a CSV file, select the **To CSV File** option. Then, go to [step 6](#).
3. For the To Clipboard in EMF option, in the Output Size (For Clipboard) area, use the Width (in.) and Height (in.) boxes to specify the size of the image.

2 Using the FreeStyle Window

Setting FreeStyle as the Default Data Visualization Application

- For the Print option, in the Print How area, select one of the following:
 - To print the graphical views in a workspace on separate pages or all the pages in a report view, select the **Separate Page** option.
 - To print the graphical views in a workspace on one page, select the **One Page** option.
- Click **OK**.
- Depending on the export type, do the following:
 - For the To CSV File option, in the Export Data dialog box, select the folder location, name the file, and click **Save**.
 - For the To Clipboard in EMF option, paste the image from the Clipboard to the appropriate document.
 - For the Print option, in the Print dialog box, select the printer, the printing preferences, the page range, and the number of copies. Then, click **Print**.

Tip To create a PDF file, select the Adobe PDF printer.

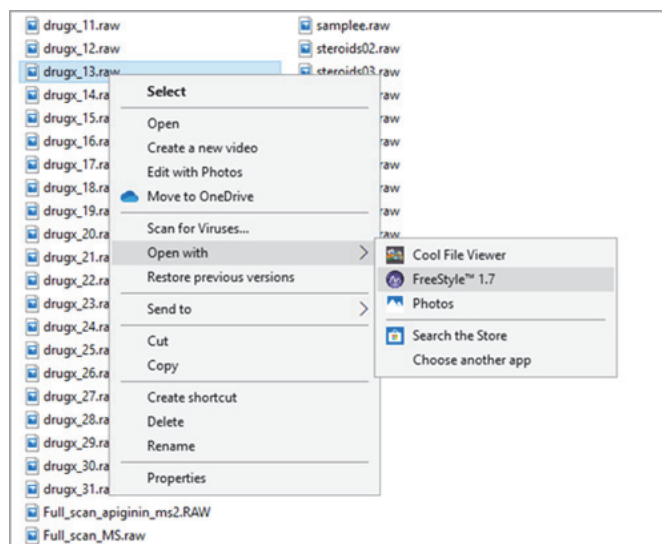
Setting FreeStyle as the Default Data Visualization Application

Currently, Thermo Fisher Scientific provides two data visualization applications: Xcalibur and FreeStyle.

❖ To specify FreeStyle as the default data visualization application

- Open Windows Explorer and browse to a RAW file.
- Right-click the file, and then choose **Open With > FreeStyle Version** (Figure 18).

Figure 18. Choosing FreeStyle as the default data visualization application



Selecting How You Open the Floating Toolboxes

You can open a floating toolbox in the Chromatogram and Spectrum (or MultiSpectrum) views to perform various functions such as averaging scans and adding chromatographic peaks. For information about these toolboxes, see [Chromatogram Toolbox](#) and [Spectrum Toolbox](#).

How you open the floating toolboxes depends on whether the Auto Show Plot Toolboxes After check box is selected on the Workspace Options page. By default, this check box is selected, and a floating toolbox opens when you pause the pointer in the Chromatogram view or point to an m/z value in the Spectrum (or MultiSpectrum) view. If you clear this check box, you can only open the toolbox by choosing Show Toolbox from the view's shortcut menu.



❖ To select the way you open the toolboxes

1. Open the Workspace Options page (see [Default Workspace Options Page](#)).
2. Do one of the following:
 - To make sure that pausing the pointer opens the toolbox, select the **Auto Show Plot Toolboxes After** check box, and then type a pause time for the pointer from **100** to **2000** ms.

With this setting, you can open the toolbox by pausing the pointer in the view or by right-clicking it and choosing Show Toolbox.

- To open the toolbox by using the Show Toolbox command only, clear the **Auto Show Plot Toolboxes After** check box.

With this setting, you can only open the toolbox by right-clicking the view and choosing Show Toolbox.

3. Click **Save**.

Toolbars

Use the FreeStyle toolbar buttons and icons to display views and to perform functions. The FreeStyle window has these toolbars:

- [Workspace Options Toolbar](#)
- [Workspace Processing Toolbar](#)
- [Display Options Toolbar](#)
- [Zoom Options Toolbar](#)
- [Text and Graphic Annotation Toolbar](#)
- [Sequence Toolbar](#) (in the [Reviewing Chromatographic Data](#) chapter)

Tip The FreeStyle window displays only the toolbars and toolbar buttons that are appropriate for the selected view. For example, the Peak Detection buttons in the Workspace Processing toolbar are only available when a Chromatogram view is selected, and the Library Search buttons in this toolbar are only available when a Spectrum or MultiSpectrum view is selected.

The Sequence toolbar appears when you select a file on the Sequence File page.

Workspace Options Toolbar

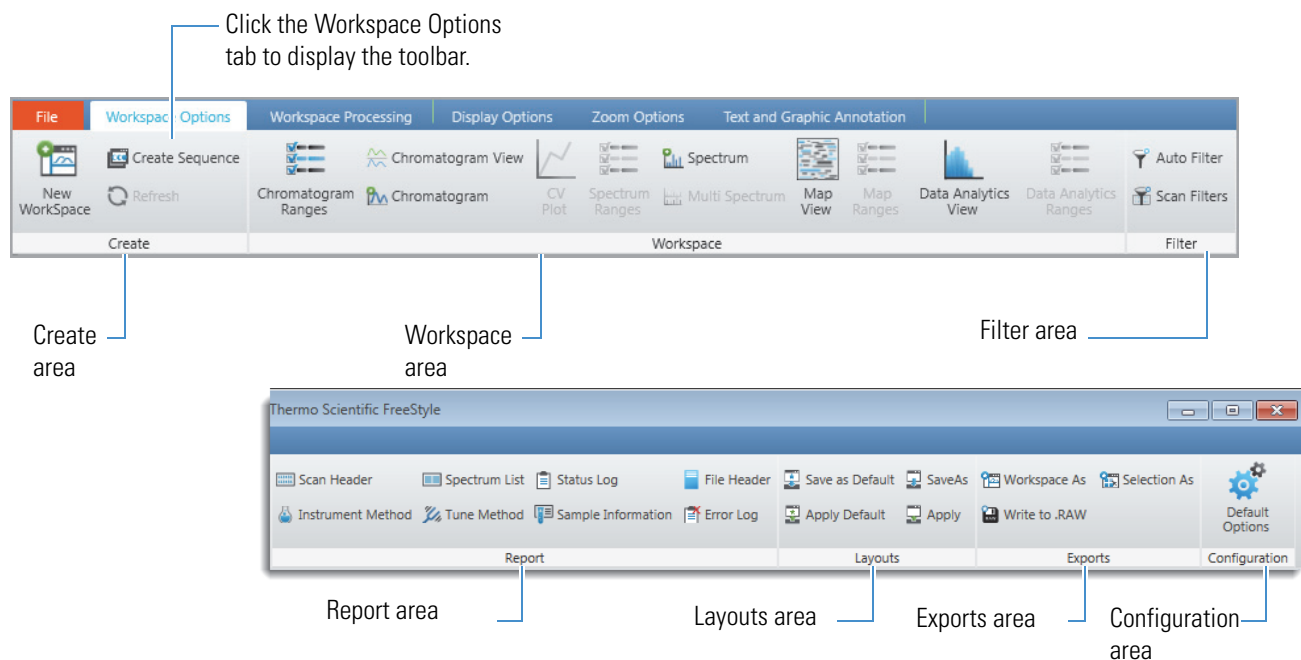
Use the buttons in the Workspace Options toolbar to create, customize, and refresh workspaces.

❖ To display the Workspace Options toolbar

Click the **Workspace Options** toolbar tab.

[Figure 19](#) shows the Workspace Options toolbar buttons that are available when a Chromatogram view is selected.

Figure 19. Workspace Options toolbar for the Chromatogram view (in two parts, left and right)



Depending on the active view, these additional buttons—Spectrum Ranges, Multi Spectrum, CV Plot, and Write to .RAW—become available as follows:

- Selecting a Spectrum view or a MultiSpectrum view enables the Spectrum Ranges and Multi Spectrum buttons.
- Opening a raw data file (RAW) with FAIMS data enables the CV Plot button.
- Selecting a Chromatogram view, a Spectrum view, or a MultiSpectrum view enables the Write to .RAW button.

Table 6 describes the buttons in the Workspace Options toolbar, from left to right.

Table 6. Workspace Options toolbar buttons (Sheet 1 of 4)

Button	Description
Create	
New WorkSpace	Displays the Open Raw File dialog box. You create a workspace by opening a raw data file or a sequence of raw data files. A workspace can display selected information from one or more raw data files. The workspace tab includes the file name of the raw data file when the Workspace displays information from only one raw data file.
Create Sequence	Opens the Create Sequence dialog box. Use the Create Sequence dialog box to build a sequence from raw data files.

Table 6. Workspace Options toolbar buttons (Sheet 2 of 4)

Button	Description
Refresh	Updates a view that is showing a chromatogram or a map from a raw data file that is currently being acquired. The data system expands the display range to show the full range of the data acquired.
Workspace	
Chromatogram Ranges	Opens the Chromatogram Ranges view where you can set up the chromatogram traces to display.
Chromatogram View	Adds a Chromatogram view to the Workspace.
Chromatogram	Adds another chromatogram trace within the selected Chromatogram view. Use the Chromatogram Ranges view or the Auto Filter button to select the chromatograms of interest.
CV Plot	Available for raw data files with FAIMS data. Opens the CV Plot Traces dialog box, where you select the CV plots of interest.
Map View	Adds a Map view of the active chromatogram trace. Displays a map of the currently selected raw file in the map view where you can view the retention time (RT) and m/z values.
Spectrum Ranges	Available when a Spectrum view or a MultiSpectrum view is selected and these views do not include a library spectrum. Displays the Spectrum Ranges dialog box, where you select the spectral trace or traces to display in a Spectrum view or MultiSpectrum view, respectively.
Spectrum	Displays a Spectrum view, where you view the spectrum for the chosen retention time (RT) and scan number.
Multi Spectrum	Displays the MultiSpectrum view, where you can group multiple spectra together in one view.
Filter	
Auto Filter	Populates the Chromatogram view with an unfiltered plot (No Filter) and a plot for every scan filter up to the number specified on the Workspace Options page of the Default Options Configuration dialog box. Also populates the Chromatogram Ranges view with the specified number of scan filters. See Adding Chromatogram Traces with the Auto Filter Feature .
Scan Filters	Opens the Scan Filters page in the Info Bar (see Using the Scan Filters Page to Display a Filtered Chromatogram).

Table 6. Workspace Options toolbar buttons (Sheet 3 of 4)

Button	Description
Report	
Scan Header	Displays the Scan Header view, where you view the scan header of the active raw data file. Click the chromatogram trace to display the scan header for the chosen retention time and scan number. See Scan Headers and Scan Header Abbreviations .
Spectrum List	Displays the Spectrum List view, where you view spectral peak information in a table. Click the chromatogram trace to display the spectrum list for the chosen retention time and scan number.
Status Log	Displays the Status Log view, where you view instrument readback parameters. Click the chromatogram trace to display the status log for a specific retention time and scan number.
Sample Information	Displays the Sample Information view, where you view sample-specific information.
Instrument Method	Displays the Instrument Method view, where you view the instrument method parameters that the instrument used to obtain the raw data file.
Tune Method	Displays the Tune Method view, where you view the tune method parameters that the instrument used to obtain the raw data file.
File Header	Displays the File Header view, where you view information from the acquisition sequence, the autosampler, and the mass spectrometer.
Error Log	Displays the Error Log view, where you view a list of error messages generated during data acquisition.
Layouts	
Save As Default	Saves the current layout as the default template file, DefaultTemplate.xml. See Saving and Applying Layout Templates .
Save As	Saves the current layout as an XML template file. See Saving and Applying Layout Templates .
Apply Default	Applies the default layout template, DefaultTemplate.xml. See Saving and Applying Layout Templates .
Apply	Opens the DataBrowser Templates dialog box, where you select a previously saved layout template. See Saving and Applying Layout Templates .
Exports	
Workspace As	Opens the Copy to Clipboard/Export dialog box, where you specify settings to print all the workspace views, export them to a CSV file, or copy them to the Clipboard (see Exporting or Printing the Contents of a View or Workspace).

Table 6. Workspace Options toolbar buttons (Sheet 4 of 4)

Button	Description
Selection As	Opens the Copy to Clipboard/Export dialog box, where you specify settings to print the selected view, export it to a CSV file, or copy it to the Clipboard (see Exporting or Printing the Contents of a View or Workspace).
Write to .RAW	<p>For a chromatogram trace, saves the MS scans for the displayed time range of the selected chromatogram trace to a new .raw file and appends _FS to the file name (see Exporting the Scans for a Filtered Chromatogram to a New RAW File). The new .raw file only includes the MS scans for the filtered chromatogram.</p> <p>For a spectrum plot, saves the selected spectrum to a new raw file and appends Scan [<i>scan number</i>] to the file name (see Exporting a Selected Spectrum to a New RAW File).</p>
Configuration	
Default Options	Opens the Default Options Configuration dialog box, where you set up the default settings for the application (see FreeStyle Default Settings).

Workspace Processing Toolbar

The available buttons in the Workspace Processing toolbar depend on whether a Chromatogram view or a Spectrum (or MultiSpectrum) view is selected.

Note The following buttons are always available—Elemental Composition, Isotope Simulation, Xtract, Peptide Fragments, and Library Manager.

For information about the Workspace Processing toolbar, see these topics:

- [Chromatogram – Workspace Processing Toolbar](#)
- [Spectrum – Workspace Processing Toolbar](#)
- [Map View - Workspace Options Toolbar](#)
- [Data Analytics View - Workspace Options Toolbar](#)

Display Options Toolbar

For information about specific Display Options toolbars, see these topics:

- [Chromatogram – Display Options Toolbar](#)
- [Spectrum – Display Options Toolbar](#)
- [Spectrum List – Display Options Toolbar](#)
- [Map View - Display Options Toolbar](#)
- [Data Analytics View - Display Options Toolbar](#)

Zoom Options Toolbar

Use the buttons in the Zoom Options toolbar to adjust the display of chromatograms or spectra.

Note The Zoom Options toolbar tab is available when a Chromatogram view, a Spectrum view, a map view, or a MultiSpectrum view is selected.

Figure 20 shows the Zoom Options toolbar, and Table 7 describes the toolbar buttons.

Figure 20. Zoom Options toolbar

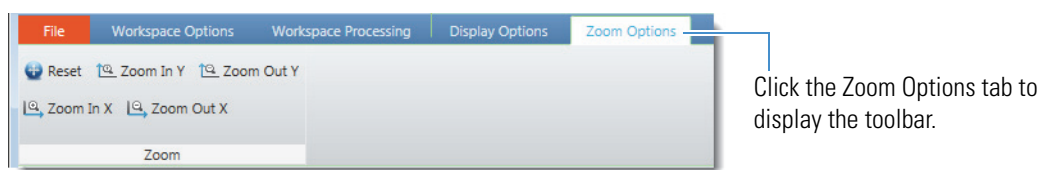


Table 7. Zoom Options toolbar buttons (Sheet 1 of 2)

Button	Description
Reset	Restores the data display to the full range of the x axis and y axis.
Zoom In Y	Zooms in on the y axis by a factor of two from the current baseline to show more detail. For example, you can change the y -axis range from 0–100 to 0–50.

Table 7. Zoom Options toolbar buttons (Sheet 2 of 2)

Button	Description
Zoom Out Y	Zooms out from the <i>y</i> axis by a factor of two from the current baseline to show more data. For example, you can change the <i>y</i> -axis range from 0–25 to 0–50.
Zoom In X	Zooms in on the <i>x</i> axis by a factor of two from the center to show more detail. For example, change the <i>x</i> -axis range from 0–20 to 5–15. Note The first time you click Zoom In X, the application zooms in on the full time range of the active chromatogram plot. The plot's time range depends on the selected scan filter and can be shorter than the data acquisition time for the raw data file. A horizontal scrollbar automatically appears below the <i>x</i> axis of the graphical view when you zoom in.
Zoom Out X	Zooms out from the <i>x</i> axis by a factor of two from the center to show more data. For example, change the <i>x</i> -axis range from 7.5–12.5 to 5–15.

Text and Graphic Annotation Toolbar

Use the buttons and icons in the Text and Graphic Annotation toolbar to annotate chromatograms and spectra with text, lines, and boxes. You can also select color, text alignment, and various text options.

For information about working with the Text and Graphic Annotation toolbar, see [Adding Text, Graphic, and Structure Annotations to a Graphical View](#).

Note The Text and Graphic Annotation toolbar tab is available when a Chromatogram view, a Spectrum view, or a MultiSpectrum view is selected.

Figure 21 shows the Text and Graphic Annotation toolbar, and Table 8 describes the toolbar buttons and icons.

Figure 21. Text and Graphic Annotation toolbar

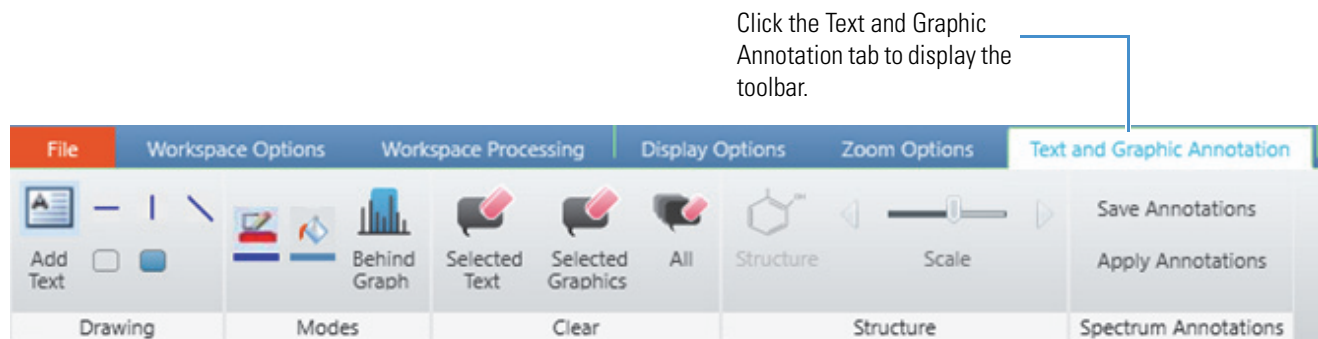


Table 8. Text and Graphic Annotation toolbar buttons and icons (Sheet 1 of 2)






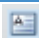



Button/Icon	Description
Drawing	
Horizontal Line 	Draws a horizontal line when you drag the mouse pointer horizontally.
Vertical Line 	Draws a vertical line when you drag the mouse pointer vertically.
Diagonal Line 	Draws a diagonal line when you drag the mouse pointer diagonally.
Box 	Draws an empty box when you drag the mouse pointer diagonally.
Filled Box 	Draws a box filled with the fill color when you drag the mouse pointer diagonally.
Add Text 	Annotates with a text label (see To add a text label).
Modes	
 Line	Opens the color box to select the line color.
 Fill	Opens the color box to select the fill color and the text color.
 Behind Graph	Places text, lines, and boxes behind the chromatogram traces or mass spectra.
Clear	
Selected Text	Deletes text annotations when you drag the pointer over them.
Selected Graphics	Deletes graphic annotations when you drag the pointer over them.
All	Deletes all text and graphic annotations.
Structure	

Table 8. Text and Graphic Annotation toolbar buttons and icons (Sheet 2 of 2)

Button/Icon	Description
Structure	Displays the structure annotations of the selected spectrum.
Spectrum Annotations	
Save Annotations	Saves the text annotation for the selected trace in XML format.
Apply Annotations	Imports the text annotation to the selected trace.

Views

A view is a pane in the Workspace. Use the FreeStyle views to display results and to open lists for entering parameters or selecting options. You can drag views to arrange them within the workspace of the application (see [Arranging Views](#)).

The FreeStyle window displays these individual views (by type in alphabetical order) inside the Workspace view. In addition, [Table 9](#) lists related topics for detailed information and the type of data the view contains for export.

Table 9. Views (Sheet 1 of 2)

View	Topic	Copy as
Primary		
Chromatogram	Chromatogram View	Image
Chromatogram Ranges	Chromatogram Ranges View	Tabular text
Multi Spectrum	MultiSpectrum View	Image
Spectrum	Spectrum View	Image
Map View	Working with the Map View	Image
Data Analytics View	Adding Data Analytics View to the Workspace	Image
Reports		
Error Log	Error Log View	Tabular text
File Header	File Header View	Tabular text
Instrument Method	Instrument Method View	Tabular text
Sample Information	Sample Information View	Tabular text
Scan Header	Scan Header View	Tabular text
Spectrum List	Spectrum List View	Tabular text
Status Log	Status Log View	Tabular text
Tune Method	Status Log View	Tabular text

Table 9. Views (Sheet 2 of 2)

View	Topic	Copy as
Data processing – Elemental composition		
Elemental Composition Results	Elemental Composition Results View	Tabular text
Data processing – Chromatographic peak detection and integration		
Peaks List	Peaks List View	Tabular text
Data processing – Library searches		
Chemical Structure	Chemical Structure view from mzVault or NIST	Image
Compounds list	mzVault Search Results View	Tabular text
Compounds list	NIST Search Results View	Tabular text
Data processing – Xtract deconvolution		
Deconvolved Spectrum	Deconvolved Spectrum View	Image
Xtract Results	Xtract Results View	Tabular text

Info Bar Pages

On the far left of the FreeStyle window, use the pages of the Info Bar to select detector, sequence, or filter options. You can also set the parameters for peak detection, isotope simulation, elemental composition, peptide fragments analysis, searching mass spectrum libraries, or applying the Xtract deconvolution algorithm.

The following pages appear in the Info Bar:

- Chromatogram-range pages: Only one of these two pages is available at any time.
 - [Detector Type Page](#)
 - [Trace Type Page](#)
- [Scan Filters Page](#)
- [Elemental Composition Page](#)
- [Isotope Simulation Page](#)
- [Modifying an mzVault Search from the mzVault Search Page](#)
- [MSn Browser Page](#)
- [Peptide Fragments Info Bar](#)
- [Modifying a NIST Search from the NIST Search Page](#)
- [Peak Detection Page](#)

2 Using the FreeStyle Window

Info Bar Pages

- [Avalon Peak Detection Page](#)
- [Genesis Peak Detection Page](#)
- [ICIS Peak Detection Page](#)
- [PPD Peak Detection Page](#)
- [Sequence File Page](#)
- [Xtract Page](#)

To rearrange or hide the Info Bar pages, follow these topics:

- [Rearranging the Info Bar Pages](#)
- [Hiding the Info Bar Pages](#)

Rearranging the Info Bar Pages

❖ To move an Info Bar page

Drag its tab left or right.

Hiding the Info Bar Pages

❖ To hide an Info Bar page

Right-click its tab and choose **Auto Hide**.

❖ To hide all the Info Bar pages

Right-click the Info Bar title bar and choose **Auto Hide**.

Reviewing Chromatographic Data

To review the chromatographic data, follow these topics.

Contents

- [Adding Chromatogram Views](#)
- [Closing Chromatogram Views](#)
- [Adding Chromatogram Traces](#)
- [Deleting Chromatogram Traces](#)
- [Displaying Mass Defect Filtered Chromatogram Traces](#)
- [Setting up Instrument Status Traces](#)
- [Displaying an EIC Trace by Using the Spectrum Toolbox](#)
- [Using the Scan Filters Page to Display a Filtered Chromatogram](#)
- [Setting Up the Display Options for a Chromatogram Trace](#)
- [Changing the Zoom Level of a Chromatogram](#)
- [Automatically Detecting and Integrating Chromatographic Peaks](#)
- [Selecting the Manual Noise Region for the Genesis and ICIS Algorithms](#)
- [Manually Adding, Undoing, and Deleting Chromatographic Peaks](#)
- [Importing Components from a Processing Method](#)
- [Exporting the Scans for a Filtered Chromatogram to a New RAW File](#)
- [Working with Sequences](#)
- [Chromatogram-Specific Toolbars](#)
- [Chromatogram-Specific Views](#)
- [Chromatogram Toolbox](#)
- [Chromatogram-Specific Pages in the Info Bar](#)

Note For information about adding images and annotations to and copying an image of Chromatogram view, see [Chapter 2, “Using the FreeStyle Window.”](#) .

Adding Chromatogram Views

You can add multiple Chromatogram views to a Workspace.

❖ To add a Chromatogram view

In the Workspace Options toolbar, click **Chromatogram View**.

Result:

- If the Workspace includes one or more Chromatogram views, a copy of the selected or active Chromatogram view (as indicated by the tilde in its title bar) appears.
- If the Workspace does not include a Chromatogram view, a new Chromatogram view appears with the default chromatogram (based on the raw data file) and display settings. The chromatogram trace is linked to the MSn Browser page and any other opened view.

Note If the Workspace does not include a Chromatogram view, you can also add the first Chromatogram view by clicking Auto Filter on the Workspace Options toolbar (see [Adding Chromatogram Traces with the Auto Filter Feature](#)).

Closing Chromatogram Views

A Workspace must include at least one of the following views: a Spectrum view, a Chromatogram view, a Map view, or a report view. When you close all the Chromatogram views, the remaining views and the MSn Browser page retain information based on the last selected chromatogram trace.

❖ To close a Chromatogram view

Click the **Close** icon.

The effect of closing a Chromatogram view depends on whether it is the only Chromatogram view in the Workspace, whether any other views are open, and the linkage state of any spectra ([Table 10](#)).

Table 10. Effect of closing a Chromatogram view

When you close ...	The application does the following ...
The last Chromatogram view in the Workspace	<ul style="list-style-type: none"> Releases any spectrum that is linked to the chromatogram. The text appended to the Spectrum view's title bar changes from CxTy to Released. Closes the Chromatogram Ranges view if it is open. Closes the CV Plot view if it is open. Closes the Workspace if the Chromatogram view was the last view in the Workspace.
One of the Chromatogram views in the Workspace	<ul style="list-style-type: none"> Makes any spectrum that was linked to a chromatogram in the closed view available for linking to another chromatogram. <ul style="list-style-type: none"> The text appended to the Spectrum view's title bar changes from CxTy to 'Select Trace to Link'. Clicking another Chromatogram view automatically links the spectrum to the selected chromatogram. Links an opened Chromatogram Ranges view to the active Chromatogram view. Closes the CV Plot if it is open.

Adding Chromatogram Traces

The factory default Workspace includes a Chromatogram view with a single trace and a Spectrum view that is linked to the trace (see [Opening Raw Data Files or Sequence Files](#)). You can add multiple chromatogram traces to the Chromatogram view and multiple Chromatogram views to the Workspace.

By default, when you add multiple traces to a Chromatogram view, the traces are stacked. As you add more traces, each successive trace appears in a different color until the view contains eight traces. The color order is as follows: (1) black, (2) brick-red, (3) green, (4) blue, (5) light orange, (6) magenta, (7) blue-green, and (8) gray. As you add more traces, this color pattern repeats.

The application supports the following types of chromatogram traces for MS data ([Table 11](#)). To select the TIC, BPC, Mass Range, Base Peak, Neutral Fragment, or Mass Defect Filter trace type, use the Chromatogram Ranges view.

Table 11. Supported chromatogram traces

Trace type	Description
Total ion current (TIC)	A trace of the signal of all ions combined as a function of the retention time or scan number.
Base peak (BPC)	A trace of the most intense ion signal as a function of the retention time or scan number.
Mass range	A trace of the ion signal of all ions in one or more mass ranges as a function of the retention time or scan number.
Extracted ion (EIC or XIC)	A trace of the signal of a single ion as a function of the retention time or scan number.
Neutral fragment	A trace of the ion signal of all ions that produce a specific neutral fragment as a function of the retention time or scan number.
Mass Defect Filter	A trace of the ion signal of all ions in one or more mass defect ranges as a function of the retention time or scan number.

Follow these topics as needed:

- [Adding Chromatogram Traces Manually](#)
- [Adding Chromatogram Traces with the Auto Filter Feature](#)

Adding Chromatogram Traces Manually

To manually add chromatogram traces to a Chromatogram view, follow these procedures as needed:

- To add a chromatogram trace by using the **Chromatogram** button
- To add a chromatogram trace by using the shortcut menu command
- To add chromatogram traces by using the **Chromatogram Ranges** view
- To add a chromatogram trace from a raw data file in a sequence list

Note To automatically populate a Chromatogram view with traces from multiple scan filters, see the next topic “[Adding Chromatogram Traces with the Auto Filter Feature.](#)”

❖ To add a chromatogram trace by using the **Chromatogram** button

1. Click a Chromatogram view or a specific trace in the Chromatogram view to select it.
2. Click the **Workspace Options** toolbar tab.
3. In the Workspace Options toolbar, click **Chromatogram**.

A copy of the selected chromatogram appears at the bottom of the view.

❖ **To add a chromatogram trace by using the shortcut menu command**

Right-click a Chromatogram view and choose **Insert Chromatogram**.

A copy of the selected chromatogram appears at the bottom of the view.

❖ **To add chromatogram traces by using the Chromatogram Ranges view**

1. In the Workspace Options toolbar, click **Chromatogram Ranges**.

The Chromatogram Ranges dialog box linked to the active or selected Chromatogram view opens.

2. In the last row, select the check box in the Display column.

The application populates the row with a copy of the currently selected row in the list, and a duplicate chromatogram trace appears in the Chromatogram view. To fill multiple rows, see [Using the Fill Down Feature](#). To populate the Chromatogram Ranges view with ranges from a spreadsheet file, see [To specify the ranges of the chromatogram traces by using data in a spreadsheet file](#).

❖ **To add a chromatogram trace from a raw data file in a sequence list**

1. Open an existing sequence file (SLD), or create a temporary sequence list by clicking **Create Sequence** in the Workspace Options toolbar and selecting a set of raw data files.
2. To display the Sequence toolbar, click the **Sequence File** tab in the Info Bar.
3. Click **New Trace**.
4. Click a raw data file in the sequence.

The application adds the raw data file to the chromatogram ranges list and displays the new chromatogram trace in the selected Chromatogram View.

Adding Chromatogram Traces with the Auto Filter Feature

Use the Auto Filter button to repopulate an existing Chromatogram view or add the first Chromatogram view with these possible characteristics:

- A plot showing the chromatogram without any scan filters
- Plots for each scan filter applied to the chromatogram up to the number of scan filters in the data file or the maximum number specified on the Default Workspace Options page, whichever is fewer

Note To manually add and delete traces in a Chromatogram view, see [Adding Chromatogram Traces](#).

❖ **To specify the maximum number of chromatogram traces to display**

1. Open the Workspace Options page (see [Default Workspace Options Page](#)).

3 Reviewing Chromatographic Data

Deleting Chromatogram Traces

2. In the #Number of Auto Filter box, type an integer from **1** to **500**.

The default maximum number of traces is 8.

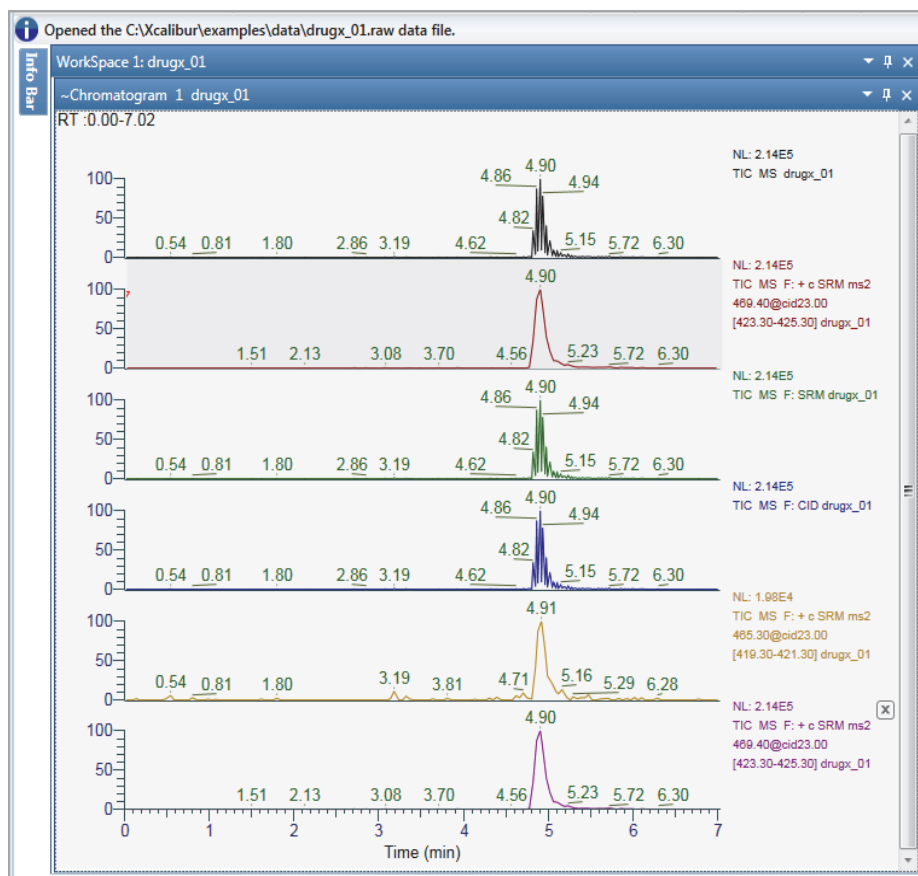
❖ **To automatically add plots for all the scan filters to a Chromatogram view**

In the Workspace Options toolbar, click **Auto Filter**.

The application populates the Chromatogram view with an unfiltered plot (No Filter) and a plot for every scan filter up to the number specified on the Default Workspace Options page of the Default Options Configuration dialog box.

Figure 22 shows the result of clicking the Auto Filter button for drugx_01.raw in the *drive:\Xcalibur\examples\data* folder.

Figure 22. Chromatogram view for a data file with five scan filters and one unfiltered TIC



Deleting Chromatogram Traces

To delete chromatogram traces or undo the most recent deletion of a chromatogram trace in a Chromatogram view, follow these procedures as needed:

- To delete a trace by clicking the delete icon

- To delete a trace by using the Chromatogram Ranges view
- To undo the most recent deletion of a chromatogram trace

Note Deleting a chromatogram trace makes any spectrum that is linked to it available for linking to another trace—that is, the text appended to the Spectrum view’s title bar changes from CxYy to ‘Select Trace to Link’. Deleting all the traces in a chromatogram view closes the view.

❖ **To delete a trace by clicking the delete icon**

1. In the Chromatogram view, point to the name of the trace that you want to delete.

A delete icon appears.


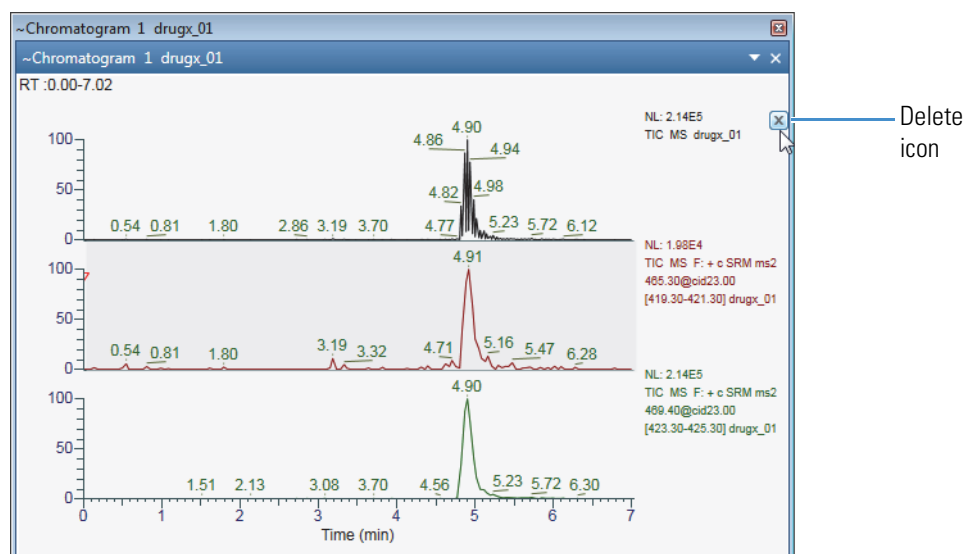

Figure 23 shows a Chromatogram view with three stacked traces. The pointer is on the delete icon, , for the first trace.

Figure 23. Chromatogram view with three stacked traces



2. Click the delete icon, , to delete the trace.

❖ **To delete a trace by using the Chromatogram Ranges view**

- Clear the trace’s associated check box in the Display column.

—or—

- a. Select the row, and then press the **Delete** key.
- b. At the prompt, click **Yes**.

❖ **To undo the most recent deletion of a chromatogram trace**

Right-click the Chromatogram view and choose **Undo Delete Chromatogram**.

Defining a Chromatogram Trace from the Chromatogram Ranges View

The following parameters define the range for a chromatogram trace—raw data file, detector, scan filter, trace type, and mass range for the Mass Range trace type. After you open a raw data file (see [To open a raw data file or a sequence file](#)), you use the Chromatogram Ranges view to select the parameter settings for each chromatogram plot.

You can copy the rows in the Chromatogram Ranges view to the Clipboard, and you can copy the data from a CSV file into the Chromatogram Ranges view.

To define the chromatogram ranges, follow these topics as needed:

- [Manually Defining Chromatogram Ranges](#)
- [Defining Chromatogram Ranges by Using a Spreadsheet File](#)

Manually Defining Chromatogram Ranges

To specify the range for each chromatogram trace by manually entering the parameter settings in the Chromatogram Ranges view, follow these procedures:

- [To manually specify the ranges of the chromatogram traces](#)
- [To specify smoothing for a plot](#)
- [To specify reference \(fixed\) plots](#)

For information about setting a delay time or using the plot operators, see [Chromatogram Ranges View](#). For information about adding multiple rows with the same settings, see [Using the Fill Down Feature](#).

❖ To manually specify the ranges of the chromatogram traces

1. In the Workspace, select the Chromatogram view to define.
2. In the Workspace Options toolbar, click **Chromatogram Ranges**.

The Chromatogram Ranges view opens. Its title bar specifies the linked Chromatogram view.

Note Each row in the Chromatogram Ranges view defines one trace in the selected Chromatogram view.

3. For each trace to display in the Chromatogram view, select another check box in the Display column.
4. In the File Name column, select the raw data file for each trace.
The available detector types depends on the experimental data in the raw data file.
5. In the Detector Type column, select the detector.

Other selections, such as Filter and Trace Type, depend on the selected detector.

Detector	Data type
MS	Mass spectrometry data The raw data files generated by an LC/MS or LC/MS/MS experiment typically contain multiple scan filters to optimize data collection. Use the remaining columns to define the trace range.
MS Trending	Instrument status readings for the mass spectrometer Use the Trace Type column to specify the specific status readback. For more information, see Setting up Instrument Status Traces .
UV	Data from an analog detector Use the Trace Type column to define the trace.
PDA	Wavelength scan data from a photodiode array detector
A/D Card	Data from an analog-to-digital converter

6. For MS data, select a scan filter from the Filter list or from the Scan Filter page.

Tip To display the Scan Filter page in the Info Bar, in the Workspace Options toolbar, click **Scan Filters**.

7. In the Trace Type column, do one of the following:

- For UV data, select the channel.
- For PDA data, select the trace type to determine the signal intensity for each measured time point.

Trace type	Effect
Total Scan	Plots the total signal.
Wavelength Range	Plots the total signal for the wavelength range that you specify in the Ranges box. Continue at step 10 .
Spectrum Maximum	Plots the maximum signal.

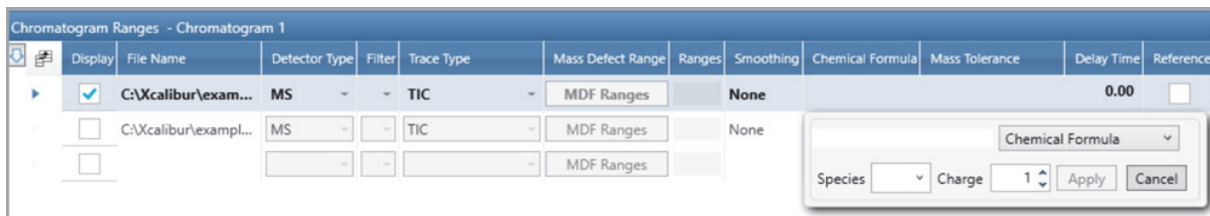
- For A/D data, select the card.
- For MS data, select **TIC, Mass Range, Base Peak, Neutral Fragment, or Mass Defect Filter**.

Note Entering a chemical formula or peptide sequence in the Chemical Formula column automatically changes the trace type to Mass Range and populates the Range column with a calculated m/z value.

3 Reviewing Chromatographic Data

Defining a Chromatogram Trace from the Chromatogram Ranges View

- For MS Trending data, select the status readback.
8. For MS data and the Mass Range, Base Peak, or Neutral Fragment trace type, specify the mass range by doing one of the following:
- In the Ranges box, type the mass range. Use a hyphen to define the mass range.
- or–
- a. In the Chemical Formula column, click the table cell to open a dialog box.



- b. Select **Chemical Formula** or **Peptide**.
- c. Type a chemical formula or a peptide sequence.

Use the IUPAC nomenclature (periodic table symbols) for chemical formulas. Use the one-letter amino acid abbreviations for peptide sequences. See [Appendix C, “One- and Three-Letter Abbreviations for Amino Acid Residues.”](#)

- d. Select an adduct species and a charge.
- e. Click **Apply**.

The application populates the Chemical Formula column with the specified formula or peptide sequence. It also populates these columns:

- Trace Type: Mass Range
- Ranges: m/z value of specified ion
- Mass Tolerance: 0.05 amu for ion trap data or 5 ppm for FTMS data

The EIC trace for the specified mass range appears in the Chromatogram view.

9. For the Mass Range trace type, in the Mass Tolerance column, specify the mass tolerance and the units.
10. For PDA data and the Wavelength Range trace type, in the Ranges box, type the wavelength range. Use a hyphen to define the wavelength range.

❖ To specify smoothing for a plot

1. In the Smoothing column, select one of these smoothing algorithms: **None**, **Gaussian**, or **Moving Mean**.
2. For Gaussian and Moving Mean smoothing, select the smoothing level from a list of odd integers: **3**, **5**, **7**, **9**, and **11**.

The smoothing level increases as the selected value increases.

❖ To specify reference (fixed) plots

Note A reference plot is a fixed plot—that is, the application does not replace a reference plot when you compare raw data files in a sequence set.

Select the check box in the Reference column.

Defining Chromatogram Ranges by Using a Spreadsheet File

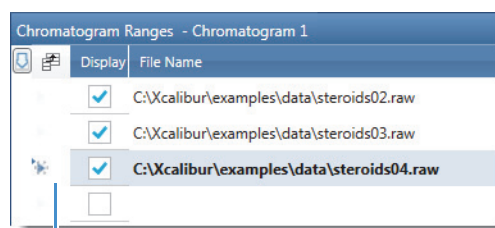
For information about editing the columns in the Chromatogram Ranges view, see [Chromatogram Ranges View](#).

To add defined chromatograms by using the data in a spreadsheet file, follow this procedure.

❖ To specify the ranges of the chromatogram traces by using data in a spreadsheet file

1. Copy rows from the Chromatogram Ranges view to a spreadsheet file as follows:
 - a. In the Workspace Options toolbar, click **Chromatogram Ranges**.
 - b. Click in the area to the left of a row to select the row to copy ([Figure 24](#)).

Figure 24. Example pointer location to copy a row



Click this area.

Tip Use the CTRL key to select non-adjacent rows. Use the SHIFT key for adjacent rows, or drag the pointer across a group of rows.

- c. Press CTRL+C.
- d. Open the Excel file and select the starting cell that you want to paste into.
- e. Press CTRL+V, and then save the file.

In the Excel file, the pasted rows appear to the right and down from the starting cell.

Tip To copy the entire contents of the Chromatogram Ranges view to a CSV file, click **Selection As** in the Workspace Options toolbar.

2. Open the CSV file and edit its contents. Do not change the column format.
3. Paste rows from a spreadsheet file into the Chromatogram Ranges view, as follows:

3 Reviewing Chromatographic Data

Displaying Mass Defect Filtered Chromatogram Traces

- a. In the spreadsheet file, select the rows that you want to copy. Make sure that this block is highlighted and has the same number of columns as in the Chromatogram Ranges view.
- b. Click **Copy** or press CTRL+C.
- c. In the Chromatogram Ranges view, select the row where you want to paste the data.
- d. Press CTRL+V.

The product-name application overwrites the selected row and the rows below it with the copied rows.

Displaying Mass Defect Filtered Chromatogram Traces

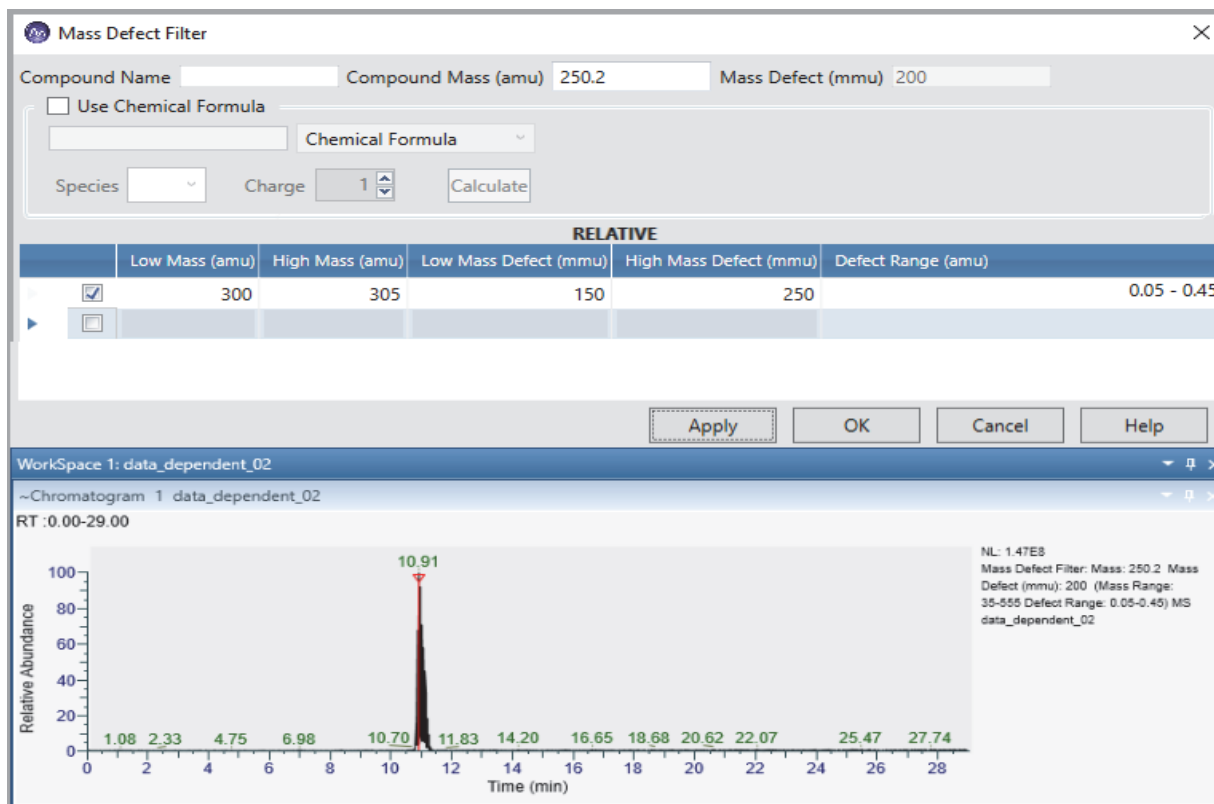
Mass defect is the difference between the accurate mass of an element (or compound) and its closest integer value (nominal mass). It might be positive (larger than the nominal mass) or negative (smaller than the nominal mass). To use MDF, you must use high mass accuracy data from high resolution analyses.

The Multiple Mass Defect Filter (MMDF) in FreeStyle is a data filtering method that removes ions with the mass defect outside the defined mass defect window. The filtering of ions can be based on the mass defect of the parent compound and its modifications. In FreeStyle, you can combine the results from multiple mass defect filters.

Mass defect filter is applied to remove the ions that fall outside the defined molecular weight range, and those that lies within the range but exceeds the expected mass defect range. This method lets you to focus on the analysis of potential drug metabolites.

[Figure 25](#) shows an example of the Mass Defect Filtered (MDF) Chromatogram Trace.

Figure 25. Mass Defect Filtered Chromatogram



The mass defect filtered chromatogram displays the sum of intensities of multiple centroid masses within the specified mass defect range.

Table 12 displays the parameters for calculating the mass defect filtered chromatogram trace.

Table 12. Mass Defect Filter Parameters (Sheet 1 of 2)

Parameter	Description
Compound Name	Specifies the name of the parent compound.
Compound Mass (amu)	Specifies the exact mass or chemical formula. You can also leave the column empty for calculation. <ul style="list-style-type: none"> • Format: 9999.9999 • Compound Mass: 0.0000 and/or empty. If you add the chemical formula, the Compound Mass box becomes non-editable.
Mass Defect (mmu)	Displays the first three decimal values of the compound mass, rounded to the third decimal place.

3 Reviewing Chromatographic Data

Displaying Mass Defect Filtered Chromatogram Traces

Table 12. Mass Defect Filter Parameters (Sheet 2 of 2)

Parameter	Description
Use Chemical Formula	Select to calculate the mass defect based on the chemical formula.
Chemical Formula	Specify the chemical formula used for calculation.
Species	Specify the element.
Charge	Specify the charge.
Low Mass (amu)	Displays a low mass defect range. If the compound mass is given, the value is relative to the compound mass. If the compound mass is empty or zero, the value is an absolute value used for the mass defect filter calculation.
High Mass (amu)	Displays high mass defect range. If the compound mass is given, the value is relative to the compound mass. If the compound mass is empty or zero, the value is an absolute value used for the mass defect filter calculation.
Low Mass Defect (mmu)	Three digits in mmu. The valid range is 0-500. Default: 50 mmu If the compound mass is empty or zero, the low mass defect value is an absolute value. If the compound mass is given, the value is a relative value used for calculation.
High Mass Defect (mmu)	Three digits in mmu. The valid range is 0-500. Default: 50 mmu If the compound mass is empty or zero, the High Mass Defect value is an absolute value. If the compound mass is given, the value is a relative value used for calculation.
Defect Range (amu)	Low Mass Defect -- High Mass Defect

Applying Mass Defect filter in a Chromatogram

❖ To apply mass defect filter

1. Select the Chromatogram view, and then click **Chromatogram Ranges**.

-or-

2. In the Workspace Options toolbar, click **Chromatogram Ranges**.
3. In the Trace Type list, select **Mass Defect Filter**.
4. In the Mass Defect Range, click **MDF Ranges**.
5. In the Mass Defect Filter dialog box, enter the following parameters:
 - Compound Name: Specify the compound name or leave the column empty.
 - Compound Mass (amu): Specify the exact mass of the compound or leave the column empty.
 - Use Chemical Formula: Select if you want to use the chemical formula. The compound mass and mass defect automatically populates when you use chemical formula for calculation.
 - Chemical Formula: Type the chemical formula that you want to use for calculation.
 - Species: Select the species from the dropdown list.
 - Charge: Select the charge from the dropdown list.
 - Select the check box next to the first row.
 - Enter the low mass, high mass, low mass defect, and high mass defect.
 - Select the check box next to each row, and then repeat entering the multiple mass range and mass range defect parameters to create a Multiple Mass Defect Filtered trace.
6. Click **Apply**, and then click **OK**.

The mass defect filtered chromatogram displays.

Relative and Absolute Mass Range Calculation

Mass defect filtering of a chromatogram trace is calculated as either relative or absolute. If you type a compound mass value in the Mass Defect Filter dialog box or use a chemical formula, the mass defect range is calculated relative to the compound mass and its mass defect. If you type zero or leave the compound mass column empty, the mass defect value used for calculation is an absolute range based on the low and high mass, and the low and high mass defect that you have entered.

3 Reviewing Chromatographic Data

Displaying Mass Defect Filtered Chromatogram Traces

Figure 26. Relative and Absolute Mass and Mass Defect Calculation

ABSOLUTE					
	Low Mass (amu)	High Mass (amu)	Low Mass Defect (mmu)	High Mass Defect (mmu)	Defect Range (amu)
<input checked="" type="checkbox"/>	300	305	50	50	-0.05 - 0.05
<input checked="" type="checkbox"/>	300	305	50	50	-0.05 - 0.05
<input type="checkbox"/>					

❖ **To calculate the relative and absolute mass and mass defect values**

1. In the Chromatogram Ranges view, select:
 - Detector Type: **MS**
 - Trace Type: **Mass Defect Filter**
 2. In the Mass Defect Range column, click **MDF Ranges**.
- The Mass Defect Filter dialog box opens.
3. In the Compound Mass (amu) column, type the compound mass value for relative calculation

-or-

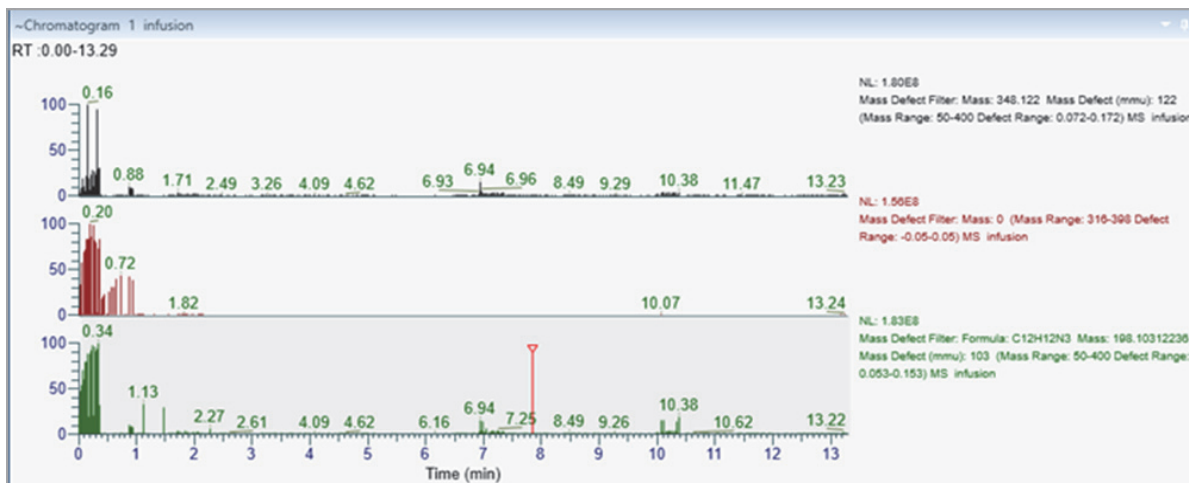
Type zero or leave the column empty for absolute calculation.

If you type a compound mass value, the Mass Defect columns displays the first three decimal values of the compound mass, rounded to the third decimal place. For example, if you type 348.122, the application automatically populates the mass defect value as 122.

4. Select each row and type the low mass and high mass range values (for example, 300-305) and the low mass and high mass defect values.
5. Click **Apply**, and then click **OK**.

Figure 27 indicates how the relative and absolute calculation of mass and mass defect values differ for different chromatogram traces in FreeStyle.

Figure 27. Relative and absolute mass defect and mass range changes for different chromatogram



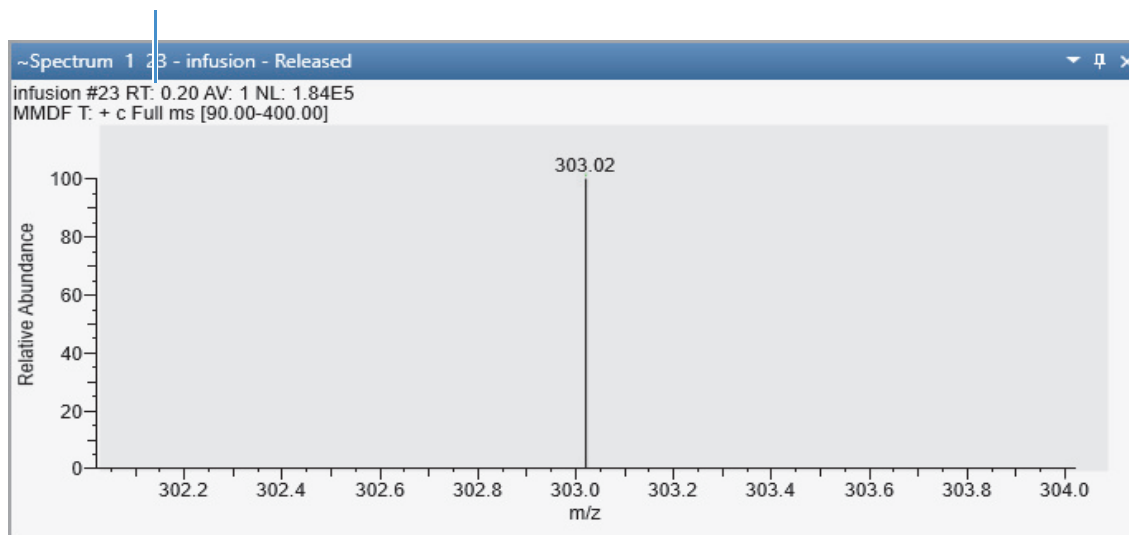
Linking a Spectrum View to an MDF Chromatogram Trace

When you select an MDF chromatogram, the Spectrum view displays mass spectral peaks corresponding to the mass defect filters applied to the chromatogram. If a single scan is selected or linked to an MDF chromatogram trace, the active or the linked spectrum:

- Displays the spectrum title as MMDF T: +c d Full ms2...
- Updates the filtered mass data corresponding to the active MDF chromatogram trace.
- Updates the Spectrum list with the filtered mass data corresponding to the active MDF chromatogram trace.
- Enables the Write to .RAW feature.
- Deactivates the NIST search and mzVault search options.

Figure 28. Spectrum linked to the MDF Chromatogram trace

Linked to MDF chromatogram trace



❖ To link a spectrum view to an MDF chromatogram trace

1. In the Chromatogram Ranges view, select:
 - Detector Type: **MS**.
 - Trace Type: **Mass Defect Filter**.
2. In the Mass Defect Range column, click **MDF Ranges**.

The Mass Defect Filter dialog box opens.

3. In the Mass Defect Filter dialog box, enter the low mass, high mass range, low mass defect, and high mass defect values.
4. Click **Apply**, and then click **OK**.

The title of the Spectrum view describes the linked MDF chromatogram trace in the following format:

MMDF T: +c d Full ms

Setting up Instrument Status Traces

You can set up traces for the MS detector's status readbacks and compare these traces to possible disturbances in the chromatographic data.

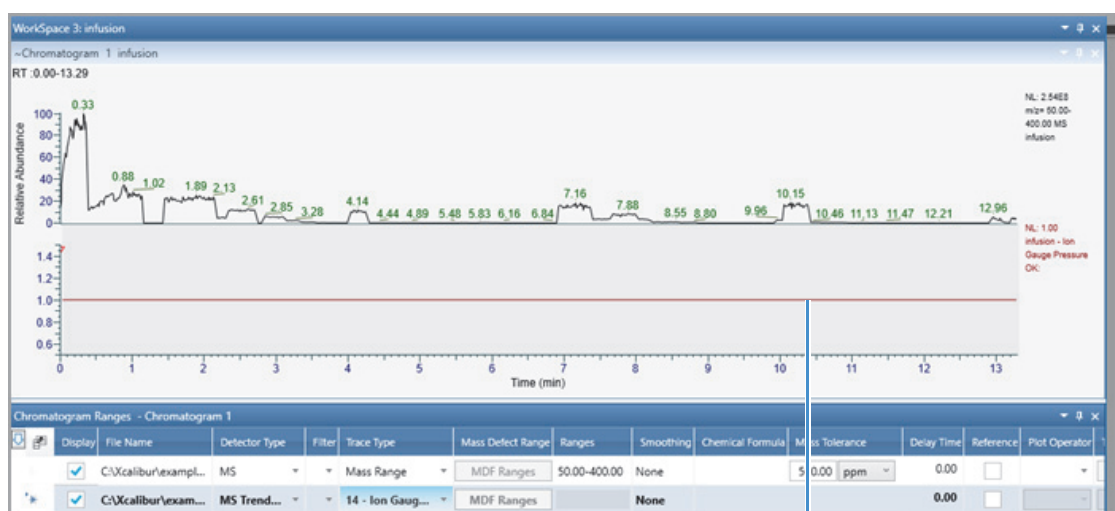
❖ To add instrument status traces to the Chromatogram view

1. Select the Chromatogram view where you want to add the trace.

2. In the Workspace Options toolbar, click **Chromatogram Ranges**.
3. For each status readback of interest, do the following:
 - a. Add a plot to the Chromatogram view (see [To add a chromatogram trace by using the Chromatogram button](#) or [To add chromatogram traces by using the Chromatogram Ranges view](#)).
 - b. In the Detector Type list of the Chromatogram Ranges view, select **MS Trending**.
 - c. In the Trace Type list, select a readback parameter.

Figure 29 shows the ion transfer tube temperature readback as a function of time.

Figure 29. Status readback displayed in the Chromatogram view



Selected readback parameter

Readback trace for the ion gauge pressure during the 13 min run

Displaying an EIC Trace by Using the Spectrum Toolbox

You can use the Spectrum toolbox to define an EIC trace when the spectrum is linked to the chromatogram.


The Spectrum toolbox has two icons for adding EIC traces to the Chromatogram view—the EIC Mass icon and the EIC Range icon. To add a plot for multiple discrete m/z values, use the EIC Mass icon. To add a trace for contiguous m/z ranges, use the EIC Range icon.

3 Reviewing Chromatographic Data

Displaying an EIC Trace by Using the Spectrum Toolbox

Tip To display an EIC trace for a single mass-to-charge value, double-click the mass-to-charge label for the peak in the Spectrum view. The Spectrum view displays the EIC trace for the selected mass-to-charge ratio ± 1 Da.

❖ To display an EIC trace for multiple discrete m/z values

1. Select the Spectrum view and open the Spectrum toolbox.
2. Click the **EIC Mass** icon, .

If the toolbox was unpinned, the application pins it so that it remains visible.

3. In the Spectrum view, click the mass-to-charge label for each peak that you want to use for the EIC trace.

A red vertical marker indicates a selected peak. [Figure 30](#) shows the selection of three mass peaks.

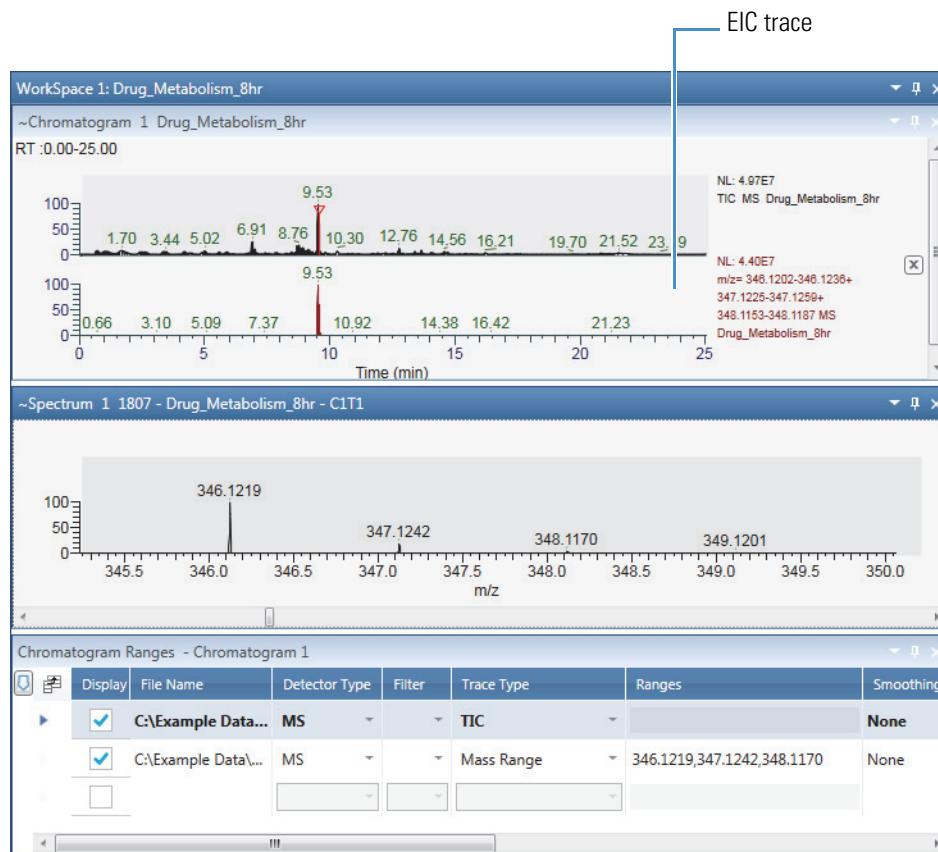
Figure 30. Spectrum view with the selection of three mass peaks



4. Click the **EIC Mass** icon again to create an EIC trace from the selected masses.

A new EIC trace appears at the bottom of the Chromatogram view ([Figure 31](#)), and the selected masses appear in the Ranges column in the Chromatogram Ranges view.

Figure 31. EIC trace created from three mass peaks



❖ **To specify the mass ranges for an EIC trace**

1. Open the Spectrum toolbox.
2. Click the **EIC Range** icon, .

If the toolbox was unpinned, the application pins it so that it remains visible.

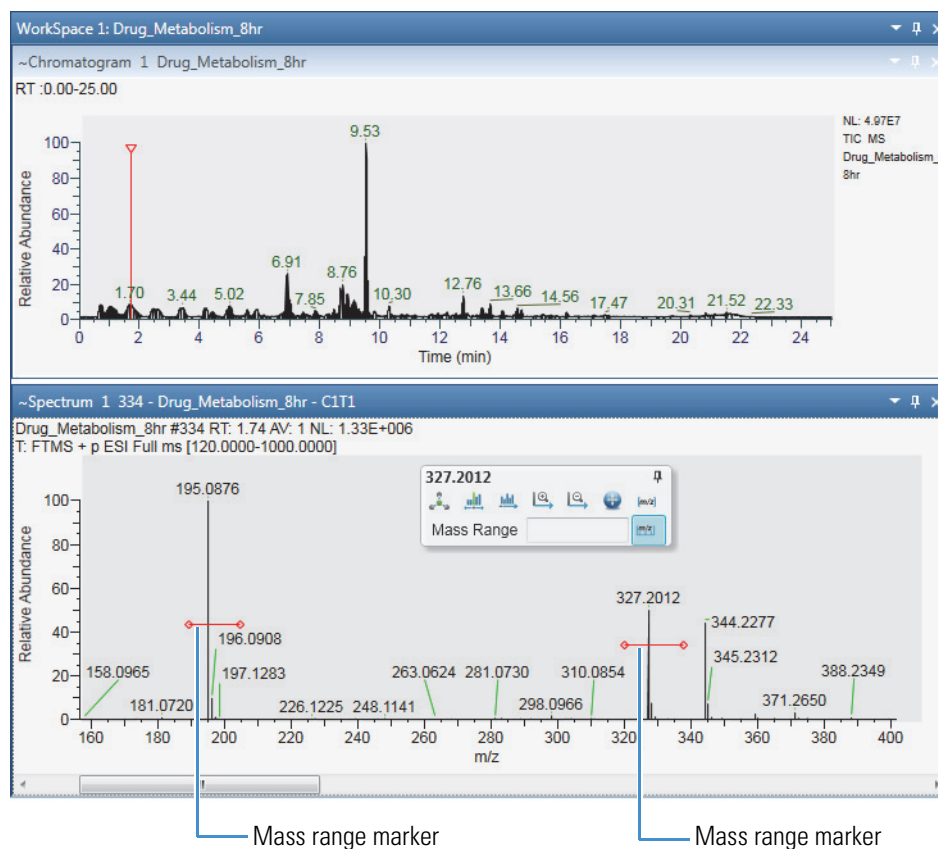
3. In the Spectrum view, drag the pointer horizontally through a contiguous or noncontiguous mass range to select the mass range that you want to use for the EIC trace.

A horizontal red line with start and end markers indicates the start and end of the selected range (Figure 32).

3 Reviewing Chromatographic Data

Displaying an EIC Trace by Using the Spectrum Toolbox

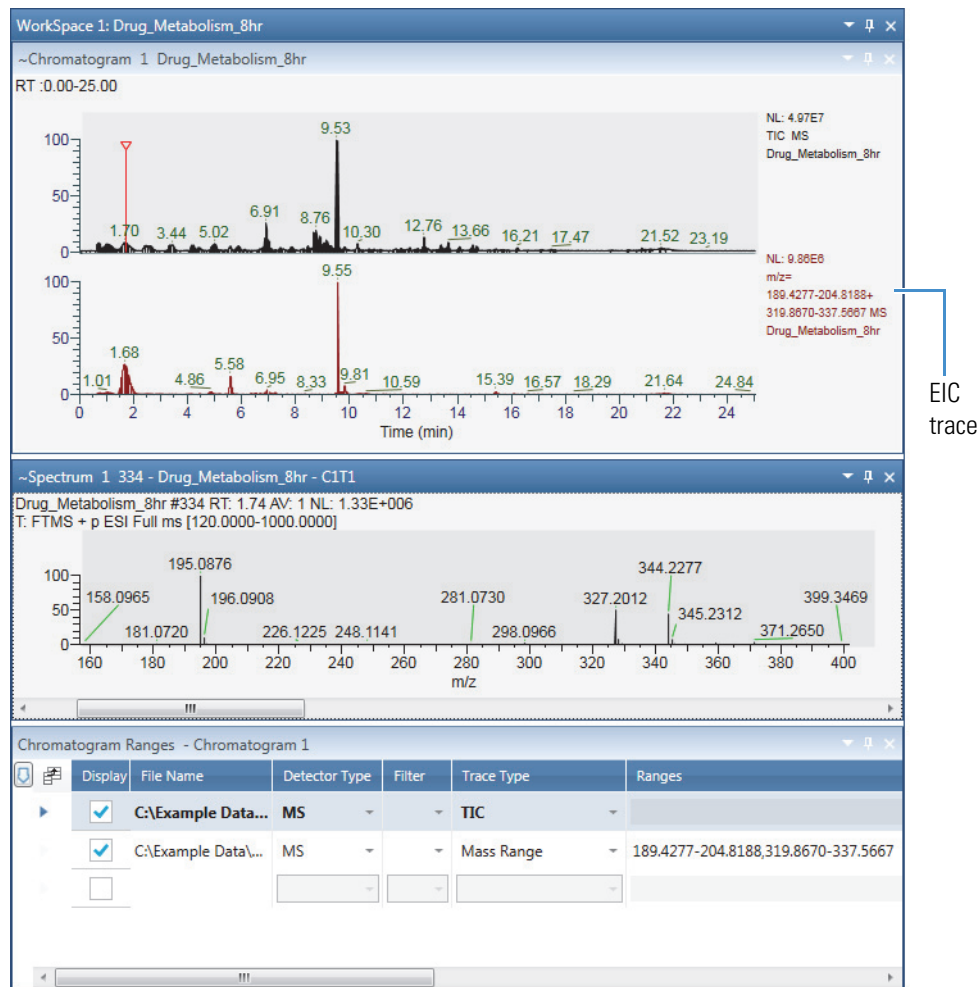
Figure 32. Two mass ranges selected in the Spectrum view



4. To create an EIC trace from the selected mass ranges, click the **EIC Range** icon again.

The selected mass ranges appear in the Ranges column of the Chromatogram Ranges view, a new EIC trace appears at the bottom of the Chromatogram view, and the application unpins the toolbox. [Figure 33](#) shows the EIC trace for two noncontiguous mass ranges.

Figure 33. EIC trace for two noncontiguous mass ranges



Using the Scan Filters Page to Display a Filtered Chromatogram

Use the Scan Filters page or the Filters column of the Chromatogram Ranges view to display filtered chromatograms.

❖ To display a filtered chromatogram

1. Open the raw data file of interest.
2. In the Workspace Options toolbar, in the Filter area, click **Scan Filters**.

The Scan Filters page appears in the Info Bar. By default, the Time Range (min) box contains an asterisk and the Track check box is clear. The asterisk means that the time range is set to the full length of the acquisition time for the raw data file.

3. (Optional) To limit the displayed time range to a portion of the acquisition time, do one of the following:

3 Reviewing Chromatographic Data

Setting Up the Display Options for a Chromatogram Trace

- In the Time Range (min) box, type the retention-time range as follows:
Start time-End time
- or–
- Select the **Track** check box, and then in the Chromatogram view, drag the pointer across the retention time of interest.
4. Select or type a filter in the filter list.

The filtered chromatogram appears in the Chromatogram view.

Setting Up the Display Options for a Chromatogram Trace

For the Chromatogram view, you can use the buttons in the Chromatogram – Display Options toolbar to stack or overlay chromatogram traces, format and label the chromatogram peaks and axes, specify how the application normalizes the chromatogram traces, and change the *y*-axis title and scale.

Note Except for the Decimal Places button, all the buttons and icons are available on the Chromatogram – Display Options toolbar. The Decimal Places button becomes available after you add the Retention Time label to your chromatograms.

Although the Peak Area, Signal To Noise, and Height buttons are always available, the application only adds these labels to integrated peaks.

To customize the display options for chromatogram plots, see these topics:

- [Formatting Chromatogram Plots](#)
- [Labeling Chromatogram Peaks or Local Maxima](#)
- [Normalizing Chromatogram Traces](#)
- [Changing the Y-Axis Title and Scale of the Chromatogram View](#)

Formatting Chromatogram Plots

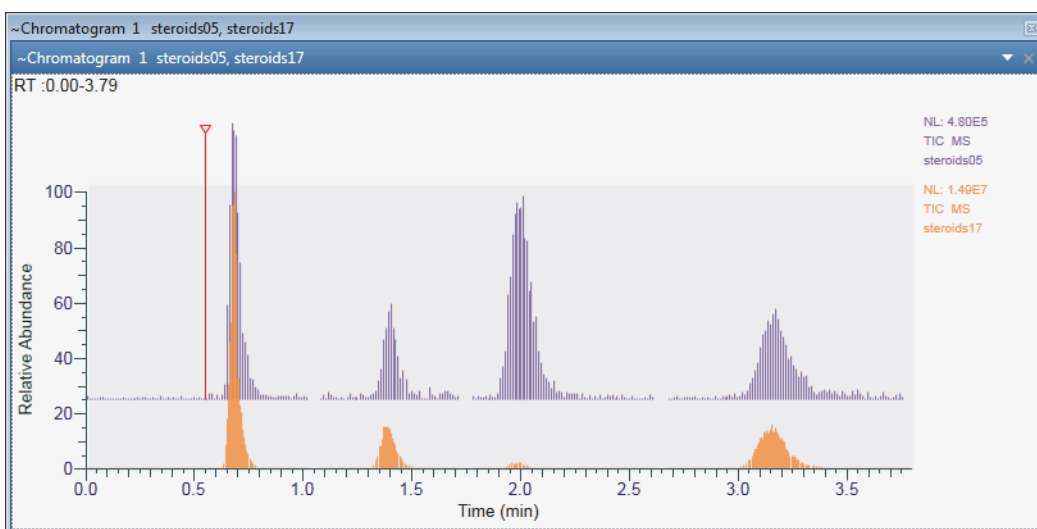
Use the Format area of the Chromatogram – Display Options toolbar to format the chromatogram plots. You can display chromatogram traces as point-to-point graphs (line graphs) or as discrete sticks for each scan. When the Chromatogram view contains multiple plots, you can stack the plots or overlay them.

With the default factory settings, the application displays multiple traces as stacked line graphs, with the top graph plotted in black. The Stack, Overlay, Point-to-Point, and Stick format options apply to all the plots. To select different colors for overlaid plots, select the color for each active plot before you overlay them. Use the Elevation slider to change the percentage overlay of the plots. Setting the slider full left overlays the traces to the same elevation on the y axis. Use the Skew slider to offset the x axis of the plots. You cannot offset the x axis when the traces are 100% overlaid to the same elevation on the y axis.

For information about all the display options, see [Setting Up the Display Options for a Chromatogram Trace](#).

Figure 34 shows two overlaid stick plots.

Figure 34. Overlaid stick plots



Labeling Chromatogram Peaks or Local Maxima

Use the Labels section of the Chromatogram – Display Options toolbar to label the chromatogram traces. When using the default factory layout, the application labels the retention time of each localized maxima in the chromatogram plots to two decimal places. When you apply a peak detection algorithm or manually draw the chromatographic baseline, the application only labels the retention time of the detected peaks.

For information about all the display options, see [Setting Up the Display Options for a Chromatogram Trace](#).

To label chromatogram traces, follow these procedures:

- To add the peak area, peak height, and signal-to-noise labels to each detected peak
- To remove a label
- To specify the format of the retention time label

3 Reviewing Chromatographic Data

Setting Up the Display Options for a Chromatogram Trace

- To change the number of decimal places for the numeric labels
 - To change the labeling threshold
- ❖ **To add the peak area, peak height, and signal-to-noise labels to each detected peak**

1. Run a peak detection algorithm or manually draw the chromatographic baseline.
2. In the Chromatogram – Display Options toolbar, click the following buttons:
 - To add the peak area label, click **Peak Area**.

Note The letters *MA* or *AA* next to the value indicate manual integration or automatic integration, respectively.

- To add the peak height label, click **Height**.

Note The letters *MH* or *AH* next to the value indicate manual integration or automatic integration, respectively.

- To add the signal-to-noise label, click **S/N Signal To Noise**.

Note The Avalon peak detection algorithm does not calculate a signal-to-noise ratio. For the Genesis and ICIS peak detection algorithms, you can choose to report the root mean square (RMS) noise values. For manual peaks, the application calculates the RMS signal-to-noise ratio.

The letters *SN* next to the value indicate that it is a signal-to-noise ratio.

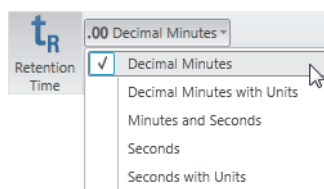
❖ **To remove a label**

In the Chromatogram – Display Options toolbar, click the associated button.

❖ **To specify the format of the retention time label**

Note By default, the Retention Time label is turned on and set to Decimal Minutes.

1. If the Retention Time label is turned off, in the Chromatogram – Display Options toolbar, click **Retention Time**.
2. Select the format for the RT label from the list to the right of the Retention Time button.



❖ **To change the number of decimal places for the numeric labels**

Click **Decimal Places** and select a value from **0** to **5** from the list.

❖ **To change the labeling threshold**

In the Label Threshold box, type a percentage from **0** to **100**.

For a value of 0%, the application labels all the trace maxima or detected peaks. At a value of 100%, the application labels only the highest data point on the trace. When the peak apex of a detected peak falls below this height, the application does not label the peak.

Filling the Chromatogram Peaks

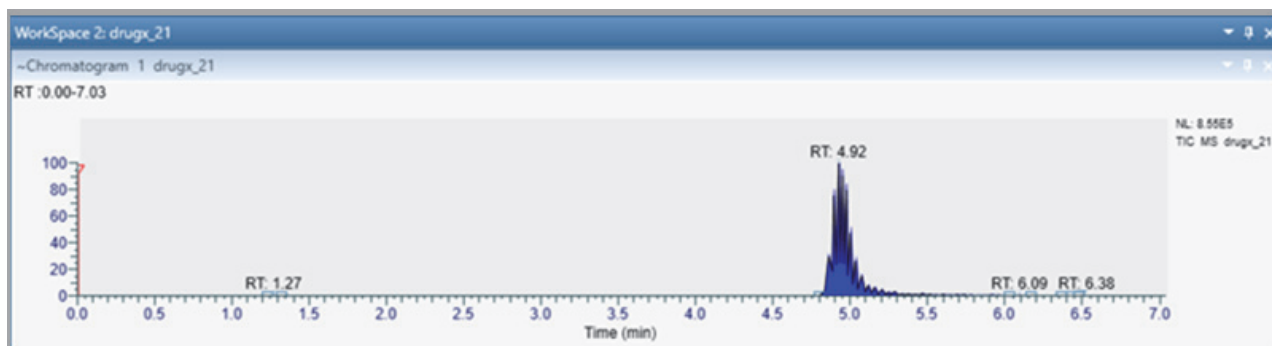
Use the Peak Fill area in the Chromatogram - Displays Options toolbar to color the chromatogram peaks. You can turn on or off the peak colors. The default color is blue.

❖ **To fill the chromatogram peaks**

1. Click the Chromatogram view to make it active.
2. In the Peak Fill area of the Chromatogram – Display Options toolbar, click the color box. The Color dialog box opens with a color palette, where you can select a preset color or customize a color.
3. Open the Peak Detection Info bar page.
4. Select the **Peak Detection Algorithm**, and then click **Apply**.

The peaks are filled with the selected color.

Figure 35. Peak Fill



5. To color the peaks of a PPD peak detection algorithm, select **Apply to PPD Peaks**.
6. To turn the color off, in the Peak Fill area, click **Fill**.

Note The Fill button is a toggle button that you can use to turn the color of the peaks on or off.

Normalizing Chromatogram Traces

Use the Normalization area of the Chromatogram – Display Options toolbar to normalize chromatogram traces. You can apply local or global normalization or turn off normalization.

Changing the normalization type changes the *y*-axis scaling:

- Local normalization scales each trace independently to its largest peak in the displayed time range.
- Global normalization scales each trace to the largest peak across the traces in the displayed time range.
- Turning off normalization scales each trace to the largest peak across the traces in the full time range, regardless of the displayed time range.

Follow these topics:

- [Applying Local Normalization](#)
- [Applying Global Normalization](#)
- [Turning Off Normalization](#)
- [Changing the Y-Axis Range for Normalized Chromatograms](#)

For information about all the display options, see [Setting Up the Display Options for a Chromatogram Trace](#).

Applying Local Normalization

When you apply local normalization (the default setting), the application normalizes the chromatogram traces so that, within the displayed time range, the most intense peak for each trace is set to 100 percent of the *y*-axis scale.

❖ To apply local normalization

1. Click the Chromatogram view to select it.
2. In the Normalization area of the Chromatogram – Display Options toolbar, click **Local**.

[Figure 36](#) shows an example of local normalization. Each trace is independently scaled to its largest peak in the displayed time range, which is the full time range.

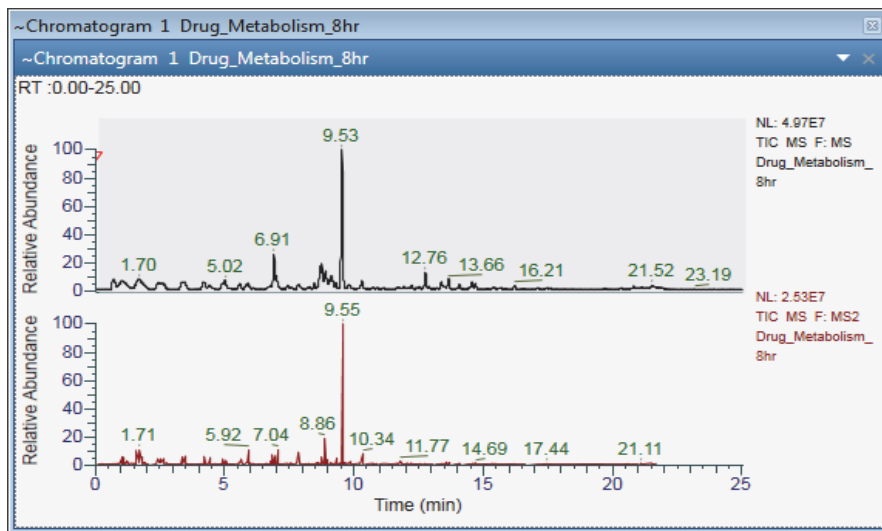
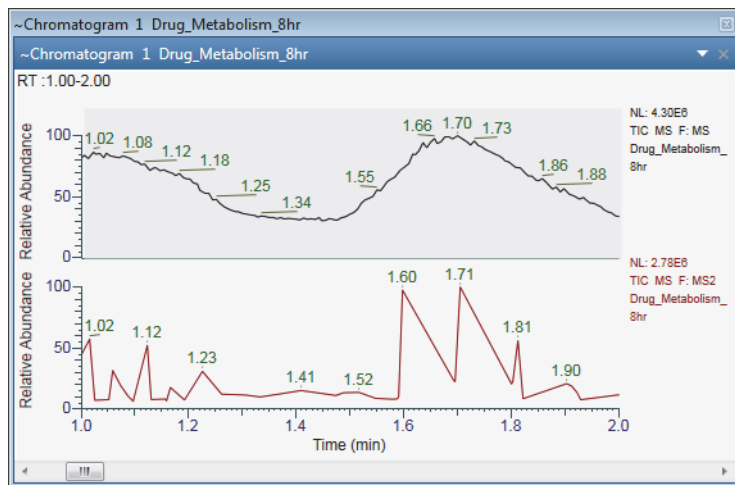
Figure 36. Local normalization with full time range displayed

Figure 37 shows the TIC traces for the same scan filters as those shown in Figure 36, but only the 1 to 2 min time range is displayed. In this case, each trace is independently scaled to its largest peak in the displayed time range of 1 to 2 min.

Figure 37. Local normalization for the 1 to 2 min time range

Applying Global Normalization

For global normalization, the application normalizes the chromatogram traces so that the most intense peak of all traces is scaled to 100% of the y axis.

❖ To apply global normalization

1. Click the Chromatogram view to select it.
2. In the Normalization area of the Chromatogram – Display Options toolbar, click **Global**.

3 Reviewing Chromatographic Data

Setting Up the Display Options for a Chromatogram Trace

Figure 38 shows an example of global normalization. Both traces are scaled to the largest peak across the traces, rather than being independently scaled to the largest peak in each trace.

Figure 38. Global normalization with the full time range displayed

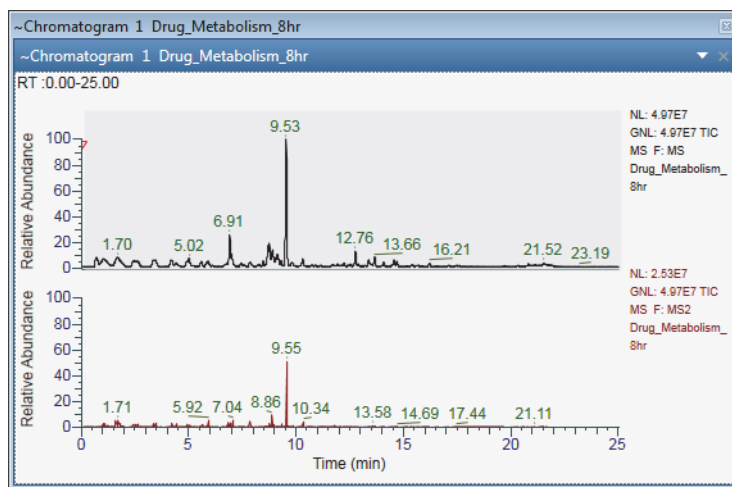
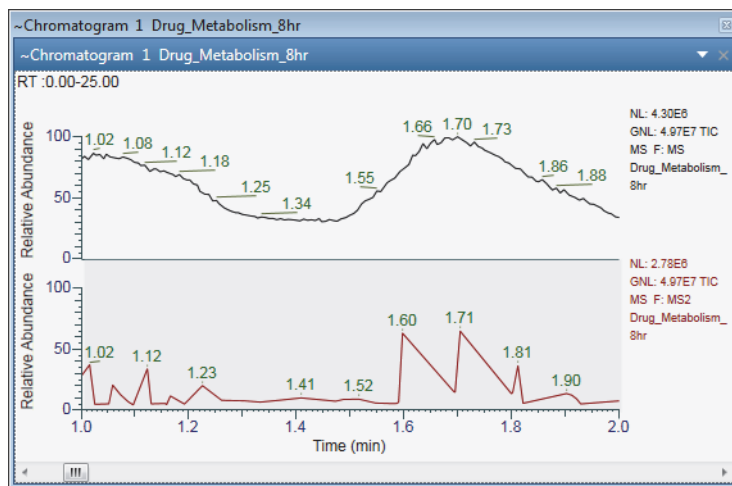


Figure 39 shows the TIC traces for the same scan filters as those shown in Figure 38, but only the 1 to 2 min time range is displayed. In this case, each trace is scaled to the largest peak across the traces in the displayed time range of 1 to 2 min.

Figure 39. Global normalization with the 1 to 2 min time range displayed



Turning Off Normalization

When you turn off normalization, the application scales all the traces by using the maximum and minimum absolute values in the full time range, regardless of the time range displayed.

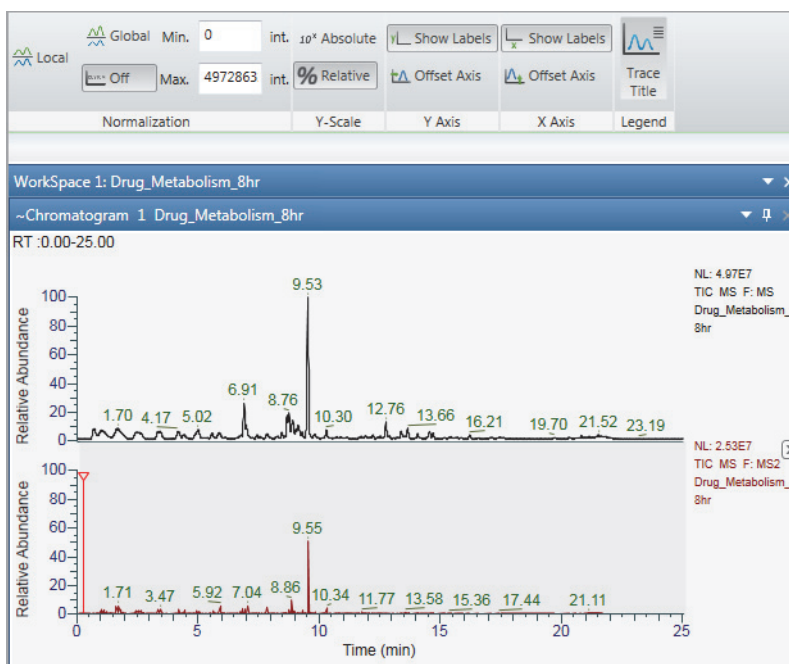
❖ To turn off normalization

1. Click the Chromatogram view to select it.

2. In the Normalization area of the Chromatogram – Display Options toolbar, click **Off**.

A comparison of [Figure 40](#) and [Figure 41](#) shows the effect of turning off normalization. Regardless of the time range displayed, the traces are scaled to the largest peak across the traces in the full time range. In addition, the Max box displays the absolute intensity of the largest peak (apex height).

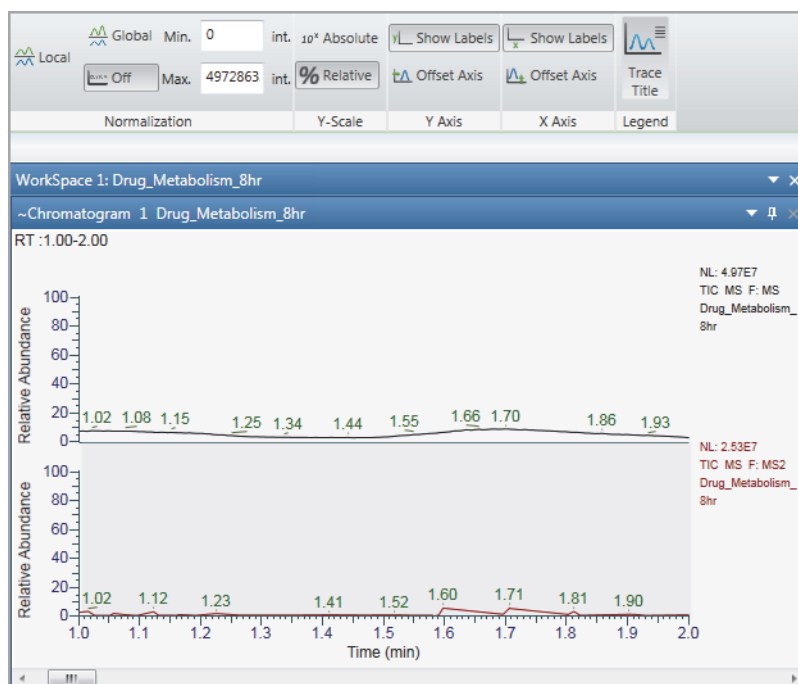
Figure 40. Normalization turned off with the full time range displayed



3 Reviewing Chromatographic Data

Setting Up the Display Options for a Chromatogram Trace

Figure 41. Normalization turned off with the 1 to 2 min time range displayed



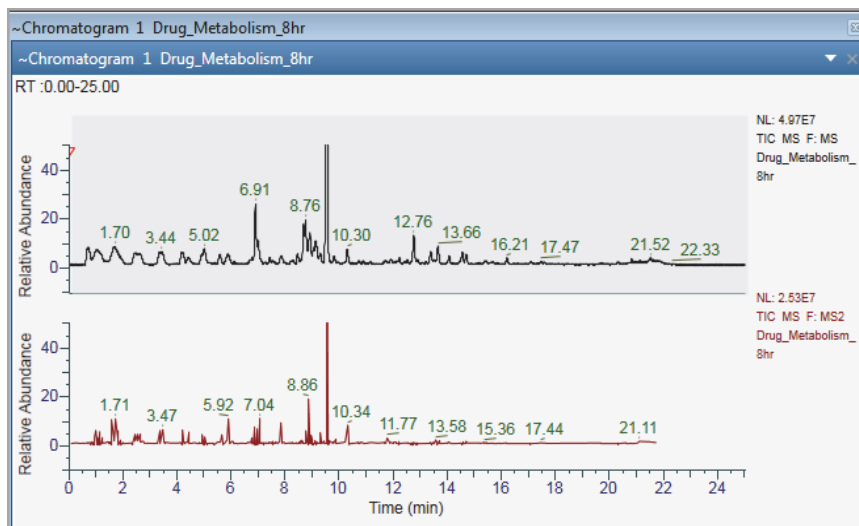
Changing the Y-Axis Range for Normalized Chromatograms

Use the Min and Max boxes in the Normalization area of the Chromatogram – Display Options toolbar to change the y-axis range from its default of full zoom (0–100%) to another zoom level.

Note Regardless of whether the y-axis scaling (Y Scale) is set to Absolute or % Relative, when normalization is turned on (local or global), the Min and Max boxes accept positive and negative values as percentages of the full y-axis scale. The Max value must be greater than the Min value.

For example, to zoom in on the –10 to 50% portion of the y axis, type –10 in the Min box and 50 in the Max box.

Figure 42 shows the result.

Figure 42. Modified y -axis range of -10 to 50% 

Changing the Y-Axis Title and Scale of the Chromatogram View

Use the y -axis and y -scale area of the Chromatogram – Display Options toolbar to change the y -axis title and scale of the Chromatogram view.

Follow these topics:

- [Changing the Y-Axis Title](#)
- [Changing the Y-Axis Scale](#)

Changing the Y-Axis Title

Enable the **Separate Axis** button in the y -axis area to change the y -axis title of the individual traces of the Chromatogram view.

❖ To change the Y-Axis title of the Chromatogram view

1. Click the Chromatogram trace to select it.
2. In the y -axis area of the Chromatogram – Display Options toolbar, click the **Separate Axis**.
3. You can type the y -axis title of your choice in the text box next to the Separate Axis.

The new title appears in the active Chromatogram trace in the Chromatogram view.

Note When you add a new Chromatogram trace, the y -axis title of the active Chromatogram trace reflects on the new trace.

Changing the Y-Axis Scale

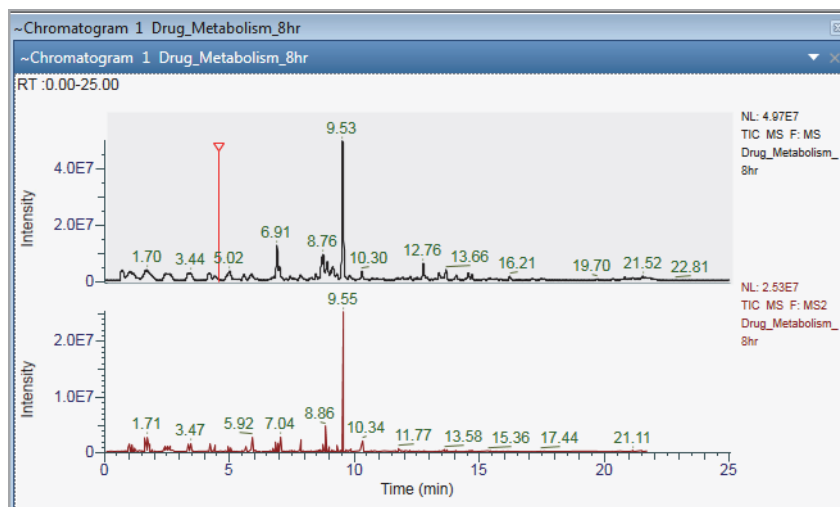
Relative abundance is the default y -axis scale (units) for the Chromatogram view. You can change the y -axis units of individual traces of the Chromatogram view by enabling the Separate Axis option in the y -axis area. Use the Default Workspace Options to change the y -scale units for each detector type. For information on how to set the units according to the detector type, see [Default Workspace Options Page](#).

❖ To change the Y-Axis units from relative abundance to absolute

1. Click the trace in the Chromatogram view to select it.
2. In the y -scale area of the Chromatogram – Display Options toolbar, click **|Y|Absolute**.
3. You can view the y -axis units in either exponent (10^x) or general form by clicking 10^x or 1.23 on the y -scale area.

[Figure 43](#) shows a Chromatogram view with the y -axis scale set to absolute (where the title is intensity) and normalization set to local.

Figure 43. Units for the y -axis set to absolute intensity



Changing the Zoom Level of a Chromatogram

You can zoom in on or out of a region of a chromatogram by using the Zoom Options toolbar, the Chromatogram toolbox, or the mouse pointer.

Follow these procedures:


- To incrementally zoom in and out on the x axis
- To zoom in on a specific section of the x axis
- To incrementally zoom in and out on the y axis

- To reset the zoom level of the *x* and *y* axes

❖ **To incrementally zoom in and out on the *x* axis**

1. Select the view.
2. To zoom in, do the following:
 - Open the Zoom Options toolbar and click **Zoom In X**.


–or–

- Open the Chromatogram toolbox and click the **Zoom In** icon, .

Each click zooms in on a 50% smaller section of the *x* axis—that is, it decreases the displayed range by a factor of 2.

3. To zoom out, do the following:
 - In the Zoom Options toolbar, click **Zoom Out X**.

–or–

- Open the Chromatogram toolbox and click the **Zoom Out** icon, .

Each click increases the displayed time range by 200%.

❖ **To zoom in on a specific section of the *x* axis**

Drag the pointer horizontally across the specific section of the *x* axis.

❖ **To incrementally zoom in and out on the *y* axis**

1. Select the view.
2. To zoom in, open the Zoom Options toolbar and click **Zoom In Y**.

Each click zooms in on a 50% smaller section of the *y* axis. For example, in a Chromatogram view where the *y* axis is set to relative abundance, clicking once zooms in on the 0 to 50% range, clicking twice zooms in on the 0 to 25% range, and so on.


3. To incrementally undo the zoom level, click **Zoom Out Y**.

Each click zooms out to a 200% larger section of the *y* axis.

❖ **To reset the zoom level of the *x* and *y* axes**

- Select the view, open the Zoom Options toolbar, and click **Reset**.
- Right-click the view and choose **Reset Scaling**.

–or–

- Select the view, open the Chromatogram toolbox, and click the **Reset Zoom** icon, .

Automatically Detecting and Integrating Chromatographic Peaks

Use the Peak Detection page in the Info Bar to select and apply any of the automated peak detection algorithms—Genesis, ICIS, Avalon, or PPD. When using the Genesis or ICIS algorithm, you can manually select the chromatographic time range that the application uses to determine the baseline noise level.

Follow these procedures:

- [To display the Peak Detection page](#)
- [To define a peak identification criteria](#)
- [To run any of the peak detection algorithms](#)
- [To run the elemental composition algorithm](#)
- [To reset the zoom level in the Chromatogram view](#)

❖ **To display the Peak Detection page**

1. Open a raw data file (RAW) or a sequence file (SLD).

The Peak Detection page appears in the Info Bar.

2. Click the **Peak Detection** tab.

❖ **To define a peak identification criteria**

1. In the Identification area, from the Returned Peaks dropdown list, select one of the following:
 - **All Peaks:** Select to identify all peaks in the chromatogram based on the selected algorithm for component identification.
 - **Nearest Peak:** Select to identify the nearest peak in the chromatogram within the specified time window.
 - **Highest Peak:** Select to identify the highest peak in the chromatogram within the specified time window.
2. If Nearest Peak or Highest Peak is selected, in the Expected Time (min) box, type or select the expected retention time for the component. The valid range is 0 to 999 minutes.
3. If Nearest Peak or Highest Peak is selected, in the Window (sec) box, type or select the retention time window for the component. The valid range is 1.0 to 999.0 seconds.
4. Specify the component name used in the processing method.

❖ **To run any of the peak detection algorithms**

1. In the list, select one of these algorithms:
 - **Genesis**—Use for mass spectrometry data.

- ICIS—Use for mass spectrometry data.
 - Avalon—Use for UV or PDA data. This algorithm detects negative chromatographic peaks more accurately than the ICIS or Genesis algorithms.
 - PPD—Use to minimize the number of optimization parameters.
2. If the Chromatogram view contains multiple plots, do one of the following:
 - To process all the plots, select the **Apply to All Plots** check box.
 - To process only the selected plot, clear the **Apply to All Plots** check box.
 3. If necessary, modify the parameters settings for the selected algorithm.

For information about the parameters settings for each algorithm, see these topics:

- [Genesis Peak Detection Page](#)
- [ICIS Peak Detection Page](#)
- [Avalon Peak Detection Page](#)
- [PPD Peak Detection Page](#)

For information about selecting the manual noise region for the ICIS or Genesis algorithms, see [Selecting the Manual Noise Region for the Genesis and ICIS Algorithms](#).

4. (Optional) To save the modified settings as the default settings for the selected algorithm, click **Save As Defaults**.

Note You can view the saved settings on the Peak Detection page of the Default Options Configuration dialog box. See [Appendix A, “FreeStyle Default Settings.”](#)

5. Depending on the algorithm, do one of the following:
 - For the Genesis, ICIS, or PPD algorithm, click **Apply**.
 - For the Avalon algorithm, click **Auto Calc Initial Events**. If necessary customize the integration by adding more events to the table. Each time you add an event, the algorithm automatically reintegrates the plots.

After the peak detection finishes, the application does the following:

- Adds the detected peaks to the peak list in the Peaks List view.
- Colors the integrated region of the detected peaks in the Chromatogram view.
- Draws a baseline under the integration region.
- Adds square integration markers at the beginning and end of the integration region. You can change the integration region by dragging the markers.

3 Reviewing Chromatographic Data

Automatically Detecting and Integrating Chromatographic Peaks

Note If the instrument method used to acquire the raw data file specified the compound names of the analytes, the application displays these names above the detected peaks. The instrument control software for the Thermo Scientific TSQ Endura™ and TSQ Quantiva™ mass spectrometers provides this feature.

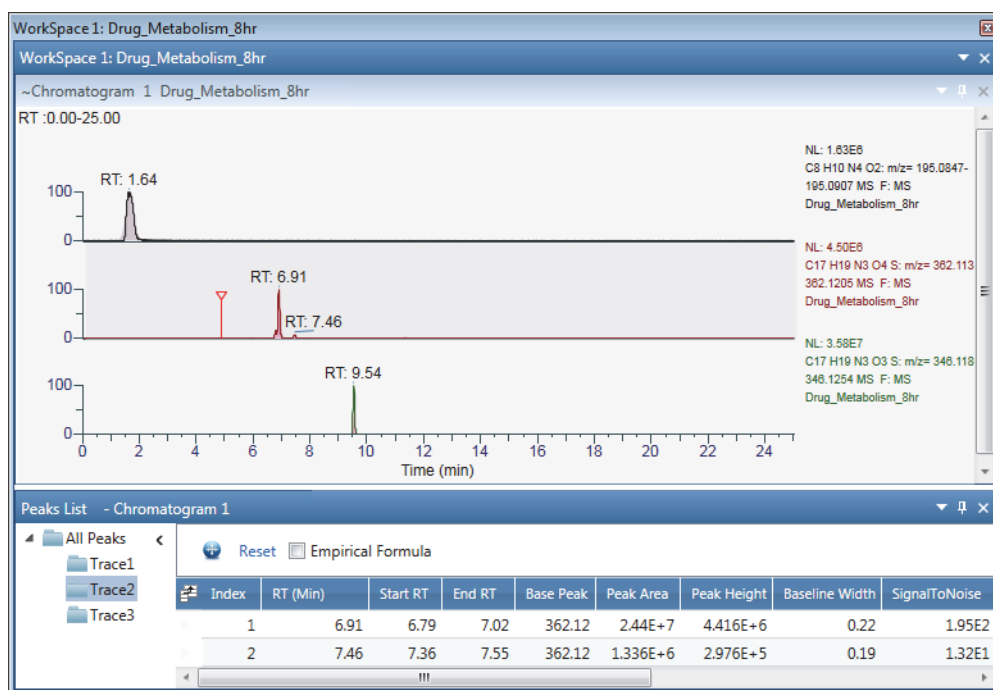
❖ To check the integration for each chromatographic peak

1. If you applied the algorithm to multiple plots, in the Peaks List view, select the plot from the trace list.

Note The Peaks List view defaults to displaying a list of all the peaks (All Peaks) detected across the chromatogram plots.

Figure 44 shows the second trace (Trace 2) selected from a list of three chromatogram plots.

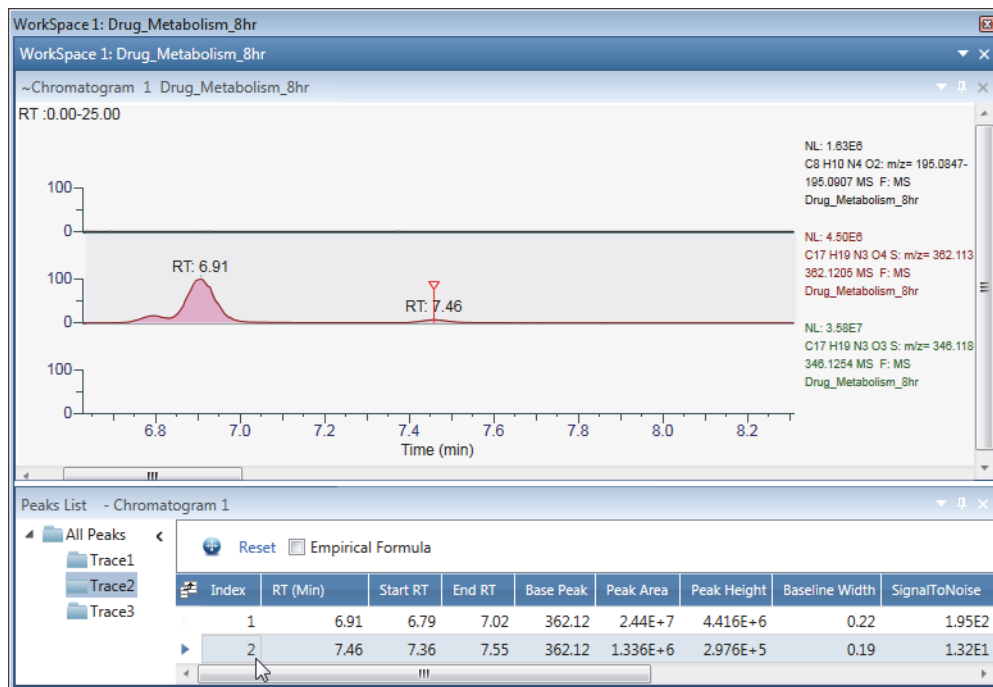
Figure 44. Peak list for Trace 2



2. To check the integration of a peak, select its row in the peak list.

In the Chromatogram view, the application zooms in on the selected peak.

Figure 45 shows the selection of the second peak in the second chromatogram plot.

Figure 45. Zoom level set to the second peak in the second chromatogram plot

❖ **To run the elemental composition algorithm**

Select the **Empirical Formula** check box.

The Empirical Formula column appears to the left of the File Name column.

❖ **To reset the zoom level in the Chromatogram view**

In the Peaks List view, click the **Reset** icon, .

Selecting the Manual Noise Region for the Genesis and ICIS Algorithms

You can manually select the noise region for the Genesis and ICIS peak detection algorithms.

❖ **To specify the manual noise region for the Genesis or ICIS peak detection algorithm**

1. Open the Peak Detection page (see [To display the Peak Detection page](#)).
2. From the peak detection algorithm list, select **ICIS** or **Genesis**, and then specify the appropriate parameter settings.
3. Click the **Workspace Processing** toolbar tab.
4. In the Peak Detection area of the toolbar, click **Select Manual Noise Range**.
5. Drag the pointer horizontally across a region of the chromatographic baseline (the x axis).

3 Reviewing Chromatographic Data

Manually Adding, Undoing, and Deleting Chromatographic Peaks

In the Chromatogram view, the application indicates the noise region with a red line and updates the peak integration. On the Peak Detection page, it automatically selects the Manual Noise Region check box and populates the RT Range box with the selected time range. In the Peaks List view, the application updates the detected peaks list.

6. Do the following:
 - To change the noise region, in the Workspace Processing toolbar, click **Select Manual Noise Range**, and then drag the pointer horizontally across a different region of the chromatographic baseline.
 - To undo the manual noise region, in the Workspace Processing toolbar, click **Clear Manual Noise Range**, and then drag the pointer horizontally across the highlighted noise region.

The application automatically updates the peak detection and integration.

Manually Adding, Undoing, and Deleting Chromatographic Peaks


You can manually add peaks to a chromatogram and the Peaks List view without running a peak detection algorithm. You can also delete peaks from the Peaks List view or undo the last peak that you added manually.

Follow these topics:


- [Manually Adding Chromatographic Peaks](#)
- [Undoing the Last Manually Added Peak](#)
- [Deleting Chromatographic Peaks](#)

Manually Adding Chromatographic Peaks

❖ To manually add peaks to the peak list

1. Click a Chromatogram view to select it.
2. Turn on the Add Peak pointer as follows:
 - a. Open the Chromatogram toolbox.
 - b. Click the **Add Peak** icon, .

Tip You can also turn on the Add Peak pointer by clicking Add Peak in the Chromatogram – Workspace Processing toolbar.

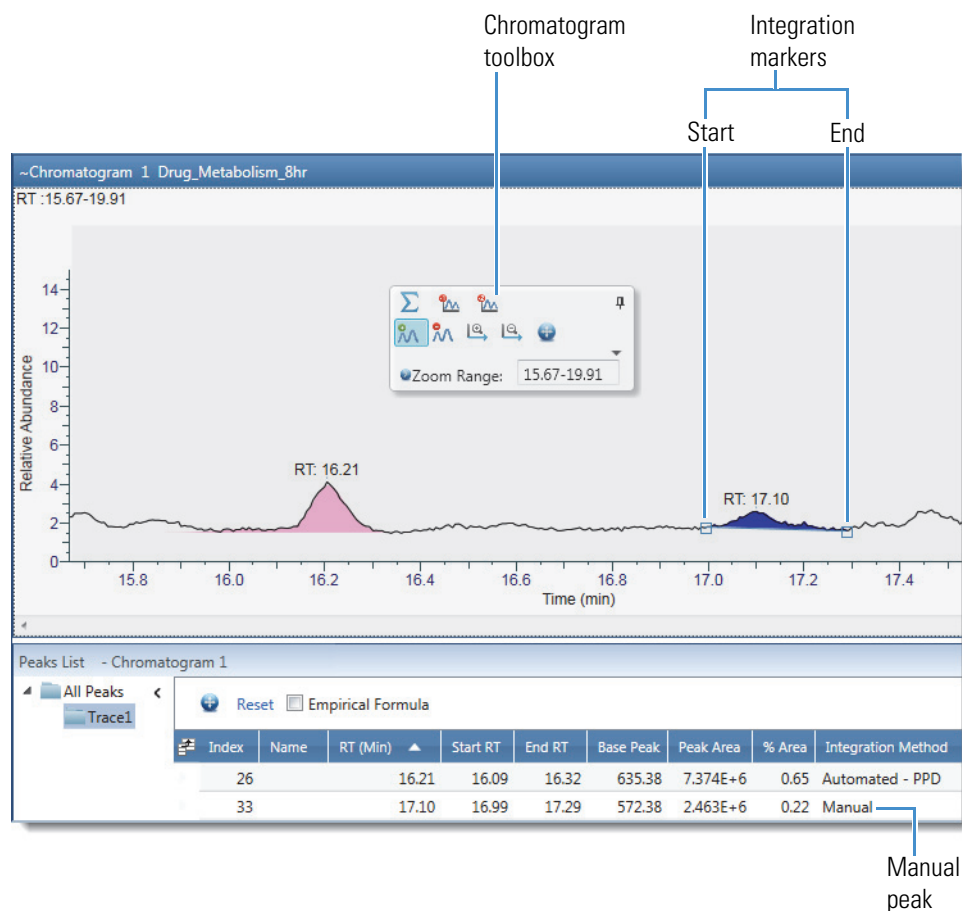
3. Drag the pointer, , along the bases of the peaks that you want to integrate and add to the peak list.

After the peak detection finishes, the application does the following:

- Adds the detected peaks to the peak list in the Peaks List view, and identifies their integration method as Manual.
- Colors the integrated region of the detected peaks blue.
- Draws a baseline under the integration region.
- Adds square integration markers at the beginning and end of the integration region. You can change the integration region by dragging the markers.

Figure 46 shows a manual peak and its integration markers.

Figure 46. Manual peak with integration markers and peaks list with a manual peak



4. To turn off the Add Peak pointer, click the **Add Peak** icon again.

Undoing the Last Manually Added Peak


- ❖ **To undo the last peak that you added manually**

Right-click the chromatogram and choose **Undo Add Peak**.

Deleting Chromatographic Peaks

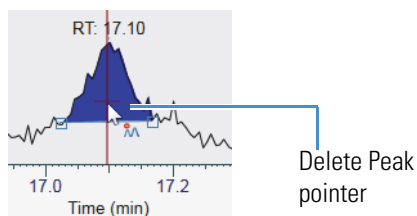
You can delete chromatographic peaks, regardless of their integration method.

❖ To delete chromatographic peaks

1. Click the Chromatogram view to select it.
2. Turn on the Delete Peak pointer as follows:
 - a. Open the Chromatogram toolbox.
 - b. Click the **Delete Peak** icon, .

Tip You can also turn on the Delete Peak pointer by clicking Delete Peak in the [Chromatogram – Workspace Processing Toolbar](#).

3. Click the peak that you want to delete.



4. To turn off the Delete Peak pointer, click the **Delete Peak** icon again.

Importing Components from a Processing Method

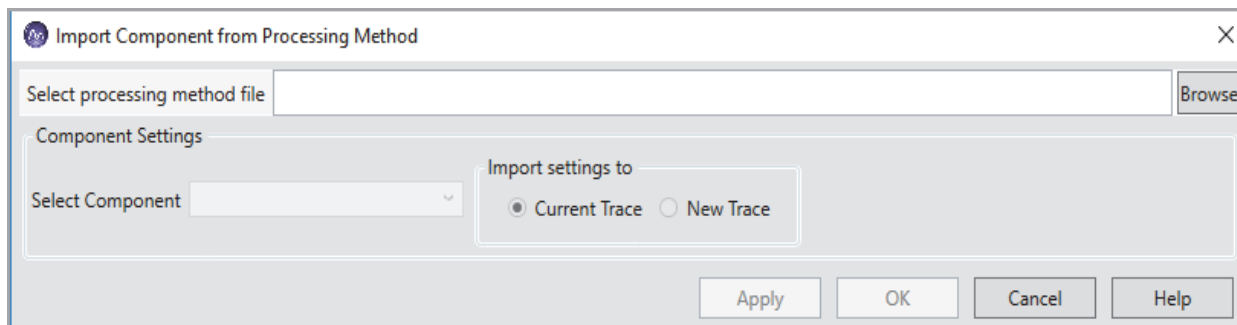
Use the Import Component option to import components from a processing method to the existing workspace. The imported processing method contains component identification parameters that you can apply to the current or new traces in the chromatogram. Processing methods are PMD files.

❖ To import components

1. Click the Chromatogram to make it active.
2. In the Workspace Processing toolbar, click **Import Component**.

The Import Component from Processing Method dialog box opens.

Figure 47. Import Component from Processing Method



3. In the Select processing method file box, click the **Browse** button to locate the .pmd file.
4. In the Open File dialog box, select the .pmd file, and then click **Open**.

The application imports the components and automatically displays the first component.

5. To change the component, click the **Select component** dropdown button, and then select another component.
6. Select the Current Trace or New Trace option to add the component to the current or new trace.
7. Click **Apply**, and then **OK**.

The peak detection is applied to the current or new trace based on the selected option. The component parameters are imported from the processing method to the Peak Detection Info bar page and the Chromatogram Ranges.

Note The Component Name box of the Peak Detection Info page displays the imported component, and you can edit the component name.

Exporting the Scans for a Filtered Chromatogram to a New RAW File

Raw data files can contain gigabytes of data. To work with only the scans related to the displayed time range of a filtered chromatogram, you can export the scans to a new RAW file, which will have a smaller file size. The new RAW file does not include the instrument method or sample information.

❖ To export the scans for a filtered chromatogram to a new RAW file

1. Select the chromatogram trace.
2. In the Workspace Options toolbar, click **Write to RAW**.

In [Figure 48](#), the scan header for the selected trace indicates that it is a filtered trace for the full MS scans (F: MS).

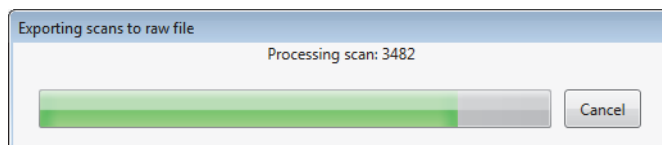
Figure 48. Floating Chromatogram view with two traces



The Export Data dialog box opens to the location of the original RAW file. By default, the application appends _FS to the file name of the original RAW file.

3. Click **Save**.

A progress bar indicates the number of scans written to the new RAW file.



Note If the chromatogram trace is filtered, the application renumbers the scans from 1 to the number written to the new RAW file.

Working with Sequences

A sequence is a list of raw data files where the file names link to the actual RAW files. You can open an existing sequence or create a new sequence from a set of raw data files.

To open an existing sequence, see [Opening Raw Data Files or Sequence Files](#).

Follow these topics:

- [Creating a New Sequence](#)

- [Comparing Chromatogram Traces in a Sequence](#)

Creating a New Sequence

You can create a sequence for any raw data file set.

❖ To create a new sequence

1. Do one of the following:
 - Choose **File > Create Sequence**.-or-
 - In the Workspace Options toolbar, click **Create Sequence**.
2. Browse to and select a set of raw data files.
3. Click **Open**.

The Sequence File page opens in the Info bar. By default, the new sequence list is named NoName.sld.

4. To save the sequence, do the following:
 - a. At the bottom of the Sequence File page, click **Save Sequence**.
 - b. Browse to an appropriate file directory, name the file, and click **Save**.

Comparing Chromatogram Traces in a Sequence

To compare one or more fixed traces against other traces, set up the fixed traces as reference traces in the Chromatogram Ranges view.

❖ To compare chromatogram traces in a sequence

1. Open a sequence (see [Creating, Positioning, Previewing, Refreshing, and Closing Workspaces](#)) or create a sequence (see [Working with Sequences](#)).
2. In the sequence list, select the raw data file that you want to use as a reference file.

The Chromatogram view displays the TIC trace for the raw data file.

3. Set up the reference trace as follows:
 - a. In the Workspace Options toolbar, click **Chromatogram Ranges** to open the Chromatogram Ranges view.
 - b. Define the trace that you want to use as a reference trace. For example, select the detector type, trace type, filter, and so on.
 - c. Copy this definition to the last row in the Chromatogram Ranges view by selecting its corresponding check box in the Display column.

3 Reviewing Chromatographic Data

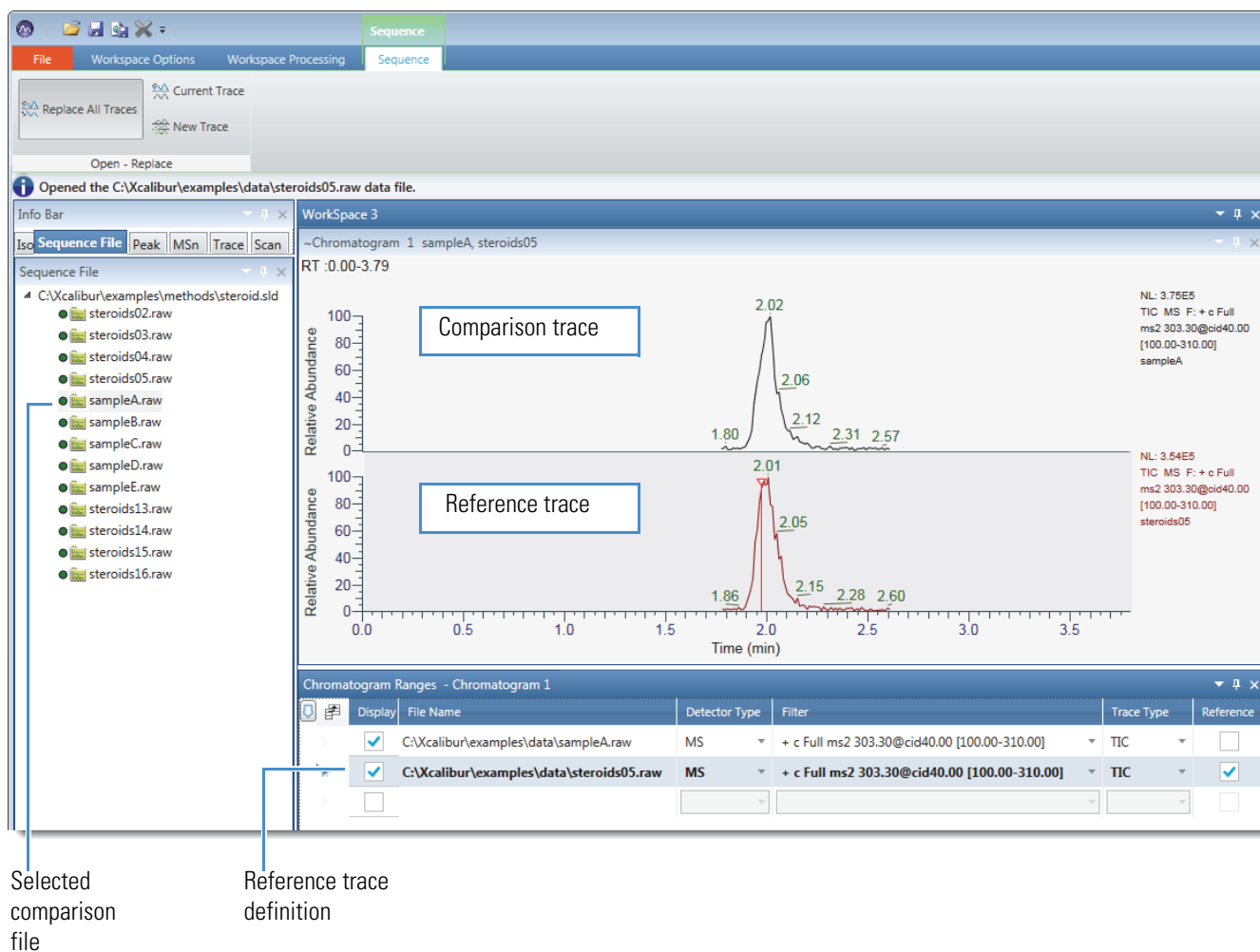
Chromatogram-Specific Toolbars

A copy of the first chromatogram trace appears below it.

- d. Select the check box in the Reference column of the row corresponding to the reference trace.
4. Click the raw data files in the sequence to compare their chromatogram traces to the reference chromatogram trace.

Figure 49 shows a Chromatogram view with two plots. The comparison trace for the selected file is on the top, and the reference trace is on the bottom.

Figure 49. Chromatogram view with a comparison trace and a reference trace



Chromatogram-Specific Toolbars

Use the following toolbars to format the Chromatogram view, detect chromatographic peaks, and work with sequence lists.

- [Chromatogram – Display Options Toolbar](#)

- Chromatogram – Workspace Processing Toolbar
- Sequence Toolbar

Chromatogram – Display Options Toolbar

Use the Chromatogram Display Options toolbar buttons, its Skew and Elevation sliders, and its Label Threshold box to customize the Chromatogram view.

❖ To display the Chromatogram – Display Options toolbar

1. Click a Chromatogram view to select it.
2. Click the **Display Options** toolbar tab.

Figure 50 shows the Chromatogram – Display Options toolbar, and Table 13 describes the toolbar buttons.

Figure 50. Chromatogram Display Options toolbar (in two parts, left and right side)

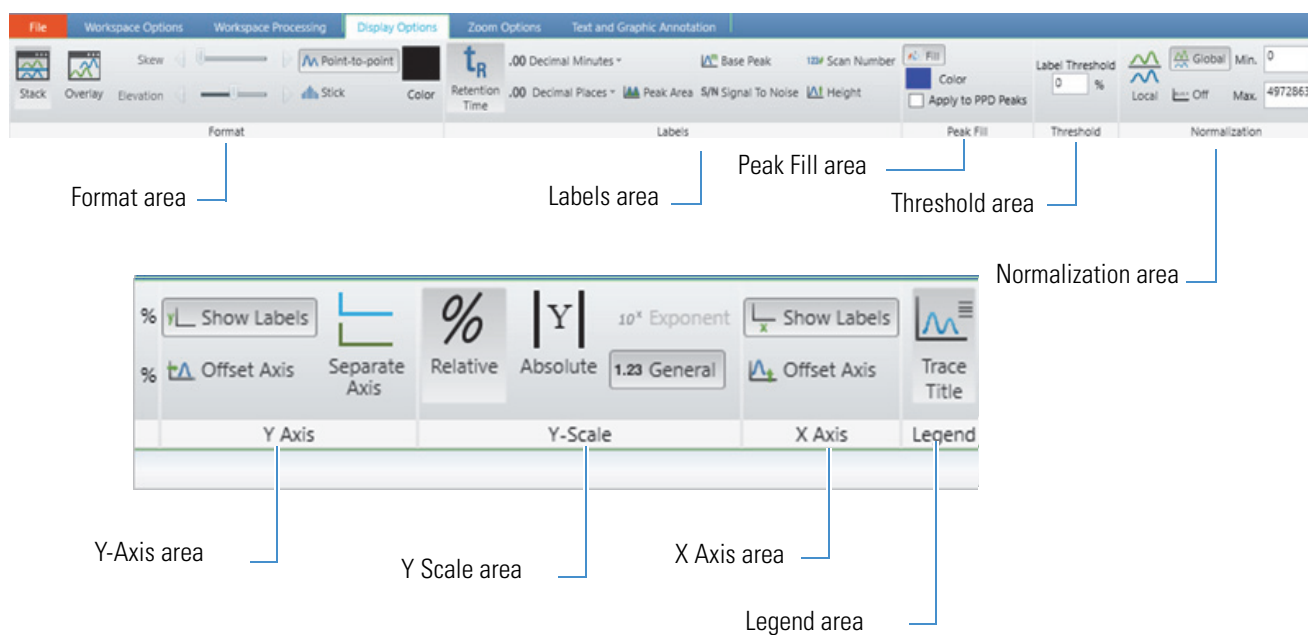


Table 13. Chromatogram – Display Options toolbar buttons (Sheet 1 of 5)

Button	Description
Format	
Use the parameters in the Format area to modify the appearance of the Chromatogram view.	
Stack	Stacks the chromatogram traces vertically.
Overlay	Overlays the chromatogram traces vertically with optional horizontal skew (time offset).

Table 13. Chromatogram – Display Options toolbar buttons (Sheet 2 of 5)

Button	Description
Skew	<p>Sets the skew angle (time offset) to a value from 0–45 degrees for an overlay arrangement of chromatogram traces.</p> <p>To set the skew, drag the Skew slider.</p>
Elevation	<p>Sets the vertical spacing for an overlay arrangement of chromatogram traces.</p> <p>To set the vertical spacing, drag the Elevation slider. Move the Elevation slider to the farthest left to overlay the plots on top of each other.</p>
Point-to-point	Connects the signal data points to form a continuous curve.
Stick	Displays the signal data points by using vertical lines.
Color	Displays the color palette, where you select the colors of the chromatogram traces.
Labels	
Use the buttons in the Labels area to annotate chromatogram peaks.	
Retention Time	<p>Adds a retention time label (in minutes) above each local maxima or chromatogram peak.</p> <p>The order of chromatogram labels for an undetected peak, from top to bottom, is retention time, base peak, and scan number. The Chromatogram view displays the retention time on all peaks that meet the selection criteria.</p> <p>The letters <i>RT</i> to the left of the value indicate a retention time.</p>
Retention time format	<p>Specifies the units and format for the retention time label.</p> <p>Selections: Decimal Minutes, Decimal Minutes with Units, Minutes and Seconds, Seconds, Seconds with Units</p>
Decimal Places	<p>Changes the number of decimal places in the retention time label.</p> <p>Range: 0–5</p>
Peak Area (of integrated peaks)	<p>Adds a peak area label above each detected chromatogram peak.</p> <p>The letters <i>MA</i> or <i>AA</i> next to the value indicate manual integration or automatic integration, respectively.</p>
Base Peak	<p>Adds a base peak mass-to-charge ratio label (in <i>m/z</i>) above each local maxima or each detected chromatogram peak.</p> <p>The letters <i>BP</i> next to the value indicate a base peak.</p>

Table 13. Chromatogram – Display Options toolbar buttons (Sheet 3 of 5)

Button	Description
Signal To Noise (of integrated peaks)	<p>Adds a signal-to-noise ratio label above each detected chromatogram peak.</p> <p>The Avalon peak detection algorithm does not calculate a signal-to-noise ratio. For the Genesis and ICIS peak detection algorithms, you can choose to report the root mean square (RMS) noise values. The application calculates the RMS signal-to-noise ratio for manual peaks.</p> <p>The letters <i>SN</i> next to the value indicate that it is a signal-to-noise ratio.</p>
Scan Number	<p>Adds the active scan number label above the chromatogram peak.</p> <p>The letter <i>S#</i> next to the value indicates the scan number.</p>
Height (of integrated peaks)	<p>Adds a peak height above each detected chromatogram peak.</p> <p>The letters <i>MH</i> or <i>AH</i> next to the value indicate manual integration or automatic integration, respectively.</p>
Label Threshold	<p>Sets the percentage of highest data point in the trace to 100%.</p> <p>The application labels only the local maxima or chromatographic peak apexes that are above the specified height threshold.</p> <p>Range: 0–100; default: 0%</p>
Peak Fill	
<p>Use the fill options in the Peak Fill area to color the chromatogram traces. See Filling the Chromatogram Peaks.</p>	
Fill	Turns on or off peak fill in the active chromatogram view.
Color	Opens the color palette, where you can select the colors to fill the integrated peaks.
Apply to PPD peaks	Fills the PPD peaks in the active chromatogram view with the selected color from the color palette.
Normalization	
<p>Use the parameters in the Normalization area to specify how the application normalizes the chromatogram traces. See Normalizing Chromatogram Traces.</p>	
Local	Normalizes each chromatogram trace with respect to the intensity of the most intense peak in that trace.
Global	Normalizes the chromatogram traces so that the most intense peak of all chromatograms is 100 percent.

Table 13. Chromatogram – Display Options toolbar buttons (Sheet 4 of 5)

Button	Description
Off	Scales all chromatograms by using the maximum and minimum absolute values. Note If you set the normalization to Off, the <i>y</i> -axis scale to Absolute, and change the trace type, the new trace might not display. To remedy this, set the normalization to Local and then to Off.
Min.	Displays the minimum of the <i>y</i> axis. Enter a value in the box to change the minimum. The application indicates whether the value is a percentage or an intensity.
Max.	Displays the maximum of the <i>y</i> axis. Enter a value in the box to change the maximum value. The application indicates whether the value is a percentage or an intensity.
Y Axis	
Use the parameters in the Y Axis area to specify how the application labels and displays the <i>y</i> axis. See Changing the Y-Axis Title and Scale of the Chromatogram View .	
Show Labels	Shows or hides the <i>y</i> -axis label.
Offset Axis	Sets the location of the displayed plot at a specified distance from the <i>y</i> axis. The <i>y</i> -axis offset moves the <i>x</i> axis slightly to the right of the <i>y</i> axis so that you can see the plot details at low <i>x</i> -axis values.
Separate Axis	Sets the Separate Axis title for the active Chromatogram trace. Note Enable the Separate Axis button to view the text box. Use this box to change the <i>y</i> axis title.
Y-Scale	
Use the parameters in the Y-Scale area to specify how the application scales the <i>y</i> axis. See Changing the Y-Axis Title and Scale of the Chromatogram View .	
Y Absolute	Scales the chromatograms traces by using the absolute values. Note If you set the normalization to Off, the <i>y</i> axis to Absolute, and change the trace type, the new trace might not display. To remedy this, set the normalization to Local and then to Off.
10 ^x Exponent	Sets the Y-Scale unit in exponent form.
1.23 General	Sets the Y-Scale unit in general form.
% Relative	Scales the chromatograms traces by using percentages.

Table 13. Chromatogram – Display Options toolbar buttons (Sheet 5 of 5)

Button	Description
X Axis	
Use the parameters in the X Axis area to specify how the application labels and displays the <i>x</i> axis.	
Show Labels	Shows or hides the <i>x</i> -axis label.
Offset Axis	Sets the location of the displayed plot at a specified distance from the <i>x</i> axis. The <i>x</i> -axis offset moves the <i>y</i> axis up slightly so that you can see the plot details at low <i>y</i> -axis values.
Legend	
Trace Title	Moves the title of the chromatogram trace from above the trace to beside it.

Chromatogram – Workspace Processing Toolbar

Use the buttons in the Peak Detection area of the Workspace Processing toolbar to detect peaks in chromatogram traces.

❖ To display the Chromatogram Workspace Processing toolbar

1. Click a Chromatogram view to select it.
2. Click the **Workspace Processing** toolbar tab.

Figure 51 shows the Peak Detection area of the Workspace Processing toolbar. The Delete Peak button becomes available after you apply a peak detection algorithm or add a peak manually.

Figure 51. Peak Detection area of the Workspace Processing toolbar

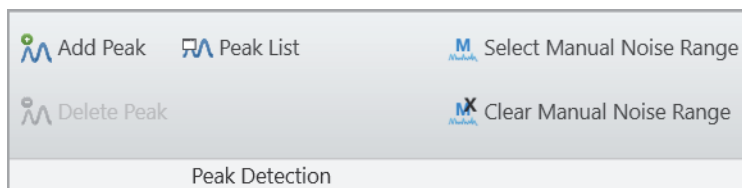


Table 14 describes the Peak Detection area buttons.

Table 14. Workspace Processing toolbar – Peak Detection area buttons

Button	Description
Detect in Active Plot	<p>Runs the Genesis, ICIS, or Avalon peak detection algorithm on the active chromatogram plot.</p> <p>Click a chromatogram plot to make it active. The Chromatogram view shades the active plot area gray. See Automatically Detecting and Integrating Chromatographic Peaks.</p>
Detect in All Plots	<p>Runs the Genesis, ICIS, or Avalon peak detection algorithm on all chromatogram plots. See Automatically Detecting and Integrating Chromatographic Peaks.</p>
Add Peak	<p>Adds a peak to the peak list. See Manually Adding Chromatographic Peaks.</p>
Delete Peak	<p>Deletes a peak from the peak list. See Deleting Chromatographic Peaks.</p> <p>Available after you apply a peak detection algorithm or manually add a chromatographic peak.</p>
Peak List	<p>Displays the Peaks List view. When you run a peak detection algorithm, the Peak List view lists the peaks that the application found in the chromatogram. You can also manually add peaks to the peak list with the Add Peak button.</p>
Select Manual Noise Range	<p>Freezes the Chromatogram view (temporarily turns off zooming with the mouse pointer) so that you can use the mouse pointer to define a noise region (see Selecting the Manual Noise Region for the Genesis and ICIS Algorithms).</p>
Clear Manual Noise Range	<p>Freezes the Chromatogram view (temporarily turns off zooming with the mouse pointer) so that you can use the mouse pointer to clear the selection of a manual noise region (see Selecting the Manual Noise Region for the Genesis and ICIS Algorithms).</p>

Sequence Toolbar

Use the Sequence toolbar buttons to determine which traces appear in the Chromatogram view as you click through the raw data files in a sequence.

❖ To display the Sequence toolbar

1. Open a sequence (see [Creating, Positioning, Previewing, Refreshing, and Closing Workspaces](#)) or create one (see [Working with Sequences](#)).

- If the Sequence File page is not open, click the **Sequence File** tab in the Info Bar.

Tip The Sequence toolbar is visible only when the Sequence File page is open.

Figure 52 shows the Sequence toolbar, and Table 15 describes the toolbar buttons.

Figure 52. Sequence toolbar

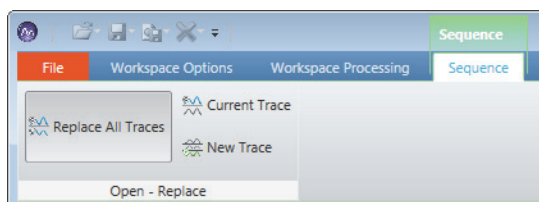


Table 15. Sequence toolbar buttons

Button	Description
Replace All Traces	<p>Replaces all the raw data files in the chromatogram ranges list—and their chromatogram traces in the Chromatogram view—with the raw data file that you select in the sequence.</p> <p>Select Replace All Traces, and then select a raw data file in the sequence on the Sequence File page for replacement.</p> <p>Note The application does not replace a trace that is set as a reference trace (its Reference check box is selected) with the selection from the sequence.</p>
Current Trace	<p>Replaces only the raw data file that you select in the chromatogram ranges list—and its chromatogram trace displayed in the Chromatogram view—with the raw data file that you select in the sequence.</p> <p>Select a trace row in the chromatogram ranges list, select Current Traces, and then select a raw data file in the sequence for replacement.</p> <p>Note The application does not replace a trace that is set as a reference trace (the box in the Reference column is selected) with the selection from the sequence.</p>
New Trace	<p>Adds new raw data files to the chromatogram ranges list and displays their chromatogram traces in the Chromatogram view.</p> <p>Select New Trace, and then select one or more raw data files in the sequence to add to the list.</p>

Chromatogram-Specific Views

For information about the chromatogram-related views, see these topics:

3 Reviewing Chromatographic Data

Chromatogram-Specific Views

- [Chromatogram View](#)
- [Chromatogram Ranges View](#)

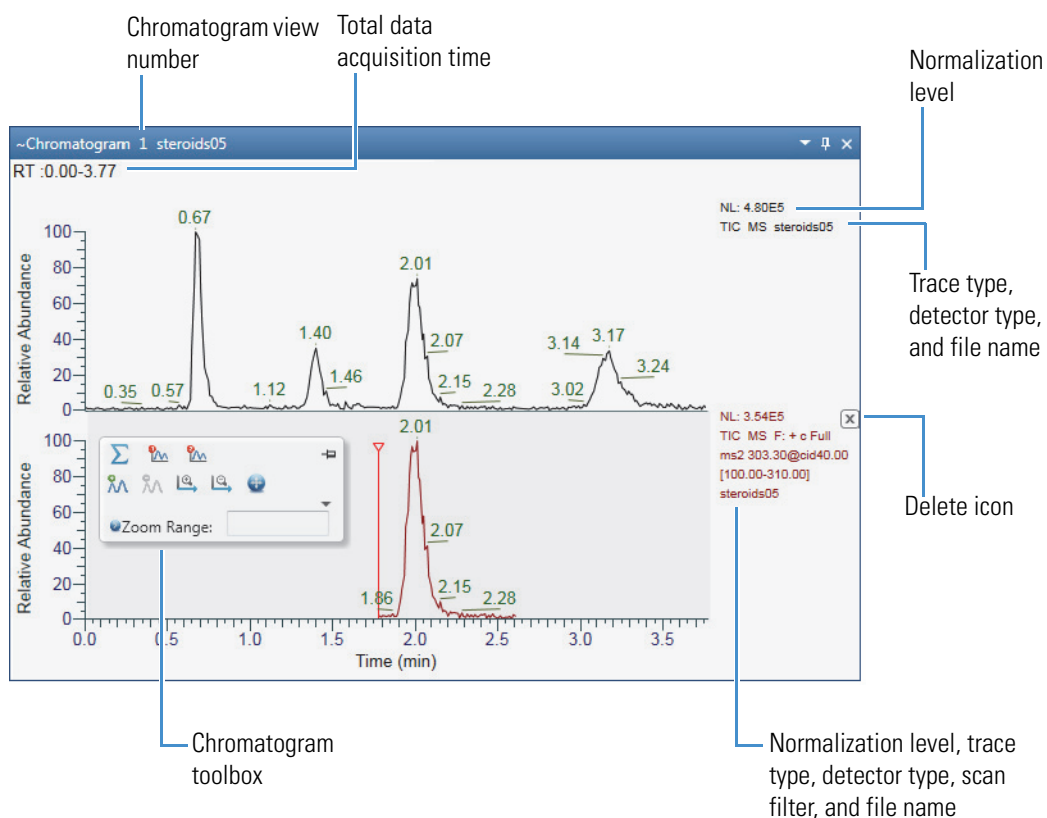
Chromatogram View

Use the Chromatogram view to display chromatogram traces. You can delete a Chromatogram view, add multiple chromatogram traces to each Chromatogram view, and add multiple Chromatogram views to the Workspace.

For information about displaying and reviewing the chromatographic data, see the list of topics in the beginning of this chapter.

[Figure 53](#) shows an example of a Chromatogram view with two traces. For information about the scan header that appears at the top right of each chromatogram trace, see [Scan Headers and Scan Header Abbreviations](#). For information about the view's toolbox, see [Chromatogram Toolbox](#).

Figure 53. Chromatogram view with two traces and the chromatogram toolbox



Note You can set a minimum trace height value, in centimeters, in the Default Workspace Options page. When you adjust the height of a Chromatogram view, if the height of the traces becomes smaller than the set minimum value, a scrollbar automatically appears to the right of the view.

Right-clicking a Chromatogram view displays a shortcut menu (Figure 54) with the commands described in Table 16.

Figure 54. Shortcut menu for a Chromatogram view

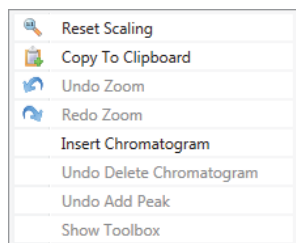


Table 16. Chromatogram view shortcut menu commands

Command	Description
Reset Scaling	Resets the scaling of all the plots in the Chromatogram view.
Copy To Clipboard	Copies an image of the Chromatogram view to the Clipboard.
Undo Zoom	Undoes the last zoom.
Redo Zoom	Reapplies the last zoom that you undid.
Insert Chromatogram	Inserts a copy of the selected chromatogram trace.
Undo Delete Chromatogram	Displays the most recently deleted chromatogram trace.
Undo Add Peak	Removes the most recently added manual peak from the Chromatogram view and the Peaks List view.
Show Toolbox	Displays the Chromatogram toolbox. To close the toolbox, click its pin icon.

These toolbars are available when a Chromatogram view is selected:

- [Chromatogram – Workspace Processing Toolbar](#)
- [Chromatogram – Display Options Toolbar](#)
- [Zoom Options Toolbar](#)
- [Text and Graphic Annotation Toolbar](#)

Related Topics

- [Reviewing Chromatographic Data](#)

Chromatogram Ranges View

Use the Chromatogram Ranges view (Figure 55) to select which chromatogram traces to display in a Chromatogram view. For instructions, see [Defining a Chromatogram Trace from the Chromatogram Ranges View](#).

Note For information about exporting the tabular data in the Chromatogram Ranges view to a CSV file, see [Exporting or Printing the Contents of a View or Workspace](#). You can paste the contents of a CSV file to the Chromatogram Ranges view if the table columns match.

❖ To open the Chromatogram Ranges view

In the Workspace Options toolbar, click **Chromatogram Ranges**.

To add, delete, and move rows in the Chromatogram Ranges view, see these topics:

- [Using the Fill Down Feature](#)
- [Moving a Row](#)
- [Deleting Rows](#)

Figure 55 shows an example of the Chromatogram Ranges view.

Figure 55. Chromatogram Ranges view showing the entry of a chemical formula

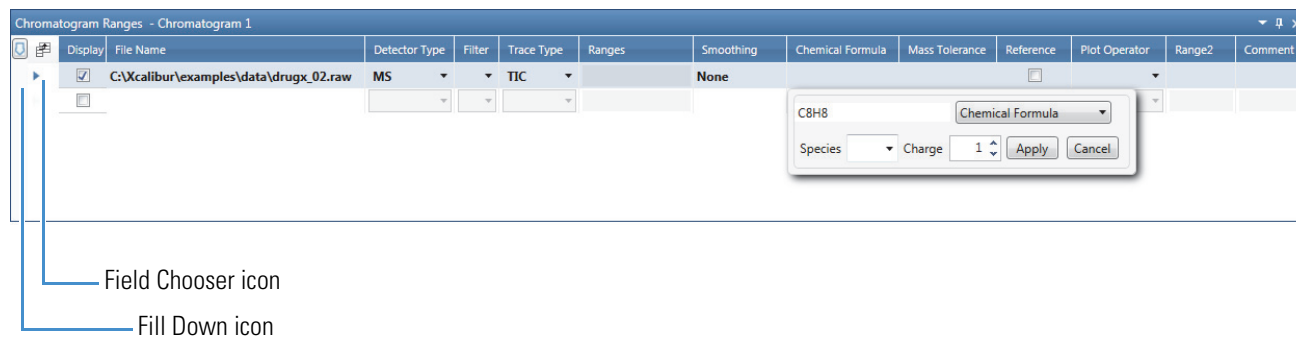


Table 17 describes the parameters and icons for the Chromatogram Ranges view. The following columns have corresponding pages in the Info Bar: Detector Type, Trace Type, and Filter.

Table 17. Chromatogram Ranges view parameters and icons (Sheet 1 of 4)


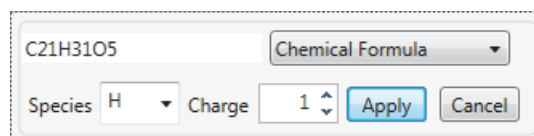
Parameter/Icon	Description
Display	Select this check box to display the chromatogram trace in a Chromatogram view. Clear this check box to remove the trace.
File Name	Displays the path and name of the raw data file. To display the Open Raw File dialog box, click  at the end of the file name.
Detector Type	Select the detector type. The application autosenses the available detector types based on the information in the raw data file. For details, see Detector Type Page .
Filter	Select or type a filter to apply to the raw data file. The application autosenses the metafilters (including SRM or compound filters) in the raw data file in this order: MS, MS2, ETD, HCD, and then the individual scan filter list. For data files with more than one MS/MS order, enter MSn to include all the MS/MS data. For grouped filters, the Filter dropdown list displays only the first filter in the group. The Spectrum view displays the actual filter for the scan. For more information, see Scan Filter Parameters and Scan Filters Page .
Trace Type	Select the trace type as follows: <ul style="list-style-type: none"> • For MS, select TIC, Base Peak, Mass Range, Neutral Fragment, or Mass Defect Filter. • For PDA, select Total Scan, Wavelength Range, or Maximum Spectrum. • For MS Trending, select one of the instrument status parameters, such as the API source voltage.
Ranges	Specify the <i>m/z</i> ranges of the displayed spectrum, such as 443.2, 534.6, 600–800. The application autosenses the default range from the raw data file. If the field is inactive, it is not applicable based on the trace type.

Table 17. Chromatogram Ranges view parameters and icons (Sheet 2 of 4)

Parameter/Icon	Description
Smoothing	Select which type of smoothing, if any, to apply to the chromatogram trace. <ul style="list-style-type: none">• Gaussian: Uses weighting coefficients corresponding to a Gaussian peak shape to average each data point with neighboring points to give the displayed value.• Moving mean: Uses equal weighting to average each data point with neighboring points to give the displayed value.
Chemical Formula	Clicking this table cell opens the following dialog box.



Select **Chemical Formula** or **Peptide**:

- For a compound, select **Chemical Formula**, use the symbols in the periodic table to specify the elements in the chemical formula, select the ionization species, and type or select the charge of the adduct ion of interest.

Example: For calcium copper titanate (which is made up of one calcium atom, three copper atoms, four titanium atoms, and twelve oxygen atoms), enter the chemical formula: CaCu₃Ti₄O₁₂.

- For a peptide, select **Peptide**, enter the peptide sequence using the one-letter symbols for the amino acid residues (see [One- and Three-Letter Abbreviations for Amino Acid Residues](#)), select the ionization species, and type or select the charge of the adduct ion of interest.




Example: For Hexarelin (a hexapeptide), use the one-letter code—HWAWFK.

When you click Apply, the Trace Type changes to Mass Range, and the mass-to-charge ratio of the specified ion appears in the Ranges column.

Table 17. Chromatogram Ranges view parameters and icons (Sheet 3 of 4)

Parameter/Icon	Description
Mass Tolerance	<p>Entering a chemical formula in the Chemical Formula column or selecting Mass Range in the Trace Type column enables this table cell.</p> <p>Specify the mass tolerance for the mass-to-charge ratio of the specified ion.</p> <p>Range: 0.00 to 10.00; default: 0.5 amu</p>
Delay Time (hidden by default)	<p>Specify a delay time (in minutes) to shift trace 1 (Trace Type column) so that it aligns with trace 2 (Trace Type2 column). A positive delay time shifts trace 1 to higher retention times by this amount.</p> <p>Range: -5.00 to 5.00; default: 0.00</p>
Reference	<p>Select this check box to make the chromatogram trace a reference trace. A reference trace is always displayed in the Chromatogram view so that you can compare other traces to it.</p>
Plot Operator	<p>Select a plot operator, either + (add) or - (subtract).</p> <ul style="list-style-type: none"> • If you select +, the Chromatogram view displays trace 1 plus trace 2. • If you select -, the Chromatogram view displays trace 1 minus trace 2. <p>For example, from a chromatogram you can use the subtract (-) plot operator to subtract contributions from a solvent or from other noise.</p> <p>For MS detectors, these are the trace 1 and trace 2 combinations:</p> <ul style="list-style-type: none"> • mass range ± mass range • mass range ± base peak • base peak ± mass range • base peak ± base peak • TIC - mass range • TIC - base peak


Table 17. Chromatogram Ranges view parameters and icons (Sheet 4 of 4)

Parameter/Icon	Description
	<p>For PDA detectors, these are the trace 1 and trace 2 combinations:</p> <ul style="list-style-type: none"> wavelength range ± wavelength range wavelength range ± spectrum maximum spectrum maximum ± wavelength range spectrum maximum ± spectrum maximum total scan – wavelength range total scan – spectrum maximum
Trace Type2 (hidden by default)	Select the type of trace to add to or subtract from trace 1.
Range2	Specify the wavelength or mass range for trace 2.
Comment	Specify a comment for the row.
Icon	
 Fill Down	<p>Opens the Fill Down dialog box for populating multiple rows with duplicate parameter settings.</p> <p>For details, see Using the Fill Down Feature.</p>
 Field Chooser	<p>Displays the Field Chooser dialog box for selecting which columns appear in the Chromatogram Ranges view.</p> <p>For details, see Selecting the Columns to Display in a View or Dialog Box with Tabular Data.</p>
Check boxes in the Display column	<p>Select the check box to add the defined chromatogram to the Chromatogram view. Selecting the check box in the last row populates the row with the settings from the previous row.</p> <p>You can edit the File Name parameter to change the new entry to that raw data file, or click  at the end of the file name to browse to and select a new raw data file.</p>

Using the Fill Down Feature

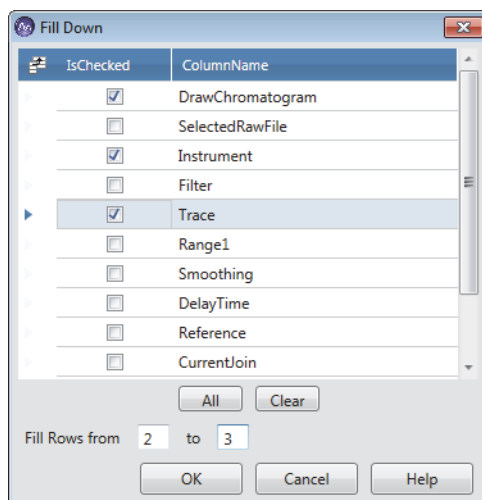
To set up several similar chromatogram traces, use the Fill Down feature in the Chromatogram Ranges view. With this feature, you can copy the selected parameter values from the currently selected row to other specified rows.

❖ **To fill down data**

1. In the Chromatogram Ranges view, select the trace row that you want to copy data from.
2. Click the **Fill Down** icon ()

The Fill Down dialog box opens ([Figure 56](#)).

Figure 56. Fill Down dialog box



3. Do one of the following:
 - Select the check box for each parameter value that you want to fill down.
 - Click **All** to select all the check boxes.

–or–

- Click **Clear** to clear all the check boxes.
4. In the Fill Rows boxes, type the starting and ending row numbers, and then click **OK**.

The application copies the selected parameter values from the currently selected row and pastes them into the specified starting and ending rows. If a value is invalid, it appears in red.

Moving a Row

You can move a row in the Chromatogram Ranges view by dragging it to the new location.

❖ **To move a row**

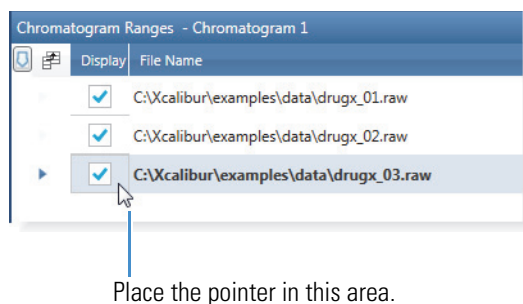
1. Place the pointer in an area somewhere in the row that you want to move, but not inside an editable field and not in the blank area to the left of the row.

For example, place the pointer in the area next to the Display check box ([Figure 57](#)).

3 Reviewing Chromatographic Data

Chromatogram-Specific Views

Figure 57. Example pointer location to move a row



2. Drag the row up or down to move it to the new location.


As you drag the row above or below another row, that other row temporarily turns red. After you drop the row, it appears above or below the red row.

Deleting Rows

You can delete rows in the Chromatogram Ranges view. For each row that you delete, the application automatically deletes the corresponding trace in the Chromatogram view.

Note When the view has only one row, you cannot delete the row.

❖ To delete rows in the Chromatogram Ranges view

1. Select the rows that you want to delete as follows:
 - To select a single row, click the arrow icon, , that appears when you point to the left of the Display column.
 - To select adjacent rows, drag the pointer across the group of rows (in the column to the left of Display column) or use the SHIFT key.
 - To select nonadjacent rows, press the CTRL key and click to the left of the Display column for each row.

Selected rows are highlighted in blue.

2. Press the DELETE key and click **Yes** in the confirmation dialog box to remove the selected rows.

Peaks List View

The Peaks List view lists the chromatogram peaks that the peak detection algorithm automatically finds and also lists the peaks that you manually add.

❖ To display the Peaks List view

- Apply the ICIS, Genesis, Avalon, or PPD peak detection algorithm (see [Automatically Detecting and Integrating Chromatographic Peaks](#)).

IMPORTANT Applying the peak detection algorithm the first time automatically opens the Peaks List view. However, if you then close this view and apply the peak detection algorithm again, the Peaks List view does not automatically reopen. To manually open this view, click the **Workspace Processing** toolbar tab, and then click **Peak List**.

- Manually add a peak (see [Manually Adding, Undoing, and Deleting Chromatographic Peaks](#)).

—or—

- Click the **Workspace Processing** toolbar tab.
- Click **Peak List**.

Figure 58 and Figure 59 show examples of the Peaks List view. In Figure 59, the Name column is populated with the compound name specified in the instrument method for the selected scan filter. The table values are read-only.

Figure 58. Peaks List view

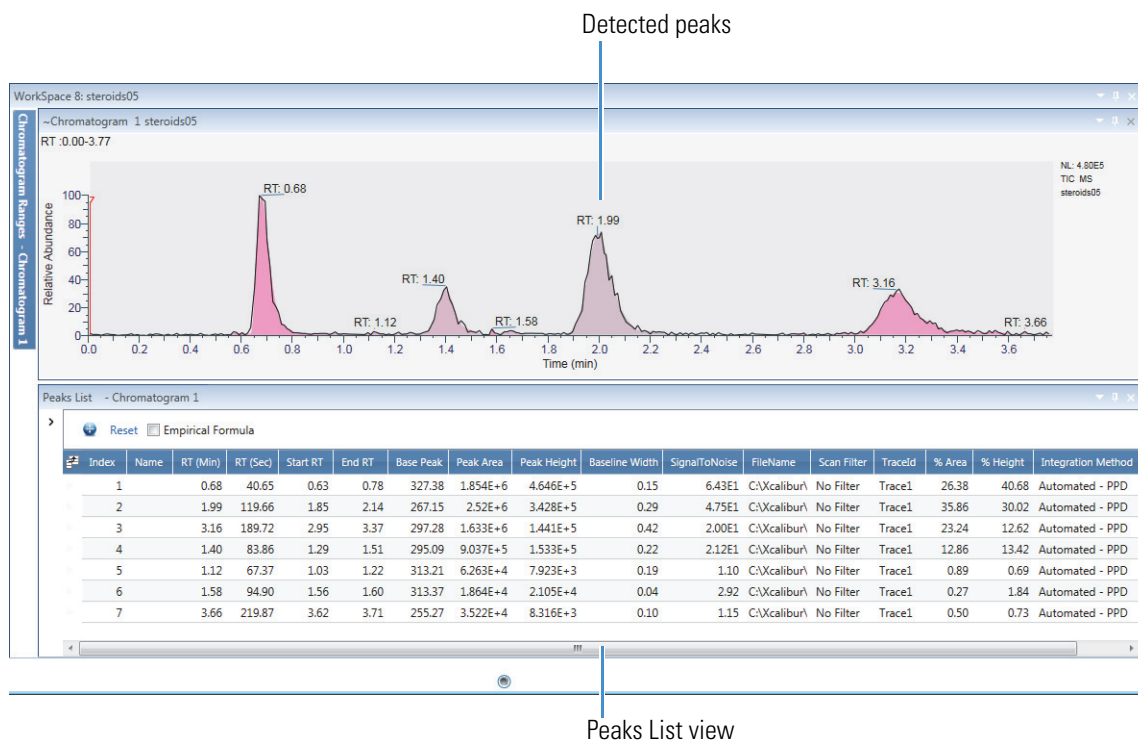


Figure 59. Peak List view with a populated Name column

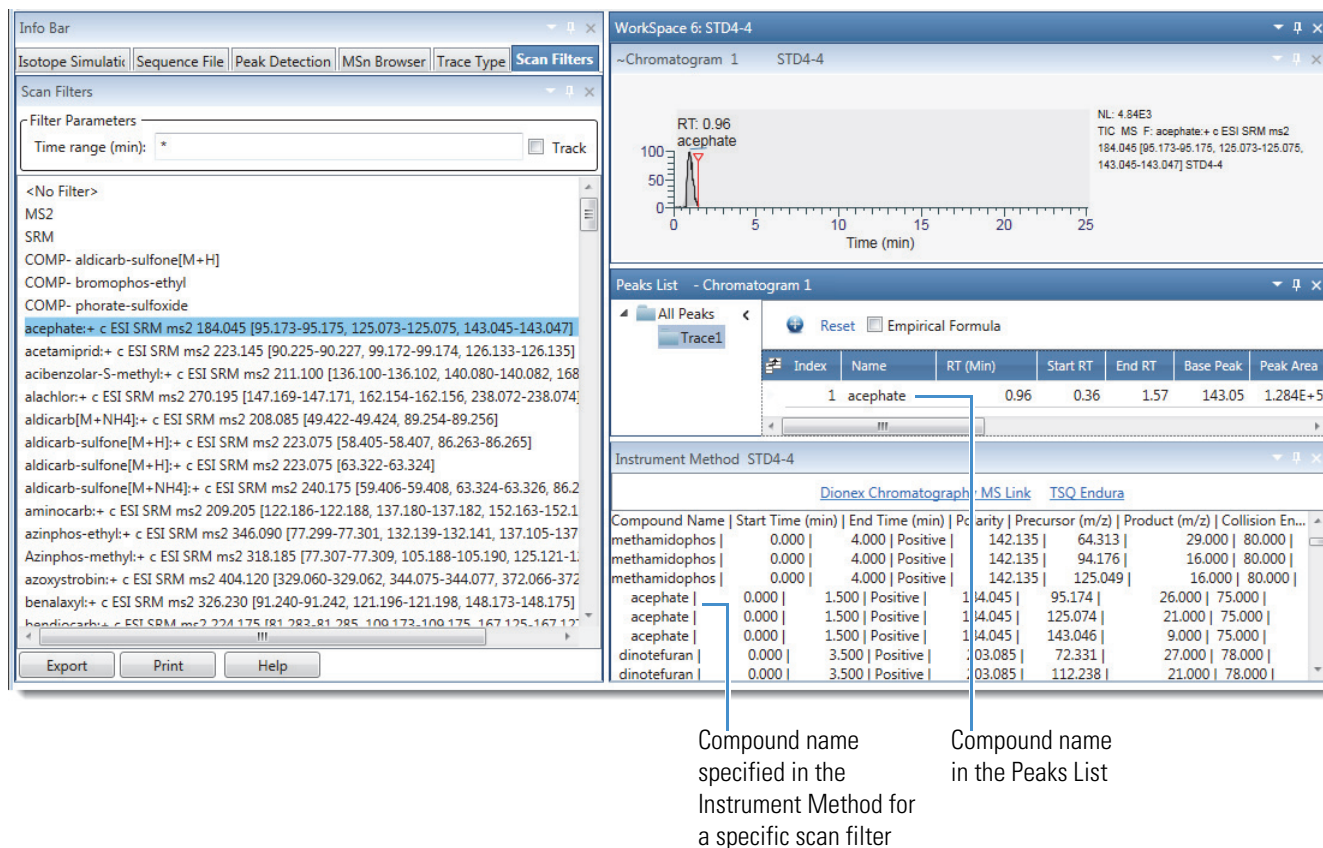


Table 18 describes the parameters and icons for the Peaks List view.

Table 18. Peaks List view parameters and icons (Sheet 1 of 3)



Parameter	Description
	Restores the data display in the Chromatogram view to the full range of the x axis and y axis after you zoom in by clicking a row in the peaks list.
	Displays the Field Chooser dialog box, where you select which fields appear in the peaks list (see Selecting the Columns to Display in a View or Dialog Box with Tabular Data).
Elemental Composition	Select to run the elemental composition algorithm. This algorithm requires full-scan data.
Field	
Index	Unique identification number for each chromatogram peak. The index increments from lowest to highest retention time in a trace and from the top trace to the bottom trace.

Table 18. Peaks List view parameters and icons (Sheet 2 of 3)


Parameter	Description
Name	Compound name from the instrument method (Figure 59). Note For some Thermo Scientific mass spectrometers, for example, the triple quadrupole product line, the instrument method that you use for data acquisition includes a table where you can associate a compound name with each scan filter. When you view the scan filters on the Scan Filters page, the associated compound names appear next to the scan filter. You cannot edit the names in the FreeStyle application.
RT (Min)	Retention time (in minutes) corresponding to the apex of the peak.
RT (Sec)	Retention time (in seconds) corresponding to the apex of the peak.
Start RT	Retention time corresponding to the start of the chromatographic peak, where the detection signal increases beyond the threshold criteria.
End RT	Retention time corresponding to the end of the chromatographic peak, where the detection signal decreases below the threshold criteria.
Base Peak	Mass-to-charge ratio of the most abundant ion at the apex of the peak. The Peak List view displays 0.00 if no mass spectral data is present.
Peak Area	Area of the peak in units of counts × seconds.
Peak Height	Number of counts at the peak apex.
Baseline Width	Difference (in minutes) between the start and end retention times.
Signal To Noise	Signal-to-noise ratio of the integrated peak.
Empirical Formula	Chemical formula determined by the elemental composition algorithm. By default, this column is hidden. Selecting the Empirical Formula check box adds the column to the table and runs the elemental composition algorithm.
File Name	Displays the path and name of the raw data file. To display the Open Raw File dialog box, click the browse icon,  , at the end of the file name.


Table 18. Peaks List view parameters and icons (Sheet 3 of 3)

Parameter	Description
Scan Filter	Select a filter to apply to the raw data file. The application autosenses the metafilters (including SRM or compound filters) in the raw data file in this order: MS, MS2, ETD, HCD, MSN, and then the individual scan filter list. For grouped filters, the Filter list displays only the first filter in the group. The Spectrum view displays the actual filter for the scan. For more information, see Scan Filter Parameters .
Trace ID	Number that identifies the chromatogram trace. The Trace ID increments from the top trace to the bottom trace.
% Area	Peak area as a percentage of the total peak area.
% Height	Peak height as a percentage of the total peak height.
Integration Method	Displays the integration method: ICIS, Genesis, Avalon, PPD, or manual.


Chromatogram Toolbox

Use the Chromatogram toolbox to create average and background subtracted spectra, add manual peaks, and zoom in or out of a Chromatogram view.

How you open the toolbox depends on your workspace setting (see [Default Workspace Options Page](#)). With the factory default setting, you can open the toolbox in two ways:

- Pause the pointer in the view. The toolbox opens with its pin icon in the unpinned (horizontal) position, . To keep the toolbox open, click its pin icon.

—or—

- Right-click the view and choose Show Toolbox. The toolbox opens with its pin icon in the pinned (vertical) position, . To close the toolbox, click its pin icon.

[Figure 60](#) shows the floating toolbox, and [Table 19](#) describes its icons and fields.

Figure 60. Chromatogram toolbox

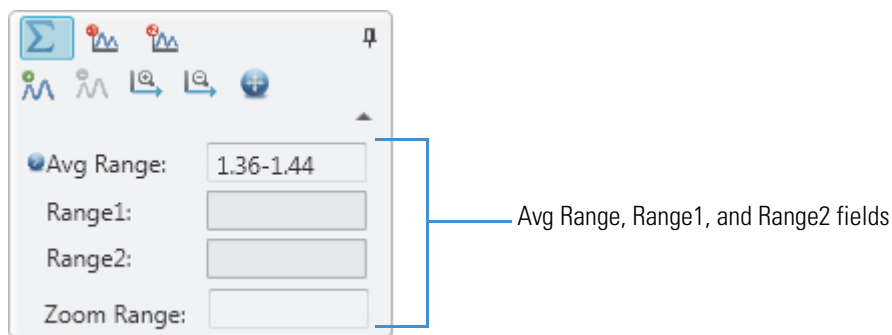
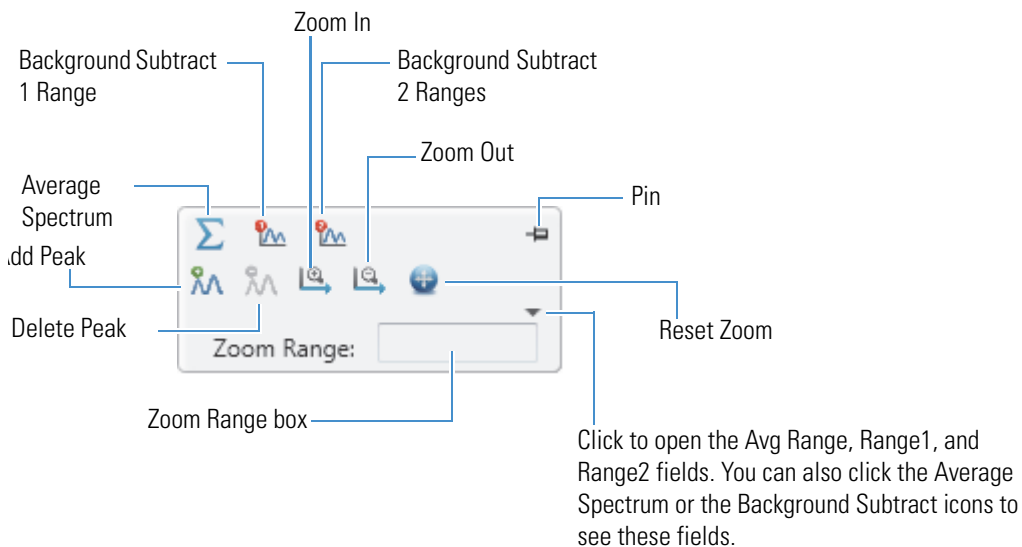


Table 19. Chromatogram toolbox icons and text boxes (Sheet 1 of 2)










Icon/ Field	Description
 Average Spectrum and the Ave Range box	Averages the mass spectra over a selected retention-time range. To create an average spectrum, see Averaging Spectra .
 Background Subtract 1 Range icon and the Range 1 box	Subtracts the mass spectrum of one retention-time range from the active mass spectrum. To create a background subtracted spectrum, see Subtracting Background Spectra .
 Background Subtract 2 Ranges and the Range 1 and Range 2 boxes	Subtracts the mass spectra of two retention-time ranges from the active mass spectrum. To create a background subtracted spectrum, see Subtracting Background Spectra .

Table 19. Chromatogram toolbox icons and text boxes (Sheet 2 of 2)

Icon/ Field	Description
 Pin	When the pin is in the horizontal position, the toolbox auto-hides when you move the pointer away from the current location. To keep the toolbox visible, click the pin icon to change it to a vertical pin.
 Add Peak	Adds a peak manually to the peak list in the Peaks List View (see Manually Adding, Undoing, and Deleting Chromatographic Peaks). Clicking the Add Peak icon again turns off the Add Peak pointer, and resumes the mouse pointer's zoom functionality.
 Delete Peak	Deletes a manual peak from the peak list (see Deleting Chromatographic Peaks). Clicking the Delete Peak icon again turns off the Delete Peak pointer, and resumes the mouse pointer's zoom functionality. Available when the peaks list includes a peak.
Tip You can also use the Add Peak and Delete Peak buttons in the Chromatogram – Workspace Processing toolbar to add and delete manual chromatographic peaks.	
 Zoom In	Displays a smaller portion of the <i>x</i> axis.
 Zoom Out	Displays a larger portion of the <i>x</i> axis.
 Reset Zoom	Resets the zoom to display the entire chromatogram trace.
Zoom Range	Enter a retention time (RT) range within the chromatogram RT limits and then press ENTER. The view rescales to display the chromatogram within the entered time range.

Chromatogram-Specific Pages in the Info Bar

Before you open a workspace in the FreeStyle window, the Info Bar contains only the Isotope Simulation page (see [Isotope Simulation Page](#)). Opening a raw data file adds the Peak Detection, MSn Browser, and Detector Type pages to the Info Bar; whereas, opening a sequence file opens these pages and the Sequence Files page. The Detector Type page lists the detectors used to acquire data for the active raw data file and the MS Trending selection if the raw data file contains mass spectrometry data.

Note If you click a row in the Filter or Trace Type columns in the Chromatogram view, the application replaces the Detector Type page with the Scan Filters page or the Trace Type page, respectively.

These are the pages in the Info Bar:

- [Detector Type Page](#)
- [Trace Type Page](#)
- [Scan Filters Page](#)
- [Peak Detection Page](#)
 - [Genesis Peak Detection Page](#)
 - [ICIS Peak Detection Page](#)
 - [Avalon Peak Detection Page](#)
 - [PPD Peak Detection Page](#)
- [Sequence File Page](#)

Detector Type Page

The detector type determines what types of traces the Chromatogram view displays. Use the Detector Type page of the Info Bar to select the detector type. The detector types shown on this page are the same as those in the Detector Type list in the Chromatogram Ranges view.

These are the possible detector types:

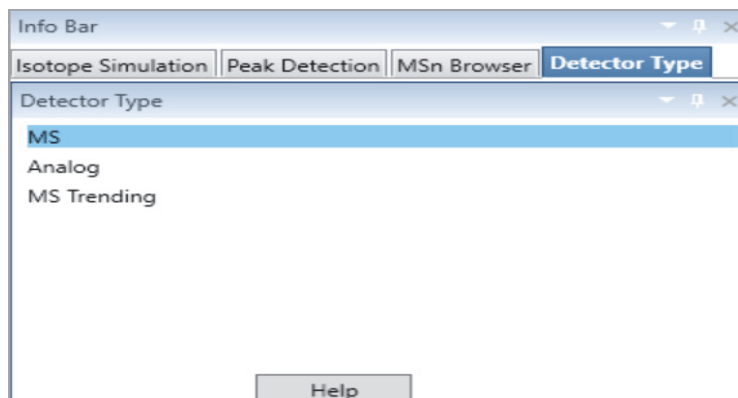
- MS—Mass spectrometry data
- A/D Card—Data from an analog-to-digital converter
- MS Trending—Instrument status readings for the mass spectrometer
- UV—Data from an analog detector
- PDA—Wavelength scan data from a photodiode array detector

❖ To display the Detector Type page

Click the **Detector Type** tab in the Info Bar.

[Figure 61](#) shows a typical Detector Type page.

Figure 61. Detector Type page



❖ **To display a plot in the Chromatogram view**

On the Detector Type page, do one of the following:

- To display an unfiltered plot of the MS data, click **MS**.
- To display a total scan plot for the PDA data, click **PDA**.
- To display a plot for a UV channel, click **UV**.
- To display a plot for the data acquired with an analog-to-digital converter, click **A/D Card**.
- To display a plot for one of the MS detector's status readbacks, click **MS Trending**, click a row in the Trace Type column in the Chromatogram Ranges view, and then, on the Trace Type page, click a trace type (instrument readback).

Trace Type Page

Use the Trace Type page to select the type of trace that the Chromatogram view displays. The available trace types depend on which detector type you select on the Detector Type page:

- For MS detectors, choose the TIC, mass range, base peak, neutral fragment, or mass defect filter trace type.
- For PDA detectors, choose the wavelength range, total scan, or spectrum maximum trace type.
- For analog detectors, choose the analog 1, 2, 3, or 4 trace type.
- For MS Trending, choose any of the instrument status selections.

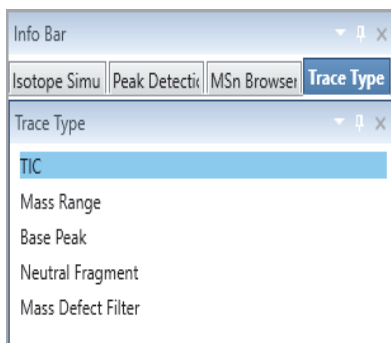
The trace types that the Trace Types page lists are the same as in the Trace Type list in the Chromatogram Ranges view.

❖ **To display the Trace Type page**

1. Open the Chromatogram Ranges view.
2. Click the dropdown arrow in the Trace Type column.

Figure 62 shows trace types for an MS detector type.

Figure 62. Trace Type page



Scan Filters Page

Use the Scan Filters page to select the type of scan filter, if any, that the application applies to the raw data file. The filter types on the Scan Filters page are the same as in the Filter box dropdown list in the Chromatogram Ranges view.

For instructions on using the Scan Filters page, see [Using the Scan Filters Page to Display a Filtered Chromatogram](#).

Note For grouped filters (MS, MS2, HCD, and so on), the Scan Filters page displays only the first filter in the group. The scan header in the Spectrum view displays the actual filter for the scan.

❖ **To display the Scan Filters page in the Info Bar**

In the Workspace Options toolbar, in the Filter area, click **Scan Filters**.

Figure 63 shows a Scan Filters page.

Figure 63. Scan Filters page

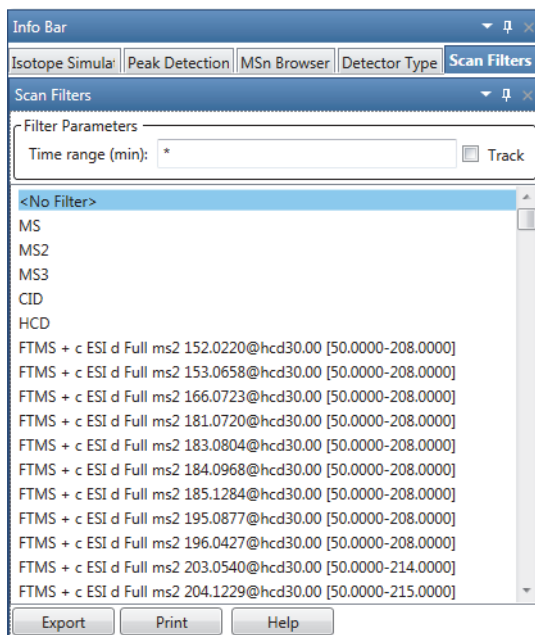


Table 20 describes the parameters on the Scan Filters page.

Table 20. Scan Filters page parameters (Sheet 1 of 2)

Parameter	Description
Filter Parameters	
Time Range (min)	Specifies the time range to display in the filter list. Filters outside the specified time range do not appear in the filter list. The specified time range must be within the retention time range for the data file. Use an asterisk to specify the full retention time range.
Track	Select to set up the time range by dragging the pointer across the Time axis in the Chromatogram view.
Filter list	
<No Filter>	Select to display an unfiltered chromatogram.
MS	Select to display only full-scan MS data.
MS _n (where n = 2 or higher)	Select to display MS/MS data of the selected MS/MS order only.
Scan filters	Select to display a chromatogram for a specific scan filter.
Buttons	
Export	Copies the file name of the raw data file and the contents of the Scan Filters page to the Clipboard.

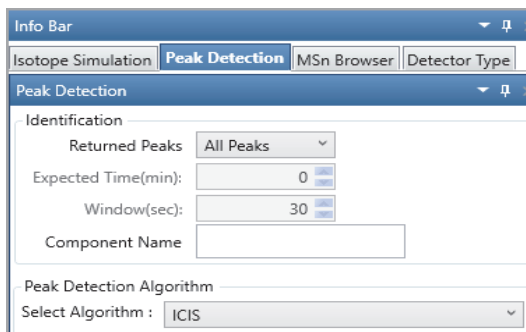
Table 20. Scan Filters page parameters (Sheet 2 of 2)

Parameter	Description
Print	Opens the Print dialog log box for selecting the printer and print preferences for printing the contents of the Scan Filters page.
Help	Opens the Help to the Scan Filters page topic.

Peak Detection Page

Use the Peak Detection Page to identify and integrate the chromatogram peaks in an active chromatogram plot. Specify a peak detection algorithm that the application applies to detect the chromatogram peaks in the active chromatogram plot. For instructions on using the Peak Detection Page, see [Automatically Detecting and Integrating Chromatographic Peaks](#). [Figure 64](#) shows the parameters for the peak identification and detection algorithms.

Figure 64. Peak Detection Info Page



[Table 22](#) describes the parameters for the peak identification and detection algorithms.

Table 21. Peak Identification and Detection algorithm parameters

Parameter	Description
Identification	
Returned Peaks	Type of peak detection results to be returned: <ul style="list-style-type: none"> All Peaks: Displays all the peaks based on the selected algorithm for component identification. Nearest Peak: Displays the peak with the nearest retention in the chromatogram within the specified retention time window for component identification. Highest Peak: Displays the peak with the highest retention in the chromatogram within the specified retention time window for component identification.

Table 21. Peak Identification and Detection algorithm parameters

Parameter	Description
Expected Time (min)	Expected retention time for the component. Enables if you select the Nearest or Highest peak.
Window (sec)	Retention time window for the component. Enables if you select the Nearest or Highest peak.
Component Name	Specifies the component name used for processing. Automatically imports the component name when a processing method is imported to the trace.
Peak Detection Algorithm	
Selected Algorithm	Specifies the peak detection algorithm to be used for processing the chromatographic peaks: <ul style="list-style-type: none">• Genesis—Use for mass spectrometry data.• ICIS—Use for mass spectrometry data.• Avalon—Use for UV or PDA data. This algorithm detects negative chromatographic peaks more accurately than the ICIS or Genesis algorithms.• PPD—Use to minimize the number of optimization parameters.

Genesis Peak Detection Page

The Genesis peak detection algorithm is the original Thermo Xcalibur peak detection algorithm and has been provided for backward compatibility with Xcalibur 1.0 data system studies. Use the Genesis Peak Detection page of the Info Bar to specify peak identification and peak integration criteria that the Genesis algorithm applies toward chromatograms in the Chromatogram view.

For information about using the Genesis peak detection algorithm, see [Automatically Detecting and Integrating Chromatographic Peaks](#).

❖ To display the Genesis Peak Detection page

1. Click the **Peak Detection** tab in the Info Bar.
2. In the list, select **Genesis**.

[Figure 65](#) shows the parameters for the Genesis peak detection algorithm.

Figure 65. Genesis Peak Detection page

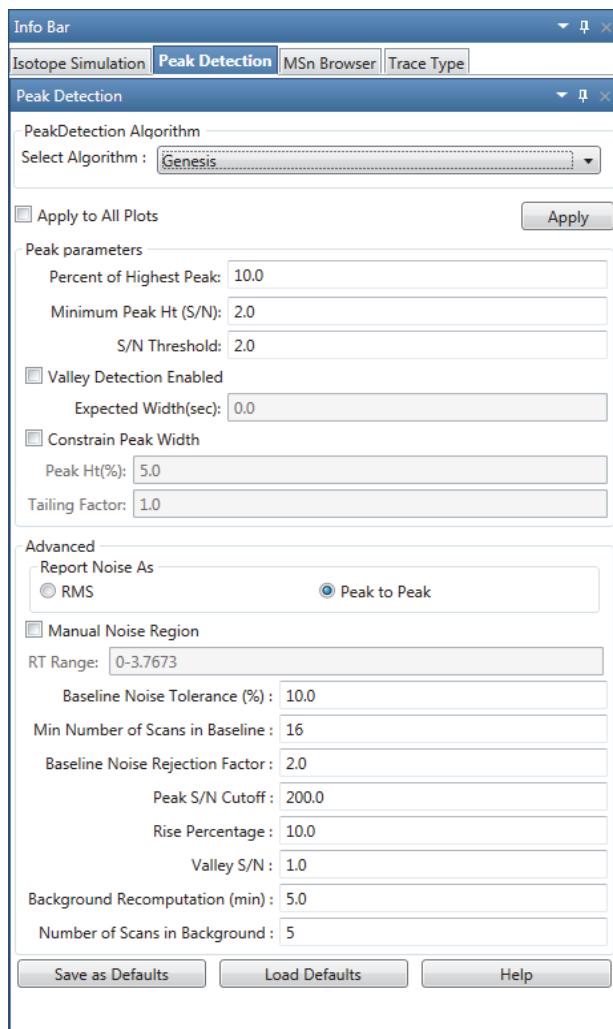


Table 22 describes the parameters for the Genesis Peak Detection page.

Table 22. Genesis peak detection parameters (Sheet 1 of 4)

Parameter	Description
Application of Settings	
Apply to All Plots	Apply the peak identification and integration settings to all displayed chromatograms. To apply the criteria to only the active chromatogram, clear this option.

Table 22. Genesis peak detection parameters (Sheet 2 of 4)

Parameter	Description
Peak Parameters	
Percent of Highest Peak	Specify a percentage threshold to limit the number of peaks to be submitted for further processing. The algorithm discards any detected peak that has an intensity less than the threshold percentage of the most intense peak. Range: 0.0–100.0; default: 10.0
Minimum Peak Ht (S/N)	Specify the peak signal-to-noise (S/N) value to equal or exceed as a criterion for peak detection. The algorithm ignores all chromatogram peaks that have S/N values less than the minimum peak height S/N value. Range: 1.0 (all peaks) to 999.0; default: 2.0
S/N Threshold	Specify the threshold for detecting peak edges. The algorithm calculates the S/N ratio by using only the baseline signal. It excludes any extraneous or minor detected peaks from the calculation. Range: 0.0–999.0; default: 2.0
Valley Detection Enabled	Select the valley detection approximation method to detect unresolved peaks. This method drops a vertical line from the apex of the valley between unresolved peaks to the baseline. The intersection of the vertical line and the baseline defines the end of the first peak and the beginning of the second peak.
Expected Width (sec)	Specify the expected peak width (in seconds). This value controls the minimum width that a peak must have if you selected the Valley Detection Enabled option. With valley detection selected, the algorithm ignores any valley points nearer than the <i>expected width/2</i> to the top of the peak. If it finds a valley point outside the expected peak width, the algorithm terminates the peak at that point. The algorithm always terminates a peak when the signal reaches the baseline, independent of the value set for the expected peak width. Range: 0.0–999.0; default: 0.0
Constrain Peak Width	Limits the peak width of a component during peak integration. To control when peak integration is turned on and off, specify a peak height threshold and a tailing factor.

Table 22. Genesis peak detection parameters (Sheet 3 of 4)

Parameter	Description
Peak Ht (%)	Specify the peak height (as a percentage) where the algorithm tests the width of target peaks. Range: 0.0–100.0; default: 5.0
Tailing Factor	Specify a factor that controls how the algorithm integrates the tail of a peak. This factor is the maximum ratio of the trailing edge to the leading edge of a constrained peak. Range: 0.5–9.0; default: 1.0
Advanced	
RMS	Calculate noise as a root-mean-square (RMS) value.
Peak to Peak	Calculate noise as a peak-to-peak value.
Manual Noise Region	Specify a region of the chromatogram that the algorithm uses to determine noise. See Selecting the Manual Noise Region for the Genesis and ICIS Algorithms .
RT Range	Specify the RT range that the algorithm uses to determine noise. The RT range must be within the chromatogram range.
Baseline Noise Tolerance (%)	Specify a percentage value that controls how the algorithm draws the baseline in the noise data. The higher the baseline noise tolerance value, the higher it draws the baseline through the noise data. Range: 0.0–100.0; default: 10.0
Min Number of Scans in Baseline	Specify the minimum number of scans that the algorithm uses to calculate a baseline. A larger number includes more data in determining an averaged baseline. Range: 2–100; default: 16
Baseline Noise Rejection Factor	Specify the baseline noise rejection factor. This factor controls the width of the RMS noise band above and below the peak detection baseline. The algorithm applies this factor to the raw RMS noise values to raise the effective RMS noise that the algorithm uses during peak detection. It responds by assigning the left and right peak boundaries above the noise and, therefore, closer to the peak apex value. This action effectively raises the peak integration baseline above the RMS noise level. Range: 0.1–10.0; default: 2.0

Table 22. Genesis peak detection parameters (Sheet 4 of 4)

Parameter	Description
Peak S/N Cutoff	Specify the S/N level that the algorithm defines as the top of the peak edge. For example, if the S/N level at the apex is 500 and the Peak S/N Cutoff value is 200, the algorithm defines the right and left edges of the peak when the S/N reaches a value less than 200. Range: 50.0–10 000.0; default: 200.0
Rise Percentage	Specify how much above the baseline the trace can rise (as a percentage) after passing through a minimum inflection point in the trace (before or after the peak). If the trace exceeds this value, the algorithm applies valley detection peak integration criteria. It applies this test to both the left and right edges of the peak. The rise percentage criterion is useful for integrating peaks with long tails. Range: 0.1–500.0; default: 10.0
Valley S/N	Specify the S/N criterion that the algorithm uses for valley detection. Range: 1.0–100.0; default: 1.0
Background Recomputation (min)	Specify the background recomputation interval (in minutes). The algorithm periodically recalculates the representative background scan it uses for background subtraction to compensate for the possibility that the composition of the background might change over the course of a run. The background recomputation interval is the time interval between these recalculations. Range: 0.5–10.0; default: 5.0
Number of Scans in Background	Specify the number of background scans the algorithm uses to determine the background. Range: 2–100; default: 5
Button	
Apply	Starts peak detection by using the Genesis peak detection algorithm.
Save As Defaults	Saves the current settings as default settings. After you save the settings as defaults, you can restore these values at any time by clicking Load Default.
Load Defaults	Restores the current default settings.
Help	Opens the product-name Help to the Genesis peak detection topic.

ICIS Peak Detection Page

The ICIS peak detection algorithm is the default algorithm that was designed for MS data. It is known for peak detection efficiency at low MS signal levels.

Use the ICIS Peak Detection page of the Info Bar to specify peak detection and integration criteria. The application applies the peak detection algorithm to the active plot or all the plots in the Chromatogram view.

For more information, see [Automatically Detecting and Integrating Chromatographic Peaks](#).

❖ To display the ICIS peak detection parameters

1. Click the **Peak Detection** tab in the Info Bar.
2. In the list, select **ICIS**.

Figure 66 shows the parameters for the ICIS peak detection algorithm.

Figure 66. ICIS Peak Detection page

The screenshot shows the 'Peak Detection' page in the Info Bar. The 'Peak Detection Algorithm' is set to 'ICIS'. The 'Apply to All Plots' checkbox is checked. The 'Peak parameters' section includes: Baseline Window (40), Area noise factor (500), Peak noise factor (10), Constrain Peak Width (checked), Peak Ht(%) (5.0), and Tailing Factor (1.0). The 'Advanced' section includes: Manual Noise Region (checked) with RT Range (0-6.97), Noise Method (INCOS Noise selected), RMS (checked) with Min peak width (3), Multiplet resolution (10), Area tail extension (5), and Area scan window (0). Buttons for 'Save as Defaults', 'Load Defaults', and 'Help' are at the bottom.

Table 23 describes the parameters for the ICIS peak detection algorithm.

Table 23. ICIS Peak Detection page parameters (Sheet 1 of 2)

Parameter	Description
Application of Settings	
Apply to All Plots	Apply the current chromatogram peak identification and integration settings to all displayed plots. To apply the criteria to only the active plot, clear this option.
Peak Parameters	
Baseline Window	Specify the number of scans, from the apex down each side of the peak, that the peak detection algorithm uses to determine the minimum baseline for the peak. A higher number of scans means a wider scan range, resulting in a lower baseline. Range: 1–500; default: 40
Area Noise Factor	Specify the noise level multiplier that the peak detection algorithm uses to determine the peak edge after the location of the possible peak. Range: 1–500; default: 500
Peak Noise Factor	Specify the noise level multiplier that the peak detection algorithm uses to determine the potential peak signal threshold. Range: 1–1000; default: 10
Constrain Peak Width	Limits the peak width of a component during the peak integration of a chromatogram. You can then set values that control when peak integration turns on and off by specifying a peak height threshold and a tailing factor.
Peak Ht (%)	Specify how much above the baseline (as a percentage) a signal must be of the total peak height before integration turns on or off. Range: 0.0–100.0; default: 5.0
Tailing Factor	Specify a factor that controls how the algorithm integrates the tail of a peak. This factor is the maximum ratio of the trailing edge to the leading edge of a constrained peak. Range: 0.5–9.0; default: 1.0
Advanced	
Manual Noise Region	Specify a region of the chromatogram that the peak detection algorithm uses to determine noise. You can enter the retention time (RT) in the RT Range box. See Selecting the Manual Noise Region for the Genesis and ICIS Algorithms .

Table 23. ICIS Peak Detection page parameters (Sheet 2 of 2)

Parameter	Description
RT Range	Specify the RT range that the algorithm uses to determine noise. The RT range must be within the chromatogram range.
INCOS Noise	Use a single pass algorithm to determine the noise level.
Repetitive Noise	Use a multiple pass algorithm to determine the noise level. In general, this algorithm is more accurate in analyzing the noise than the INCOS Noise algorithm; however, it might take longer depending on the data.
RMS	Calculate noise as root mean square (RMS). By default, the algorithm uses peak-to-peak for the noise calculation. If you determine the noise region manually, the algorithm automatically selects RMS.
Min Peak Width	Enter the minimum number of scans required in a peak. Range: 0–100; default: 3
Multiplet Resolution	Specify the minimum separation in scans between the apexes of two potential peaks. This criterion determines if two peaks are resolved. Range: 1–500; default: 10
Area Tail Extension	Specify the number of scans past the peak endpoint to use in averaging the intensity. Range: 0–100; default: 5
Area Scan Window	Specify the number of scans on each side of the peak apex to include in the area integration. Range: 0–100; default: 0
	Note Specifying zero scans means that the algorithm includes all scans in the area integration, from the peak start to the peak end.
Button	
Apply	Starts the ICIS peak detection algorithm.
Save As Defaults	Saves the current settings as the default settings.
Load Defaults	Restores the current default settings.
Help	Opens the product-name Help to the ICIS peak detection topic.

Avalon Peak Detection Page

The Avalon peak detection algorithm supports detectors other than mass spectrometers, and detects negative chromatographic peaks more accurately than the Genesis or ICIS peak detection algorithms. Use the Avalon Peak Detection page in the Info Bar to specify peak identification and integration criteria. You can apply this algorithm to the active plot or all the plots in the Chromatogram view.

❖ **To display the Avalon Peak Detection page**

1. Click the **Peak Detection** tab in the Info Bar.
2. Select the **Avalon** search algorithm.

Figure 67 shows the parameters for the Avalon peak detection algorithm.

Figure 67. Avalon Peak Detection Settings page

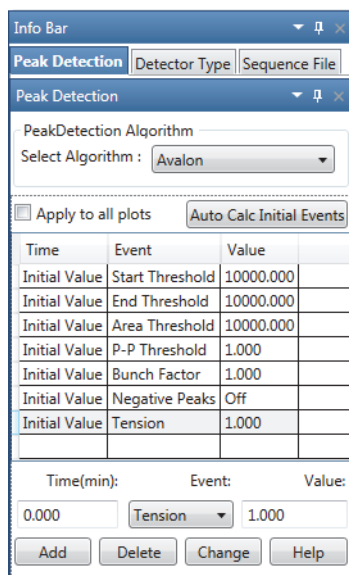


Table 24 describes the parameters for the Avalon peak detection algorithm.

Table 24. Avalon Peak Detection page parameters (Sheet 1 of 4)

Parameter	Description
Application of Settings	
Apply to All Plots	Apply the current chromatogram peak identification and integration settings to all displayed chromatogram plots. To apply the criteria to only the active plot, clear this check box.

Table 24. Avalon Peak Detection page parameters (Sheet 2 of 4)

Parameter	Description
Event List	
Event List	<p>Specify the settings for initial events and user-defined timed events in the event list.</p> <p>To calculate values for initial events, click Auto Calculate Initial Events.</p> <p>To change the settings in the event list, highlight the row and then enter the revised settings in the boxes below the list. Click Add to add a new row of entered values to the event list. Click Change to update automatically both the event list and the chromatogram display.</p> <p>There are seven permanent integration events, identified by the Initial Value setting in the Time column. These are the default integration events that the Avalon integration algorithm requires. You can change the value of an initial entry integration event, but you cannot delete it or change its time value.</p>
Event List Entry	
Time	<p>Specify the value for the Time column for the highlighted entry in the event list.</p> <p>You cannot change the time entry for initial value events.</p>

Table 24. Avalon Peak Detection page parameters (Sheet 3 of 4)

Parameter	Description
Event	<p>Select the event type from these options:</p> <ul style="list-style-type: none"> • Start/End Threshold: Directly related to the RMS noise in the chromatogram, this value is the start or end threshold, the fundamental control used for peak detection. • Area Threshold: Controls the area cutoff. The algorithm does not detect any peaks with a final area less than the area threshold. This control is in units of area for the data. • P-P Threshold: The peak-to-peak resolution threshold controls how much peak overlap must be present before two or more adjacent peaks create a peak cluster. Peak clusters have a baseline drop instead of valley-to-valley baselines. This option is specified as a percentage of the peak height overlap. • Negative Peaks: Allows or denies negative peaks. • Bunch Factor: The number of points that are grouped together during peak detection. It controls the bunching of chromatographic points during integration and does not affect the final area calculation of the peak. The Bunch Factor must be an integer between 1 and 6; a high bunch factor groups peaks into clusters. • Tension: Controls how closely the baseline must follow the overall shape of the chromatogram. A lower baseline tension traces the baseline to follow changes in the chromatogram more closely. A higher baseline tension follows the baseline less closely over longer time intervals. This option is specified in minutes. • Tangent Skim: For fused peaks that are significantly different in size, the tangent skim method allocates area to the various peaks. By default, the algorithm chooses the tallest peak in a cluster as the parent. You can also identify which peak in the cluster is the parent. The algorithm detects tangent skim peaks on one or both sides of the parent peak. Tangent skim automatically resets at the end of the peak cluster. <p>The threshold and bunch factor parameters are the most important ones in controlling peak detection.</p>
Value	<p>Specify the value for Value column for the currently highlighted entry in the event list. The range of factors allowed for each value is specific to each event.</p>

Table 24. Avalon Peak Detection page parameters (Sheet 4 of 4)

Parameter	Description
Button	
Auto Calc Initial Events	Searches for the best values of initial events that detect peaks in the data. This button is active if you have a raw data file open. The Avalon algorithm automatically estimates the initial values for the detection of peaks based on the data in the current raw data file, and then displays those initial values in the event list. Any timed event in the event list is unchanged. Automatic calculation of initial events determines initial values for the following events only: start threshold, end threshold, area threshold, P-P [resolution] threshold, bunch factor, negative peaks, and tension. Additionally, you can specify timed events for these events in the same event list.
Add	Add a row to the event list. When you click Add, both the event list and the chromatogram update automatically with the added specification in the currently selected chromatogram.
Delete	Removes a highlighted event from the event list. You cannot delete initial values.
Change	Updates a highlighted entry in the event list. When you click Change, peak detection with the updated specification occurs automatically for the currently selected chromatogram. For initial events, the algorithm changes only the values, and not the events.
Help	Opens the product-name Help to the Avalon peak detection topic.

PPD Peak Detection Page

The PPD peak detection algorithm fits the data to model peak functions.

Use the PPD Peak Detection page of the Info Bar to specify peak detection criteria. The application applies the peak detection algorithm to the active raw data file displayed in the product-name window.

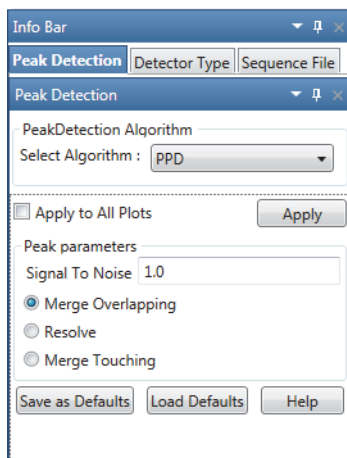
For information about using the PPD peak detection algorithm, see [Automatically Detecting and Integrating Chromatographic Peaks](#).

❖ To display the PPD Peak Detection page

1. Click the **Peak Detection** tab in the Info Bar.
2. Select the **PPD** algorithm.

[Figure 68](#) shows the settings for the PPD peak detection algorithm.

Figure 68. PPD Peak Detection page



For information about adding peak labels, see [Labeling Chromatogram Peaks or Local Maxima](#).

[Table 25](#) describes the parameters for the PPD peak detection algorithm.

Table 25. PPD Peak Detection page parameters (Sheet 1 of 2)

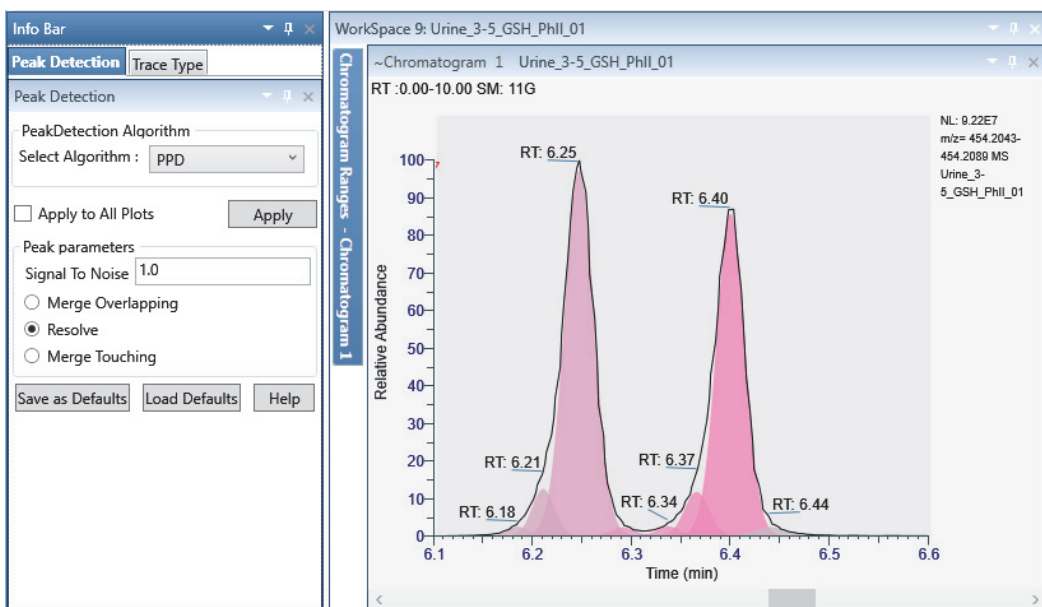
Parameter	Description
Application of Settings	
Apply to All Plots	Applies the current chromatogram peak identification and integration settings to all displayed plots. To apply the criteria to only the active plot, clear this option.
Peak Parameters	
Signal To Noise	Specifies the absolute signal-to-noise threshold. A value of 0 applies no filtering. Range: 0.00–100.00; default: 1.0
Merge Overlapping	(Default) Merges overlapping peaks into one peak. Use this option for LC/MS data to avoid reporting small shoulders as separate peaks.
Resolve	Resolves overlapping or touching peaks into two chromatographic peaks. Use this option for GC/MS data.
Merge Touching	Merges two peaks whose edges barely overlap at the 1% intensity level.
Button	
Apply	Starts the PPD peak detection algorithm.
Save As Defaults	Saves the current settings as the default settings.

Table 25. PPD Peak Detection page parameters (Sheet 2 of 2)

Parameter	Description
Load Defaults	Restores the current default settings.
Help	Opens the product-name Help to the PPD peak detection topic.

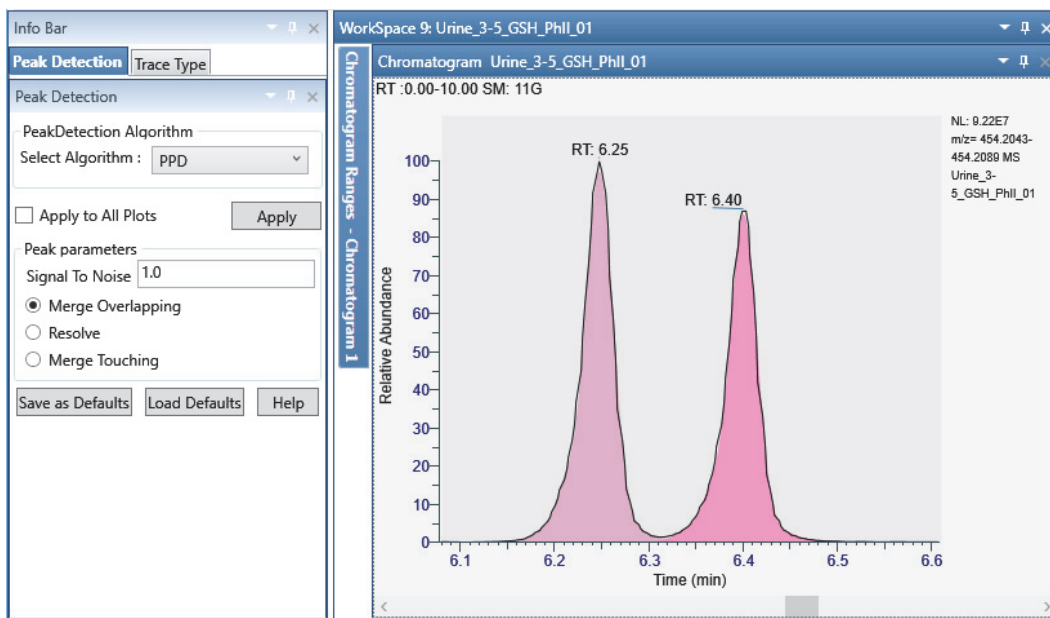
The PPD peak detection algorithm automatically resolves all the chromatographic peaks, and selecting the Resolve option displays all the detected peaks (see [Figure 69](#)).

Figure 69. PPD peak detection with the Resolve option



By default, the Merge Overlapping option is selected. For LC/MS data, use Merge Overlapping to avoid reporting small shoulders as separate peaks. Compare [Figure 70](#) where the overlapping peaks are merged and reported as two peaks to [Figure 69](#) where all the peaks are resolved and reported as seven separate peaks.

Figure 70. PPD peak detection with the Merge Overlapping option



Sequence File Page

The Sequence File page of the Info Bar displays a sequence of raw data files that you created or opened. Click a raw data file in the list to display its trace in the Chromatogram view.

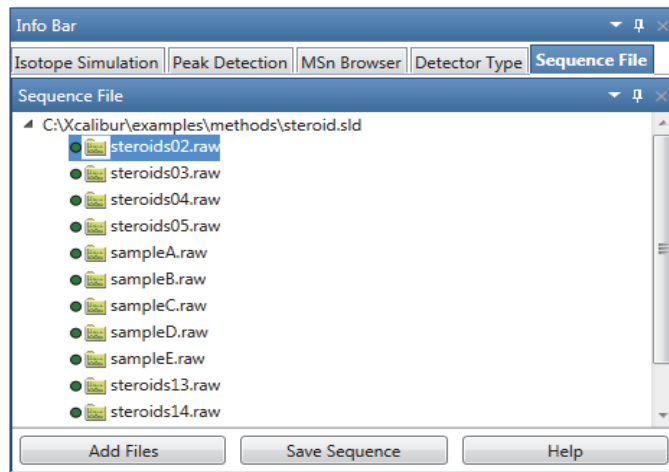
❖ To display the Sequence File page

Create or open a sequence of raw data files.

For more information, see [Creating, Positioning, Previewing, Refreshing, and Closing Workspaces](#) or [Working with Sequences](#).

Figure 71 shows a typical Sequence File page.

Figure 71. Sequence File page



Reviewing Spectral Data

Use the spectra-specific toolbars, views, and Info Bar pages of the FreeStyle application to review spectral data from the mass spectrometer or PDA detector.

To display spectra of interest, follow these topics.

Contents

- [Displaying the Scan for a Time Point in a Chromatogram](#)
- [Adding Spectrum Views to the Workspace](#)
- [Selecting Spectra from the Spectrum Ranges Dialog Box](#)
- [Creating a MultiSpectrum View and Changing Its Spectrum Plots](#)
- [Formatting a MultiSpectrum View](#)
- [Linking and Releasing Spectra to and from a Chromatogram](#)
- [Averaging Spectra](#)
- [Subtracting Background Spectra](#)
- [Exporting a Selected Spectrum to a New RAW File](#)
- [Using Precursor Markers to Open Data-Dependent Scans](#)
- [Selecting Spectra from an MSn Tree](#)
- [Setting Up the Display Options for a Spectrum View](#)
- [Changing the Zoom Level in a Spectrum View or MultiSpectrum View](#)
- [Spectra-Specific Toolbars](#)
- [Spectra-Specific Views](#)
- [Spectrum Ranges Dialog Box](#)
- [Spectrum Toolbox](#)
- [MSn Browser Page](#)

4 Reviewing Spectral Data

Displaying the Scan for a Time Point in a Chromatogram

Note For information about scan headers and custom annotations, see [Scan Headers and Scan Header Abbreviations](#) and [Adding Text, Graphic, and Structure Annotations to a Graphical View](#), respectively.

Displaying the Scan for a Time Point in a Chromatogram


Follow this procedure to display the scan for a particular time point of a chromatogram.

❖ To display the scan corresponding to a particular time point in a chromatogram

1. Display the chromatogram plots of interest as described in [Defining a Chromatogram Trace from the Chromatogram Ranges View](#).
2. Make sure that the Spectrum view or Multi Spectrum view links to the chromatogram.

Note The text appended to a Spectrum view indicates its linkage status. For example, the text CIT1 at the end of a Spectrum view's name indicates that actions performed on trace 1 in Chromatogram view 1 affect the spectrum displayed.

3. Click the chromatogram plot at the retention time or scan number that you want to view.

In the Chromatogram view, a red vertical marker, , indicates the selected data point. The Spectrum view displays the spectrum for that retention time or scan number. On the computer keyboard, use the right or left arrow keys to increment or decrement the scan number.

Note If you do not apply a scan filter, the scan number increments and decrements by one. If you do apply a scan filter, the scan number increments and decrements to the next scan number that meets the filter criteria.

Adding Spectrum Views to the Workspace

You can add multiple Spectrum views and MultiSpectrum views to the workspace.

❖ To add a Spectrum view

1. Open a raw data file (see [Opening Raw Data Files or Sequence Files](#)).

With the factory default layout, the Chromatogram view appears at the top and the Spectrum view appears at the bottom and displays scan #1.

2. To add another Spectrum view to the workspace, do one of the following:

- In the Workspace Options toolbar, click **Spectrum**.

A copy of the active spectrum appears in the new Spectrum view.

- Open the MSn Browser page and select the fragmentation spectra of interest (see [Selecting Spectra from an MSn Tree](#)).

A new Spectrum view opens for each selected spectrum.

For information about displaying a different spectrum in the Spectrum view, see these topics:

- [Displaying the Scan for a Time Point in a Chromatogram](#)
- [Selecting Spectra from the Spectrum Ranges Dialog Box](#)

Selecting Spectra from the Spectrum Ranges Dialog Box

You can use the Spectrum Ranges dialog box to select the spectrum to be displayed in a Spectrum view or the spectra to be displayed in a MultiSpectrum view. For information about creating an average spectrum or performing background subtraction, see these topics: [Creating an Average Spectrum from the Spectrum Ranges Dialog Box](#) or [Using the Spectrum Ranges Dialog Box to Subtract Background Spectra](#).

Note When a spectrum is linked to a chromatogram, you cannot select a different raw data file, detector, or scan filter from the Spectrum Ranges dialog box.

❖ To select spectra with the Spectrum Ranges dialog box

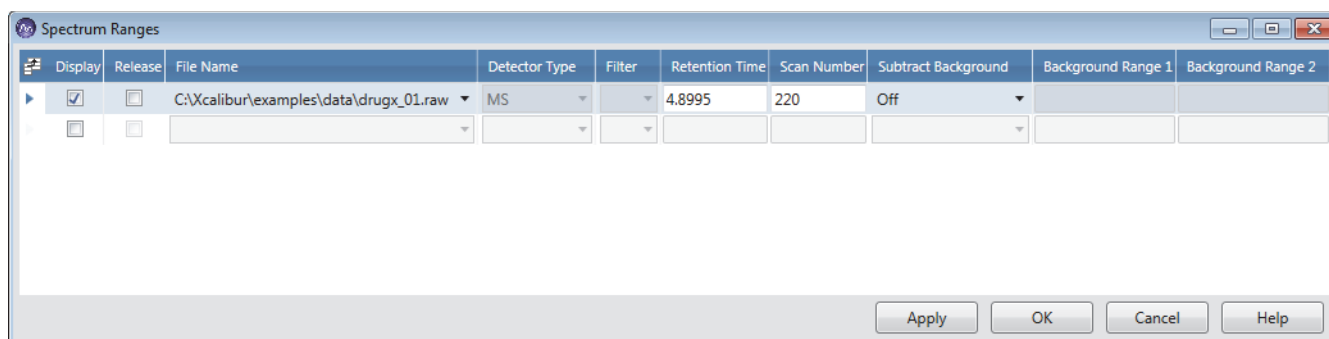
1. Do one of the following:
 - Click either a **Spectrum** view or a **MultiSpectrum** view to select it. Then, in the Workspace Options toolbar, click **Spectrum Ranges**.

–or–

- Right-click a Spectrum view or a MultiSpectrum view and choose **Ranges**.

The Spectrum Ranges dialog box opens with a list of the scans currently displayed in the selected view ([Figure 72](#)).

Figure 72. Spectrum Ranges dialog box



4 Reviewing Spectral Data

Creating a MultiSpectrum View and Changing Its Spectrum Plots

2. Do any of the following:
 - To add another spectrum plot, select the last check box in the Display column.
 - For a MultiSpectrum view, to remove a spectrum plot, clear its associated check box.
 - (Released spectrum only) From the File Name list, browse to and select a different raw data file.
 - (Released spectrum only) From the Detector Type list, select a different detector.
 - (Released spectrum only) From the Filter list, select a different scan filter.
 - In the Retention Time box, type a different retention time, and then click elsewhere in the same row.

The corresponding scan number appears in the Scan Number box.
 - In the Scan Number box, type a different scan number, and then click elsewhere in the same row.

The retention time of the selected scan appears in the Retention Time box.
3. Click **OK**.

If you add multiple plots to a Spectrum view, the Spectrum view becomes a MultiSpectrum view.

Creating a MultiSpectrum View and Changing Its Spectrum Plots

To compare spectrum plots using the same normalization level, you must use the MultiSpectrum view (see [Applying Local or Global Normalization in a MultiSpectrum View](#)).

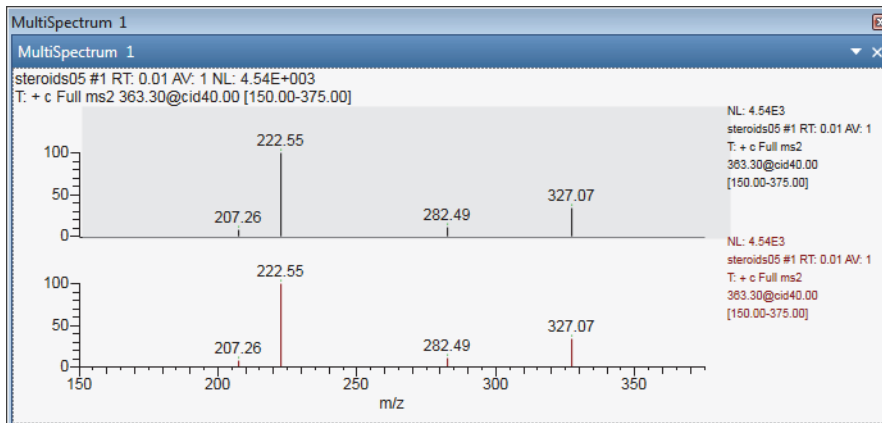
To set up the spectrum plots in a MultiSpectrum view, follow these procedures

- [To change a Spectrum view to a MultiSpectrum view](#)
 - [To add spectrum plots to a MultiSpectrum view](#)
 - [To select a different spectrum for any of the plots in a MultiSpectrum view](#)
- ❖ **To change a Spectrum view to a MultiSpectrum view**

1. Click the **Workspace Options** toolbar tab.
2. Select a Spectrum view and click **Multi Spectrum**.

The view's title bar changes to MultiSpectrum and a copy of the selected spectrum appears below the original spectrum ([Figure 73](#)).

Figure 73. MultiSpectrum view



❖ **To add spectrum plots to a MultiSpectrum view**

1. Click the MultiSpectrum view to select it.
2. In the Workspace Options toolbar, click **Multi Spectrum**.

A copy of the last plot appears at the bottom of the view.

❖ **To select a different spectrum for any of the plots in a MultiSpectrum view**

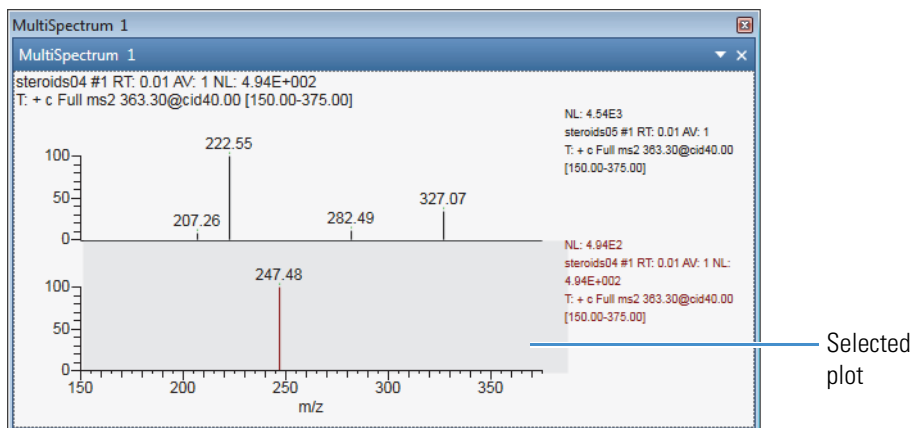
- Use the Spectrum Ranges dialog box (see [Selecting Spectra from the Spectrum Ranges Dialog Box](#)).

—or—

- a. Click the plot to select it.

An active plot has a darker background than the other plots in the view ([Figure 74](#)).

Figure 74. MultiSpectrum view with plots from different raw data files



- b. In the Chromatogram view, select a point by using the pointer or the left and right keyboard keys.

Formatting a MultiSpectrum View

For more information about the formatting options, see [MultiSpectrum – Display Options Toolbar](#).

❖ To format the spectrum plots in a MultiSpectrum view

1. Click the MultiSpectrum view to select it.

The MultiSpectrum toolbar ribbon appears.

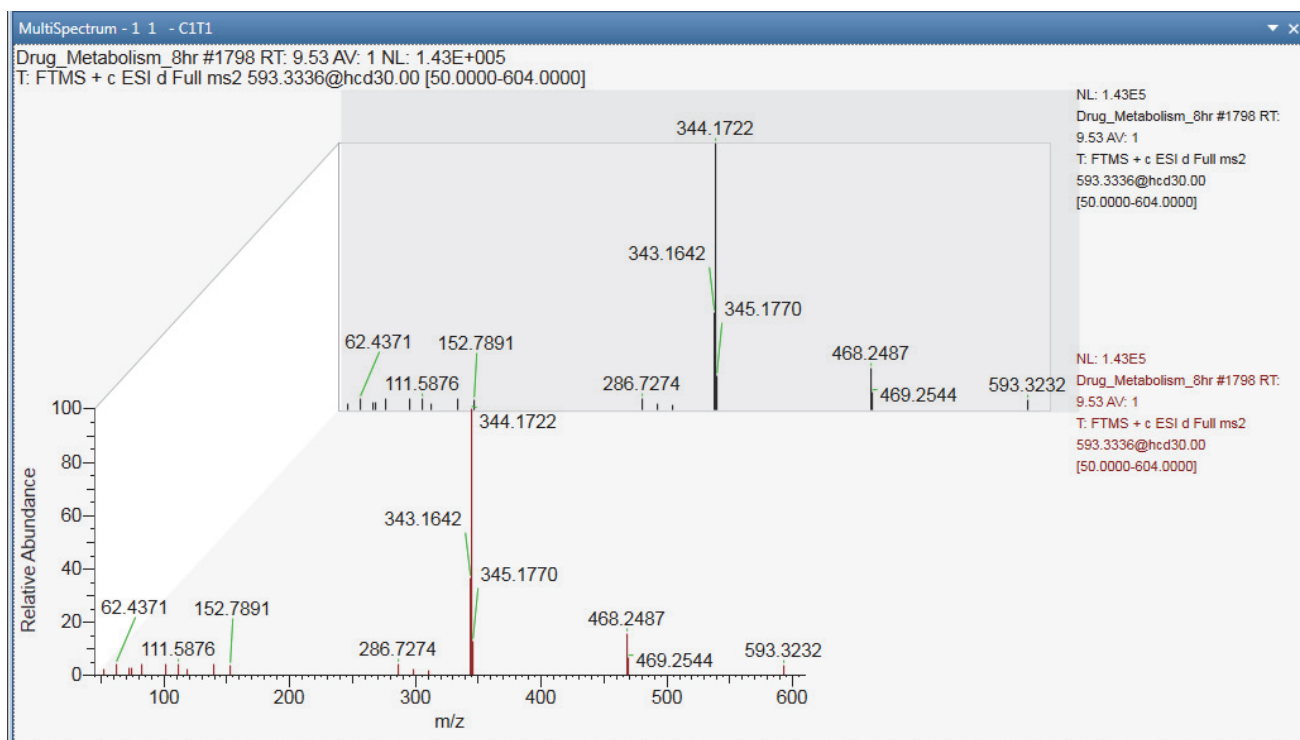
2. Click **Display Options**.

3. Do one or more of the following:

- To change the color of a spectrum plot, select the plot, and then click the **Color** icon in the Format area of the Display Options toolbar and select a color.
- To change the orientation of the plots from stacked to overlaid, click **Overlay**.
- To change the skew of the overlaid plots, use the **Skew** slider.
- To change the relative elevation of the overlaid plots, use the **Elevation** slider.
- To add a backdrop to the overlaid plots, click **Draw Backdrop**.

Figure 75 show a MultiSpectrum view with two spectrum plots with a 25% overlay, a 45° skew, and a backdrop.

Figure 75. MultiSpectrum view with two overlaid spectrum plots



Linking and Releasing Spectra to and from a Chromatogram

When you add a Spectrum view to the Workspace, the application automatically links the spectrum to the active chromatogram trace. You can release the spectrum from the trace and link it to another trace.

See these topics:

- [Linkage States for a Spectrum Plot](#)
- [Releasing a Linked Spectrum from the Chromatogram](#)
- [Linking a Spectrum View to a Chromatogram](#)

Linkage States for a Spectrum Plot

Table 26 describes the linkage states for a spectrum plot.

Table 26. Linkage states for a spectrum plot

Linkage state	Description
Linked to a chromatogram	<p>When a spectrum plot is linked to a chromatogram, clicking another time point in the linked chromatogram changes the displayed spectrum.</p> <p>The title bar of the Spectrum (or Multi Spectrum) view describes the linked chromatogram by its view number and trace number in the following format:</p> $CxTy$ <p>where</p> <p>x = the number of the Chromatogram view</p> <p>y = the number of the trace in the Chromatogram view</p>
Available for linking	<p>The following text is appended to the title bar of the Spectrum (or Multi Spectrum) view—Select Trace to Link.</p> <p>When a spectrum plot is available for linking, clicking any chromatogram in the Workspace links the plot to the selected chromatogram, updates the spectrum plot to the selected time point, and changes the appended text in the view's title bar to $CxTy$.</p>
Released	<p>When you release a spectrum plot from its chromatogram, the text appended to the Spectrum (or Multi Spectrum) view changes to Released.</p> <p>In the Released state, the spectrum plot is not affected by any actions in a Chromatogram view. To update a released spectrum plot, you must use the Spectrum Ranges dialog box.</p>

Releasing a Linked Spectrum from the Chromatogram

For a spectrum in a Spectrum view, the text appended to the view's name indicates the linkage status of the spectrum. For a MultiSpectrum view, the appended text applies to the selected spectrum.

❖ To release a spectrum from its linked chromatogram

Right-click the spectrum and choose **Release from Chromatogram**.

The application appends the following text to the view's name: Released.

Note After you release a spectrum, you can use the Spectrum Ranges dialog box to select a different detector, scan filter, or raw data file.

Linking a Spectrum View to a Chromatogram

❖ To link a spectrum to a chromatogram

1. Right-click the spectrum and choose **Link to Chromatogram**.
2. Click the chromatogram to link it.

To describe the linkage state, the application appends the following text to the view's name:

C_xT_y

where x indicates the Chromatogram view's number and y indicates the trace number

Note When a spectrum is linked to a chromatogram, you cannot use the Spectrum Ranges dialog box to select a different detector, scan filter, or raw data file.

Averaging Spectra

You can increase mass accuracy and reduce noise in a spectrum by averaging the spectrum over an appropriate scan range.

To generate an averaged spectrum, the application uses the original intensity versus frequency data from the mass spectrometer, bins the data into frequency intervals, and then uses the instrument's mass calibration file to convert the frequency data to m/z values.

You can create an averaged spectrum by using the Chromatogram toolbox or the Spectrum Ranges dialog box.


See these topics:

- [Creating an Average Spectrum from the Chromatogram Toolbox](#)
- [Creating an Average Spectrum from the Spectrum Ranges Dialog Box](#)

Creating an Average Spectrum from the Chromatogram Toolbox

When a spectrum is linked to a chromatogram trace, you can use the Chromatogram toolbox to create an average spectrum.

❖ To create an average spectrum by using the Chromatogram toolbox

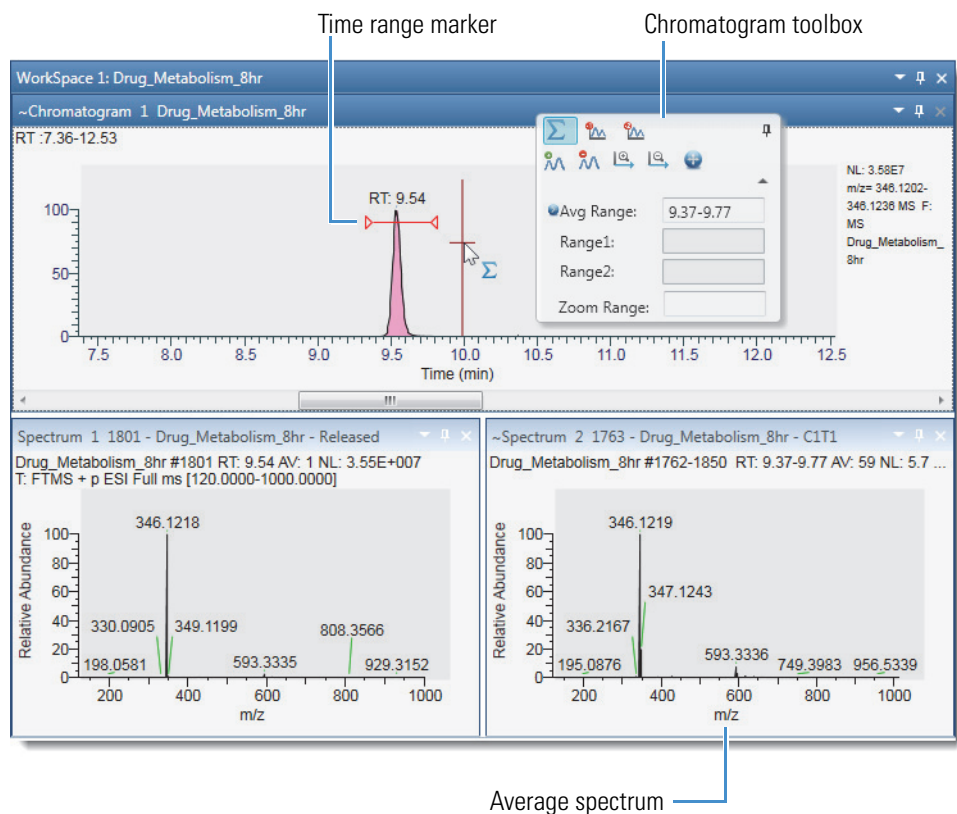
1. In the Chromatogram view that is linked to the spectrum, open the Chromatogram toolbox.
2. Click the **Average Spectrum** icon, .
3. Drag the pointer through the time range that defines the scans to average, or enter the time range in the Avg Range box.

A red line in the Chromatogram view marks the time range. You can undo spectrum averaging by clicking the Average Spectrum icon again.

[Figure 76](#) shows a Workspace view with two Spectrum views. The Spectrum view on the right shows the average spectrum for the time range defined in the toolbox.

Tip To retain the average spectrum while working with the chromatogram, release the spectrum from the chromatogram.

Figure 76. Workspace with a Chromatogram view and two Spectrum views



Creating an Average Spectrum from the Spectrum Ranges Dialog Box

❖ **To create an average spectrum by using the Spectrum Ranges dialog box**

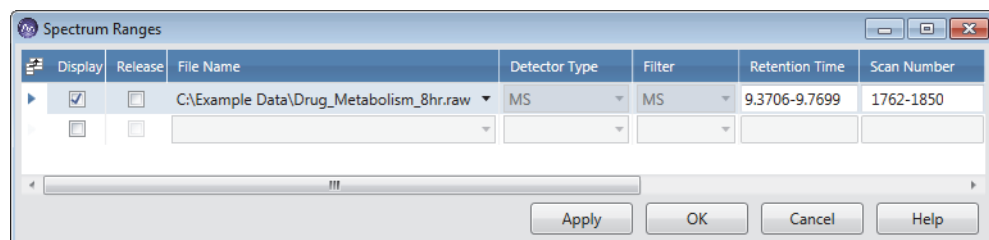
1. Right-click the Spectrum view of interest and choose **Ranges**.

The Spectrum Ranges dialog box opens.

2. In the Retention Time box, enter the time range for averaging spectra in the following format:

$$RT_{\text{start}}-RT_{\text{end}}$$

Figure 77. Spectrum Ranges dialog box with a time range for the averaged spectrum



3. Click **Apply**.

Subtracting Background Spectra

The solvent or other noise can create unwanted background spectra. You can subtract them from the spectrum for one or two ranges. The background subtraction algorithm subtracts an average of the selected scans and redraws the spectrum. The spectrum view header shows the number of subtracted scans. For example, SB: 12 indicates that the algorithm has applied background subtraction (SB) to the spectrum by using 12 scans. To the right of that is subtraction range_1 and range_2.

See these topics:

- [Using the Chromatogram Toolbox to Subtract Background Spectra](#)
- [Using the Spectrum Ranges Dialog Box to Subtract Background Spectra](#)

Using the Chromatogram Toolbox to Subtract Background Spectra

You can use the Chromatogram toolbox to subtract background spectra in a linked spectrum.

❖ To subtract background spectra from the spectrum

1. Select the Chromatogram view of interest and open the Chromatogram toolbox (see [Chromatogram Toolbox](#)).
2. Click the **Background Subtract 1 Range** icon or the **Background Subtract 2 Ranges** icon to background subtract over one or two ranges, respectively.
3. Drag the pointer through the time ranges that define the scans to subtract: once for one range or twice for two ranges. Or, enter the time ranges in the Range 1 and Range 2 boxes.

You can undo the background subtraction by clicking the Background Subtract 1 Range icon or the Background Subtract 2 Ranges icon again.

[Figure 78](#) shows the process of dragging the pointer across range 1.

Figure 78. Background subtraction process for the first time range

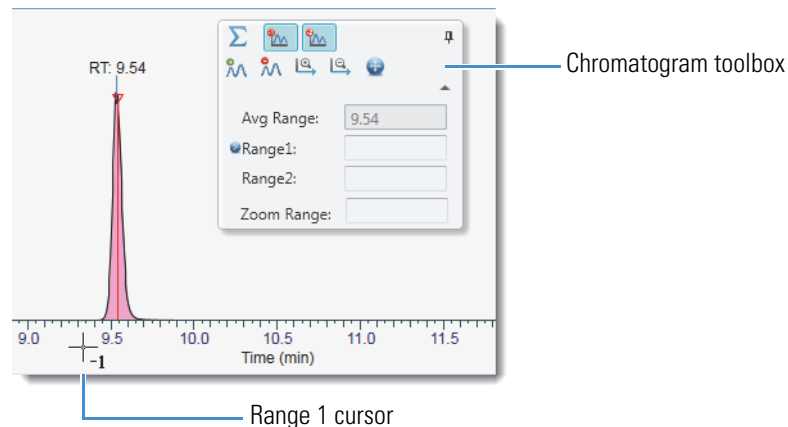


Figure 79 shows the process of dragging the pointer across range 2.

Figure 79. Background subtraction process for range 2

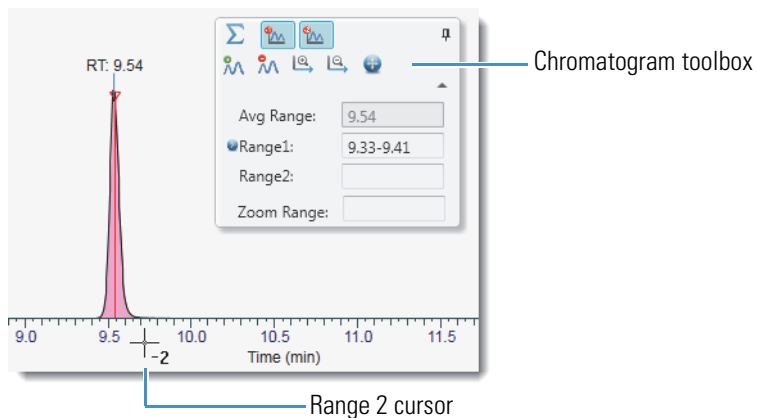
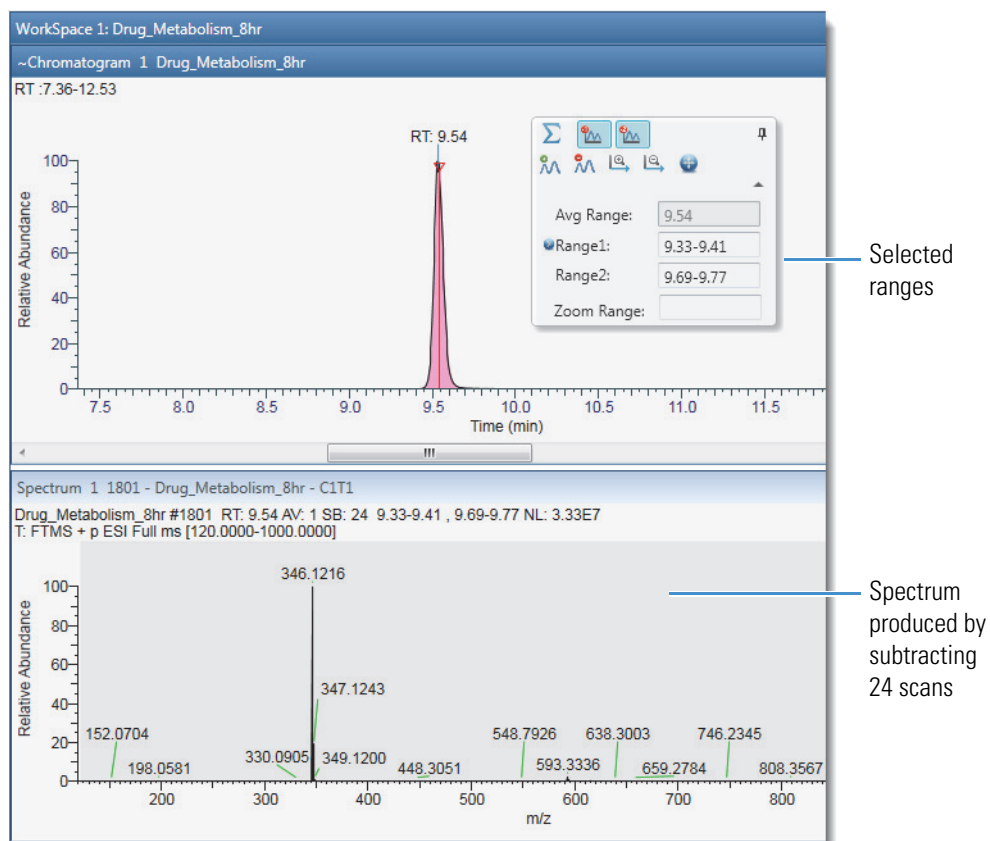


Figure 80 shows the toolbox with the selected background-subtraction ranges and the spectrum produced by subtracting 24 scans.

Figure 80. Spectrum produced by subtracting 24 scans



Using the Spectrum Ranges Dialog Box to Subtract Background Spectra

When a Spectrum view is not linked to a chromatogram, use the Spectrum Ranges dialog box to subtract background spectra.

❖ To subtract background spectra by using the Spectrum Ranges dialog box

1. Select the Spectrum view of interest.
2. In the Spectrum - Workspace Options toolbar, click **Spectrum Ranges**.
The Spectrum Ranges dialog box opens.
3. Use the File Name list and the Retention Time or Scan Number box to define the spectrum.
4. In the Subtract Background list, select **1 Range** or **2 Ranges**.
5. Enter the time range for the background spectra in the Background Range 1 and Background Range 2 boxes as appropriate.
6. Click **Apply**.

Exporting a Selected Spectrum to a New RAW File

Raw data files can contain gigabytes of data. To work with only a selected spectrum, you can export the spectrum to a new RAW file. The new RAW file will contain the selected spectrum and the sample data from the original RAW file and have a much smaller file size.

❖ To export a spectrum to a new RAW file

1. Open a raw data file.
2. Set up the spectrum of interest in a Spectrum or MultiSpectrum view.
The spectrum can be a single scan, a portion of a scan, an averaged spectrum, or a composite spectrum.
3. Click the spectrum plot of interest to select it.
4. In the Spectrum – Workspace Options toolbar, click **Write to .RAW**.
The Export Data dialog box opens. The File name box contains the original file name appended with Scan[*scan number*].
5. Browse to an appropriate storage folder and rename the file as necessary.
6. Click **Save**.

The new RAW file only contains the exported spectrum and the export date, which is listed in the File Header report as the acquisition date and time.

Using Precursor Markers to Open Data-Dependent Scans

You can add these two markers to the tops of the spectrum peaks in precursor scans: Precursor Flag and Nearby Precursors. Double-clicking these markers opens the data-dependent scans in another view.

For information about turning on these markers and using them, see these topics:

- [Using the Precursor Flag Marker](#)
- [Using the Nearby Precursors Marker](#)

Using the Precursor Flag Marker

The application can label the spectrum peaks that triggered data-dependent scans with triangles called precursor markers.

❖ To turn on the precursor flag marker

1. Click a Spectrum view to select it.
2. In the Workspace Options toolbar, click the **Display Options** tab.
3. In the Labels area of the Display Options toolbar, click **Precursor Flag**.

[Figure 81](#) shows a full MS scan with a precursor flag at m/z 195.0877.

Figure 81. Spectrum view with the Precursor Flag marker turned on

Precursor Flag label (enabled)



Precursor marker in the full MS1 scan

❖ **To display the fragmentation scan triggered by the precursor ion**

Double-click a peak's precursor marker to display its product ion spectrum in a separate Spectrum view.

Figure 82 shows a full MS1 scan with one precursor ion at m/z 195.0877 and its data-dependent product ion spectrum in a second Spectrum view.

4 Reviewing Spectral Data

Using Precursor Markers to Open Data-Dependent Scans

Figure 82. Precursor and product ion scans in separate Spectrum views

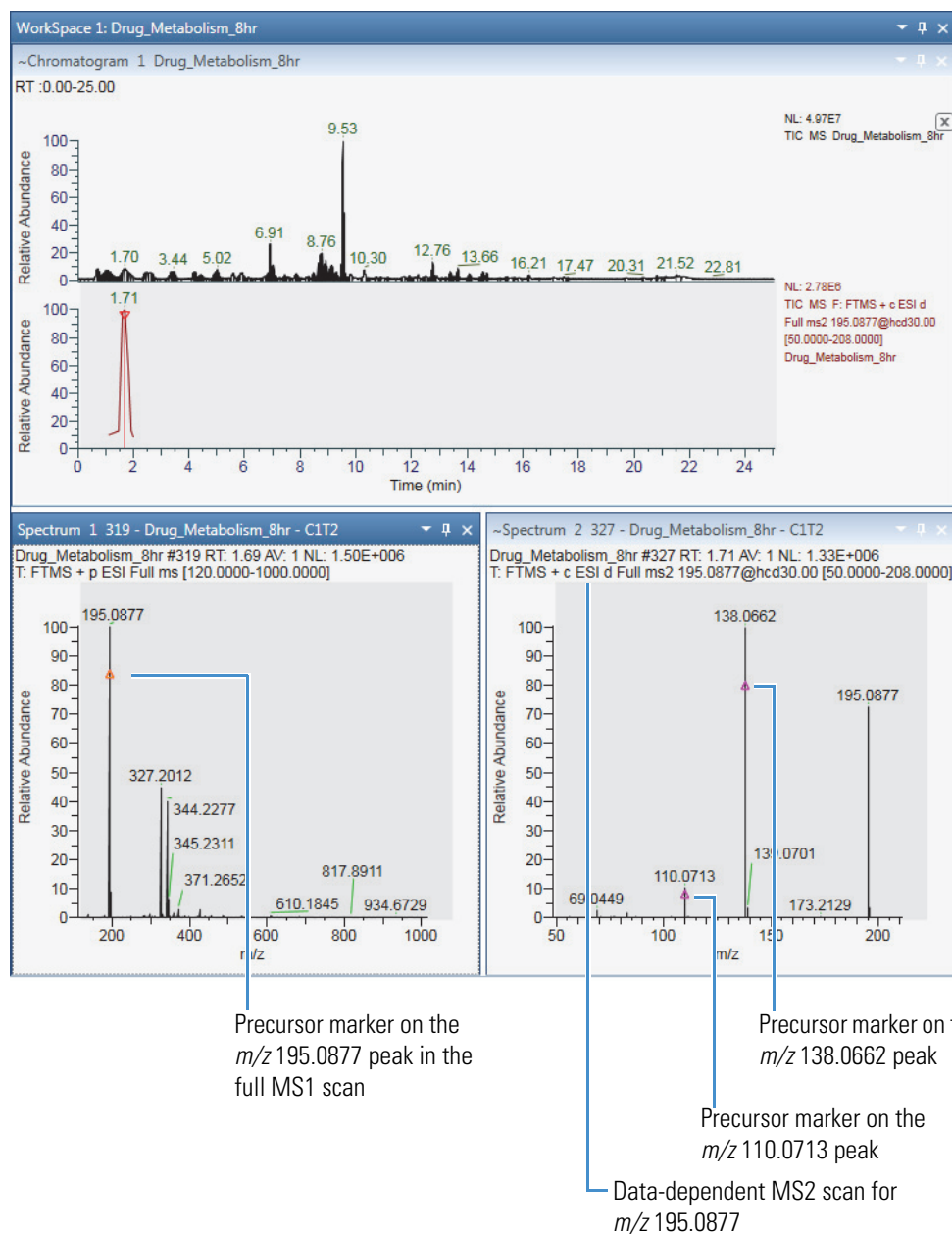
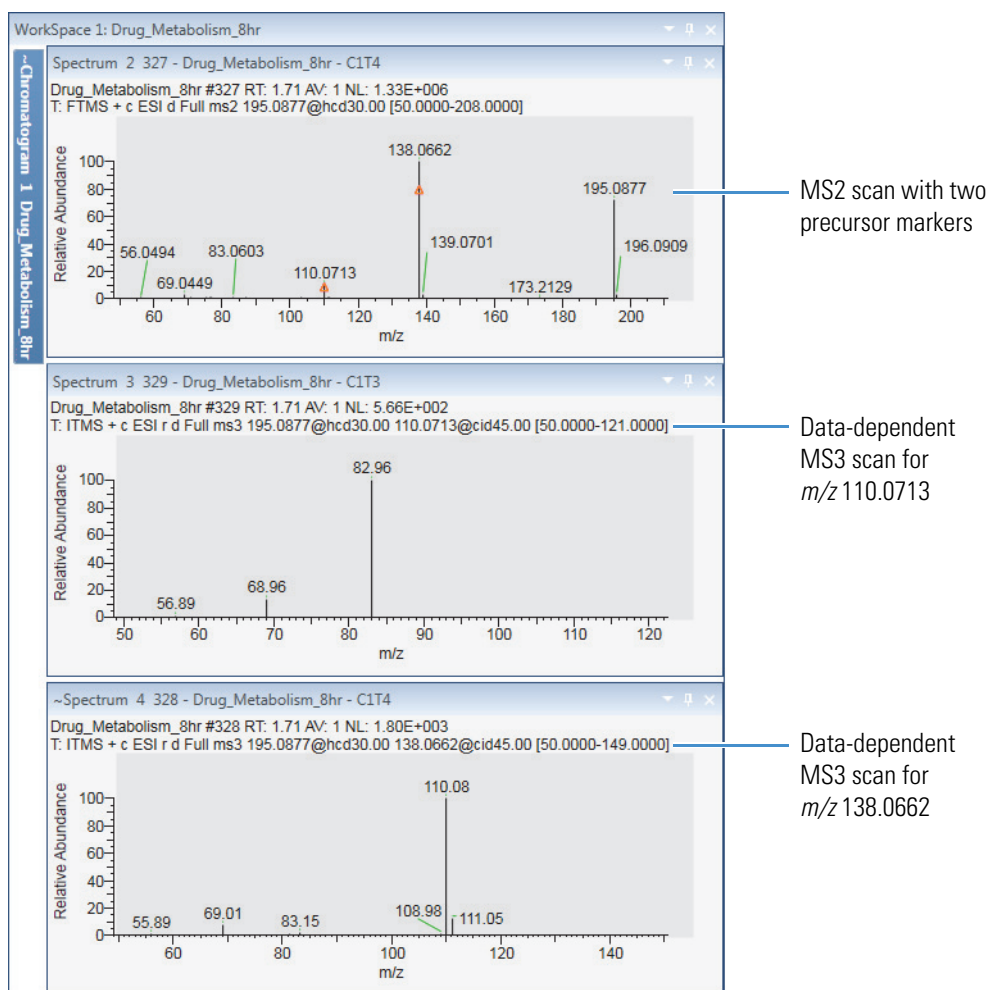


Figure 83 shows the product ion scans for the two marked precursor ions in the data-dependent MS2 scan.

Figure 83. Data-dependent MS2 scan with two precursor peaks and their MS3 product ion scans



Using the Nearby Precursors Marker

Use the Nearby Precursors marker to mark the peaks in the current MS(n-1) scan that the instrument selected for data-dependent MS_n acquisition in the current scan or another scan. The application searches for matching data-dependent scans (by the precursor ion's *m/z* value within the specified mass tolerance) within the specified time range of the current scan.

Double-clicking a marker opens the average spectrum of the data-dependent scans in a separate Spectrum view or individual spectrum plots for each data-dependent scan in a MultiSpectrum view.

You can change the default behavior for this marker on the Workspace Options page.

❖ To set up the display options for the fragmentation scans

1. In the toolbar, click **Default Options**.
The Default Options Configuration dialog box opens.
2. In the left pane, click **Workspace Options**.
3. In the Nearby Precursor area, set up the search and display options for the data-dependent scans:
 - Set the time range to search within the raw data file.

Tip Consider the experiment type when setting the time range. For infusion experiments consider setting the search range to the entire file. For chromatography experiments, consider basing the search range on the chromatographic peak width.

- Set the maximum number of hits (matching data-dependent scans) to return.
 - Set the plot type for the data-dependent scans to an average spectrum in a Spectrum view or individual plots in a MultiSpectrum view.
 - Set the mass tolerance for matching the m/z values of the peaks in the current scan to the m/z values of the precursor ions for the data-dependent scans in the next level of MS/MS activation.
- ### ❖ To turn on the Nearby Precursor marker

1. Click a Spectrum view to select it.
2. In the Workspace Options toolbar, click the **Display Options** tab.
3. In the Labels area of the Display Options toolbar, click **Precursor Flag**.

❖ To display the related data-dependent scans for a precursor ion

Double-click the peak's Nearby Precursor marker (○).

Figure 84 shows a precursor scan with Nearby Precursor markers on the left and the average spectrum for the related data-dependent scans on the right.

Figure 84. Full MS scan with Nearby Precursor markers and an average spectrum of its four related data-dependent scans (matching m/z value within specified RT)

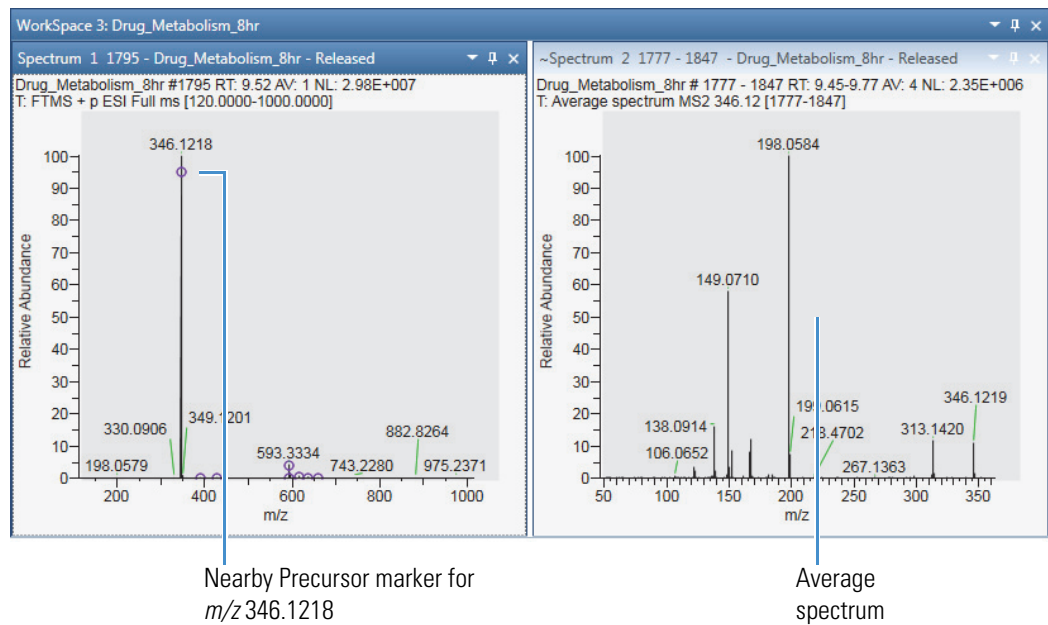
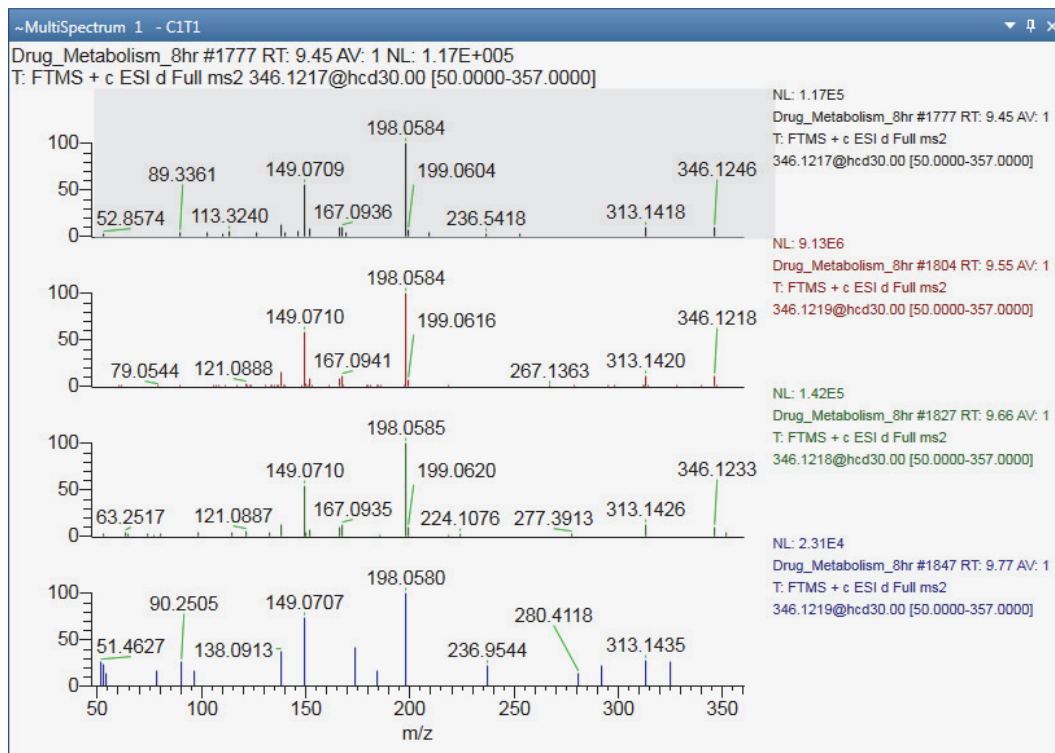


Figure 85 shows the four matching (by precursor m/z value) data-dependent scans that the application found within the specified retention time of scan #1795.

Figure 85. MultiSpectrum view showing the data-dependent scans



Selecting Spectra from an MSn Tree

On the MSn Browser page in the Info Bar view, you can select scans of interest from an MSn spectrum tree. You can also select to display an average spectrum for all the MSn scans of a specific precursor ion or a composite spectrum that includes the mass peaks from all the fragmentation scans that originate from the same precursor ion. For information about the parameter settings and the MSn tree structure, see [MSn Browser Page](#).

Follow these procedures:

- [To display the MSn Browser page](#)
- [To filter the MSn tree by a selected time range](#)
- [To filter the MSn tree by a selected mass range](#)
- [To modify the grouping of the scan data in the MSn tree](#)
- [To expand the entire MSn tree](#)
- [To display all the individual scan items](#)
- [To display an average spectrum](#)
- [To normalize a composite spectrum](#)

❖ To display the MSn Browser page

1. Open a raw data file (see [Opening Raw Data Files or Sequence Files](#)).
2. In the Info bar, click the **MSn Browser** tab.

In the MSn Parameters area, use the Time Range and Mass Range settings in the MSn parameters area to filter the spectra in the MSn Tree area. Use the Mass Tolerance setting to change the grouping of the MSn nodes.

In the MSn Tree area, the branches for each MS2 precursor appear as nested groups ordered by their increasing m/z value. Expanding an MS2 precursor node displays the MS3 precursor nodes when available, expanding the MS3 precursor nodes displays the MS4 precursor nodes when available, and so on.

When an MSn level includes more than one scan, the application displays the average spectrum for the level. If the experimental data includes multiple fragmentation levels, for example MS2 and MS3, the MS3 level includes a composite spectrum made up of the MS2 precursor scan and the MS3 fragmentation scans. An MS4 level node includes a composite spectrum made up of the MS2 and MS3 precursor scans and the MS4 fragmentation scans.

❖ **To filter the MSn tree by a selected time range**

- In the MSn Parameters area, in the Time Range (min) box, type the time range that you want to review.

–or–

- a. In the MSn Parameters area, select the **Track** check box.
- b. In the Chromatogram view, drag the pointer across the *x*-axis range that you want to review.

The selected time range appears in the Time Range (min) box, and the MSn Tree view displays the scans within the selected time range.

❖ **To filter the MSn tree by a selected mass range**

In the Mass Range box, type a beginning mass and an ending mass, separated by a hyphen (*lowest mass-highest mass*). The specified mass range must be within the mass range of the scan data.

The MSn Tree view displays only the scans where the mass of the precursor ion falls within the specified mass range.

❖ **To modify the grouping of the scan data in the MSn tree**

In the Mass Tolerance box, type an *m/z* value from **0.00** to **10.00**.

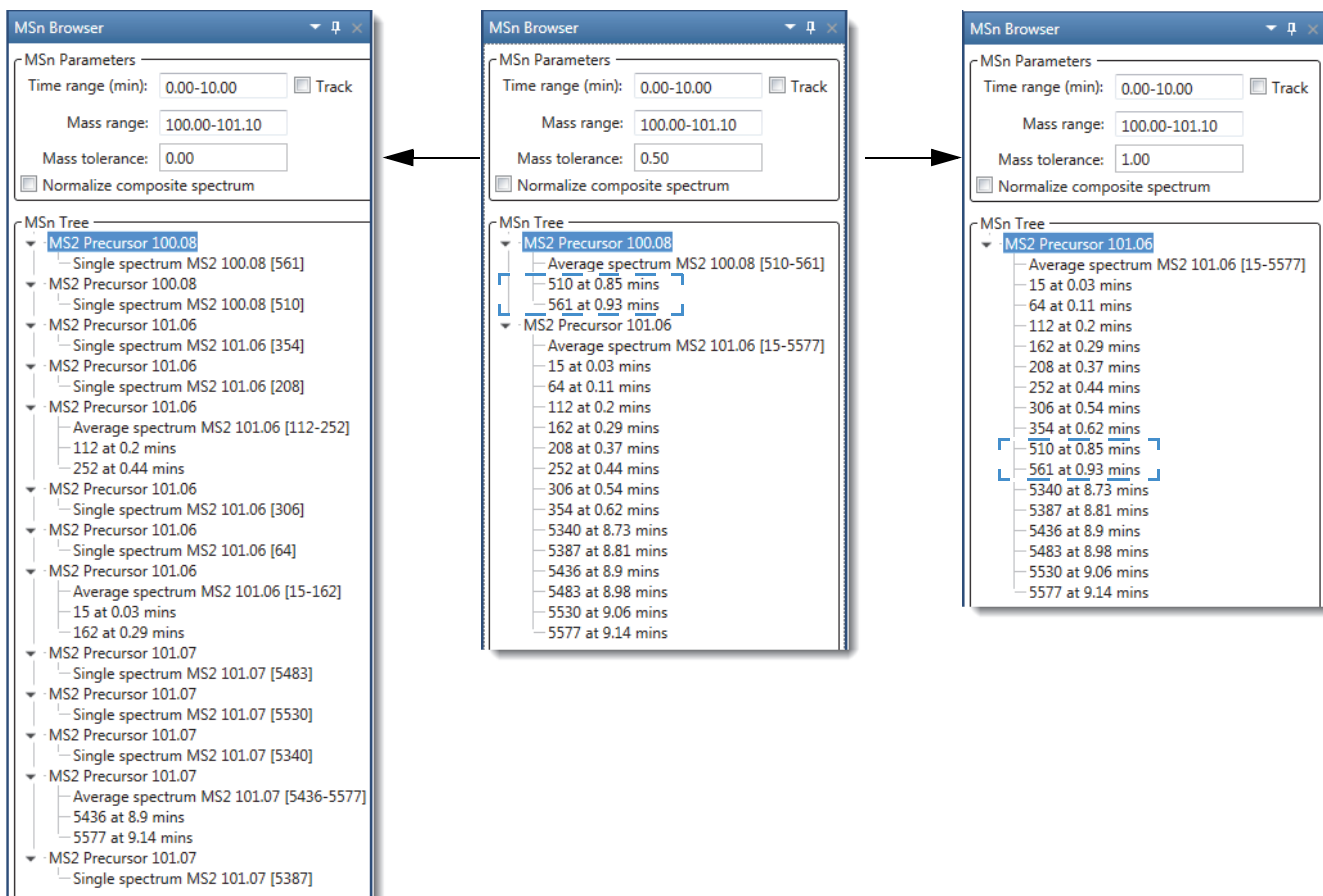
Figure 86 shows the effect of decreasing the mass tolerance from 0.50 to 0.00 and the effect of increasing the mass tolerance from 0.50 to 1.00.

- Decreasing the mass tolerance below that of the experimental data increases the number of MS2 precursor nodes.
- Increasing the mass tolerance from 0.50 to 1.00 combines the scans under the MS2 Precursor 100.08 node with the scans under the MS2 Precursor 101.06 node, as the mass difference between these nodes is less than 1.00.

Figure 86. Effect of changing the mass tolerance on the MSn tree

Decrease from 0.50 to 0.00

Increase from 0.50 to 1.00



❖ **To expand the entire MSn tree**

Right-click the MSn Tree area and choose **Expand List**.

❖ **To display all the individual scan items**

Right-click the MSn Tree area and choose **Include Individual Scans**.

❖ **To display the spectrum for a single scan**

In the MSn Tree area, double-click any of the Single Spectrum *Precursor* items or any of the *Scan Number at Retention Time* items.

A new Spectrum view appears in the Workspace.

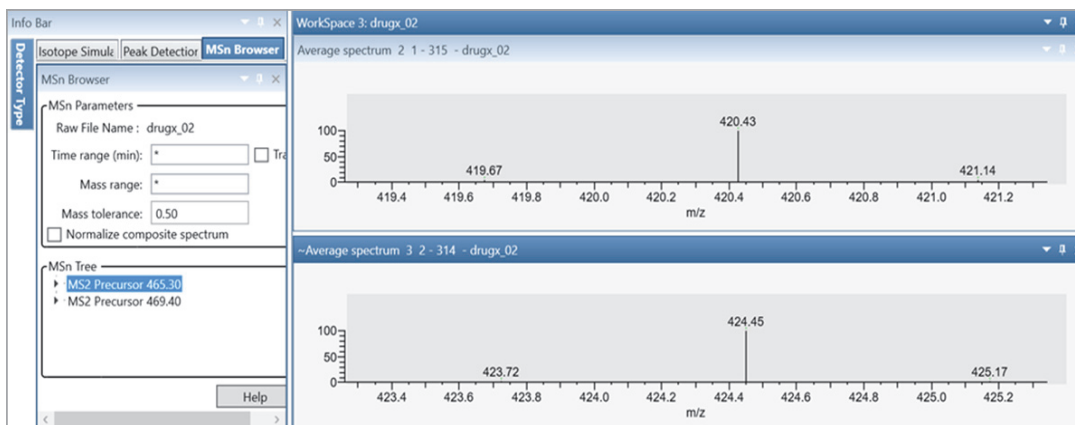
❖ **To display an average spectrum**

In the MSn Tree area, double-click an **Average Spectrum MSn precursor values (scan range)** item.

A new Spectrum view appears in the Workspace. The scan header includes the term Average Spectrum, the number of averaged scans, and the scan number range for the individual scans.

Figure 87 shows average MS2 level spectrum.

Figure 87. Average spectrum of two individual scans



❖ **To display a composite spectrum**

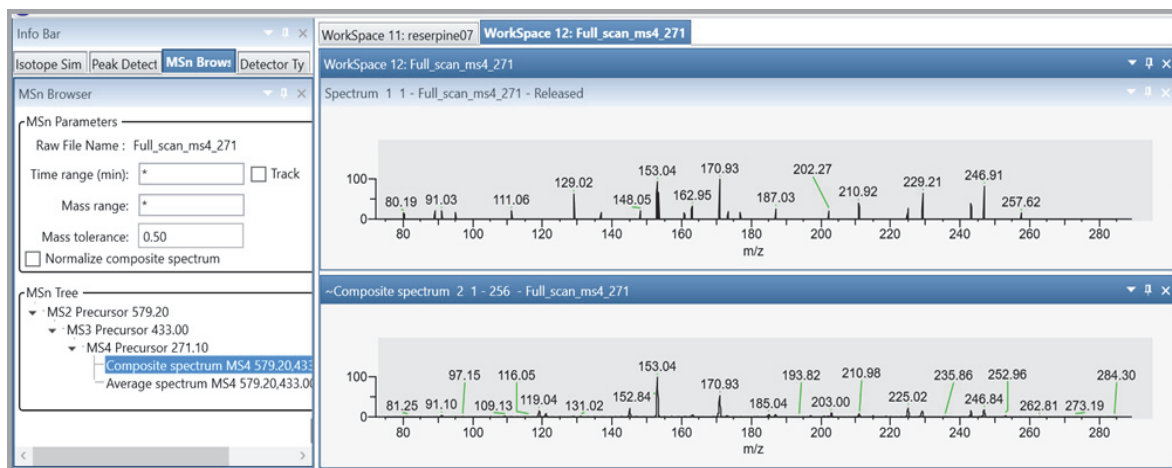
Note The MSn tree includes a composite spectrum for each branch that includes multiple fragmentation levels.

In the MSn Tree area, double-click a **Composite Spectrum MSn precursor values** item.

A new Spectrum view appears in the Workspace. The scan header includes the term Composite Spectrum and the number of averaged scans.

Figure 88 shows a composite spectrum and the two scans (MS2 and MS3 fragmentation levels) that the application combined to create it.

Figure 88. Composite spectrum and individual MS2 and MS3 scans

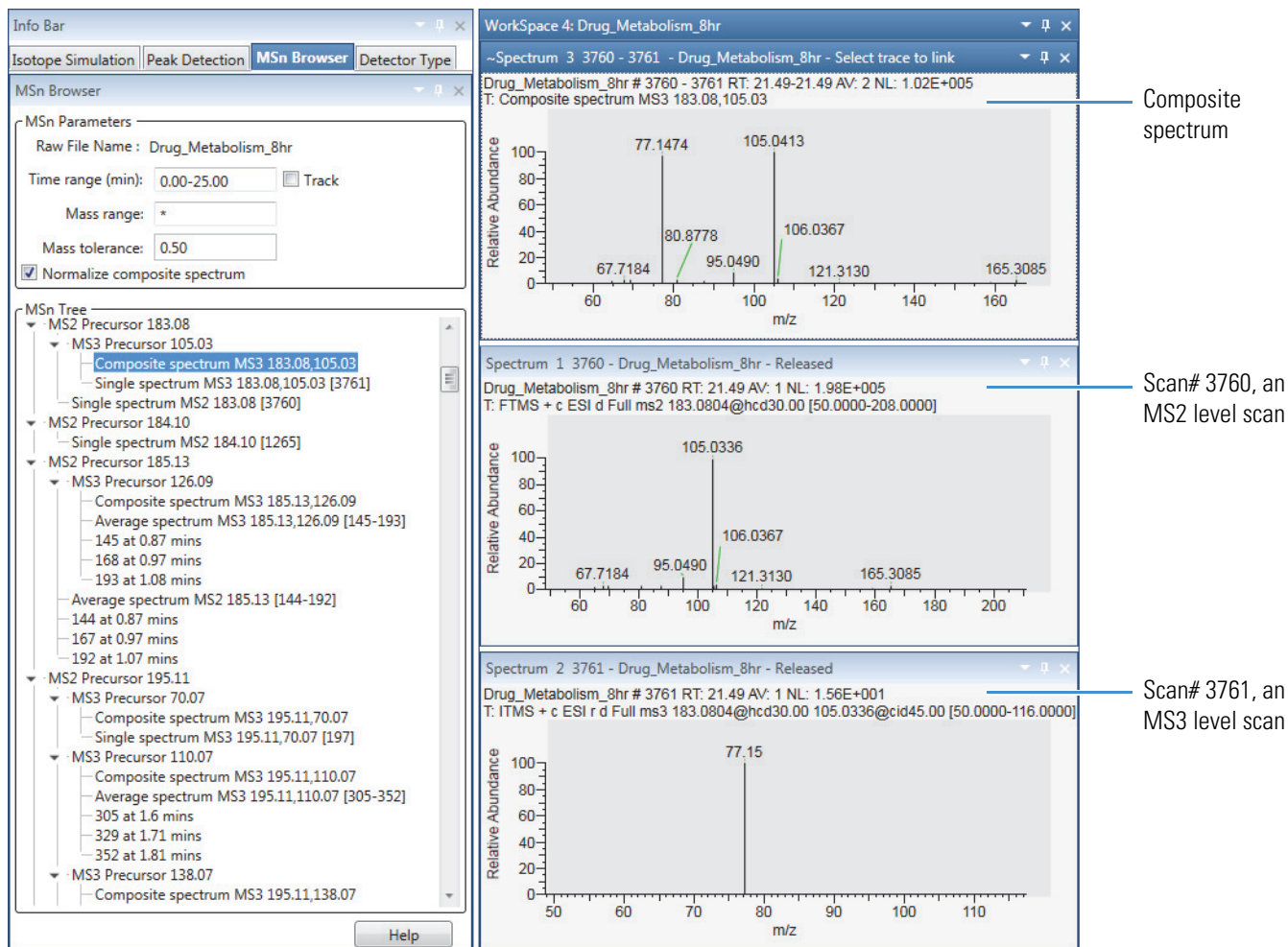


❖ **To normalize a composite spectrum**

Select the **Normalize Composite Spectrum** check box, and then double-click the composite spectrum of interest in the MSn tree.

Figure 89 shows the normalized composite spectrum for scans 3760 and 3761.

Figure 89. Normalized composite spectrum versus the individual MS2 and MS3 scans



Setting Up the Display Options for a Spectrum View

Use the Spectrum – Display Options toolbar to customize the display of the selected Spectrum or Multi Spectrum view.

To set up the display options for the Spectrum view, follow these topics:

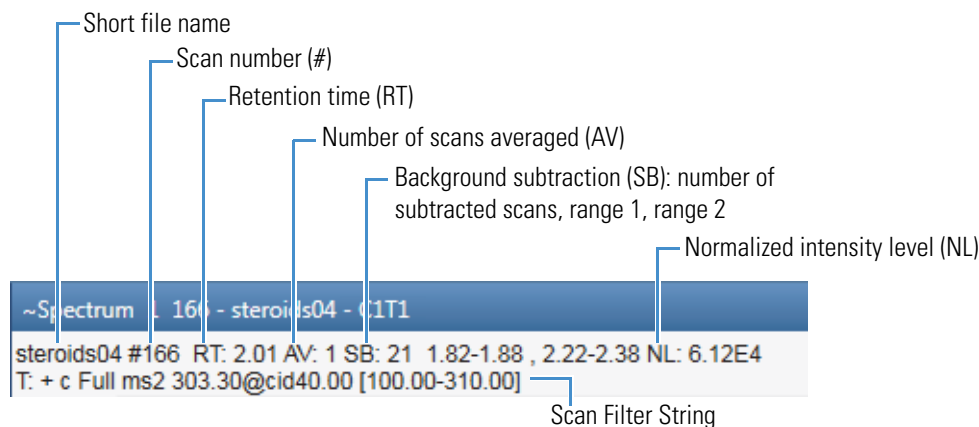
- [Modifying the Scan Header for the Spectrum View](#)
- [Applying Local or Global Normalization in a MultiSpectrum View](#)
- [Labeling Spectrum Peaks](#)
- [Labeling Mass Spectrum Peaks with Chemical Formulas](#)
- [Changing the Y-Axis Scale of a Spectrum View](#)

Modifying the Scan Header for the Spectrum View

By default, the scan header for the Spectrum view displays the following: Short File Name, Scan Number, Retention Time, Average Number of Scans, Background Subtraction Scan Numbers, Normalized Intensity, Polarity, Scan Filter String, and Mass Ranges.

Figure 90 shows a scan header for scan #166 at a retention time of 2.01 min and with background subtraction applied to two ranges.

Figure 90. Default scan header for scan #166 in steroids04.raw with background subtraction



❖ To add or delete information from the scan header for a Spectrum view

1. Open a raw data file.
2. If you are using a custom layout that does not include a Spectrum or Multi Spectrum view, in the Workspace Options toolbar, click **Spectrum** or **Multi Spectrum**.
3. Click the Spectrum (or MultiSpectrum) view of interest.

Note The selections that you make in the Scan Header dialog box affect only the selected Spectrum view (or Multi Spectrum view).

4. In the Spectrum – Display Options toolbar, click **Scan Header**.

The Scan Header dialog box opens ([Scan Headers and Scan Header Abbreviations](#)).

5. Do the following:
 - To display a field, select its corresponding check box.
 - To hide a field, clear its corresponding check box.
6. Click **OK**.

Applying Local or Global Normalization in a MultiSpectrum View

In a MultiSpectrum view, you can normalize the mass spectra to the most intense peak in all the spectra, or you can normalize the mass spectral peaks in each spectrum to the most intense peak in the spectrum.

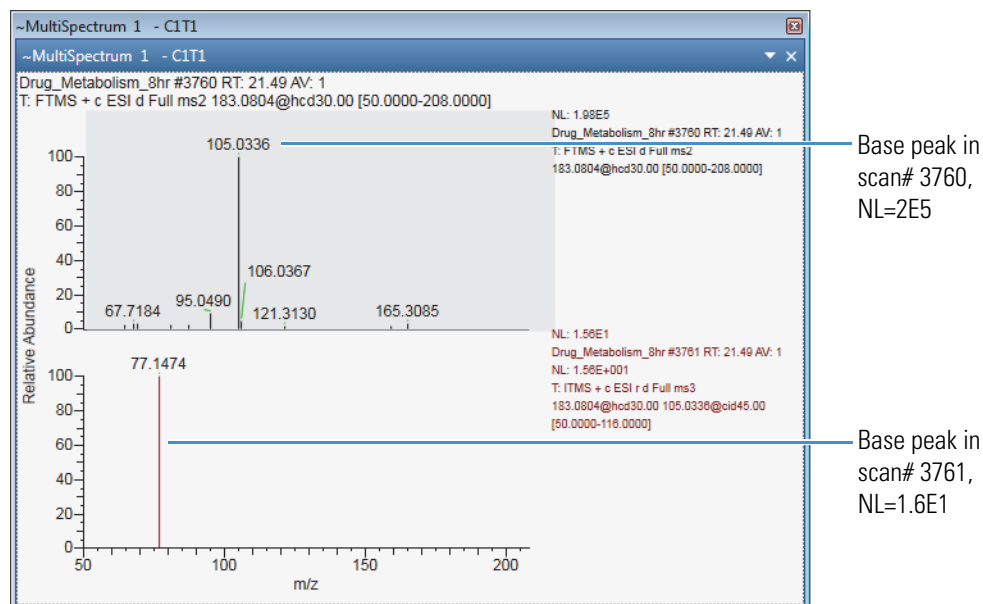
For information about displaying spectra in a MultiSpectrum view, see [Creating a MultiSpectrum View and Changing Its Spectrum Plots](#).

❖ To normalize each spectrum separately

In the Normalization area of the Multi Spectrum – Display Options toolbar, click **Local**.

Figure 91 shows a set of locally normalized spectrum plots. Each spectrum is normalized to the most intense peak in the spectrum.

Figure 91. Local normalization



❖ To normalize the spectra against the most intense peak across the spectra

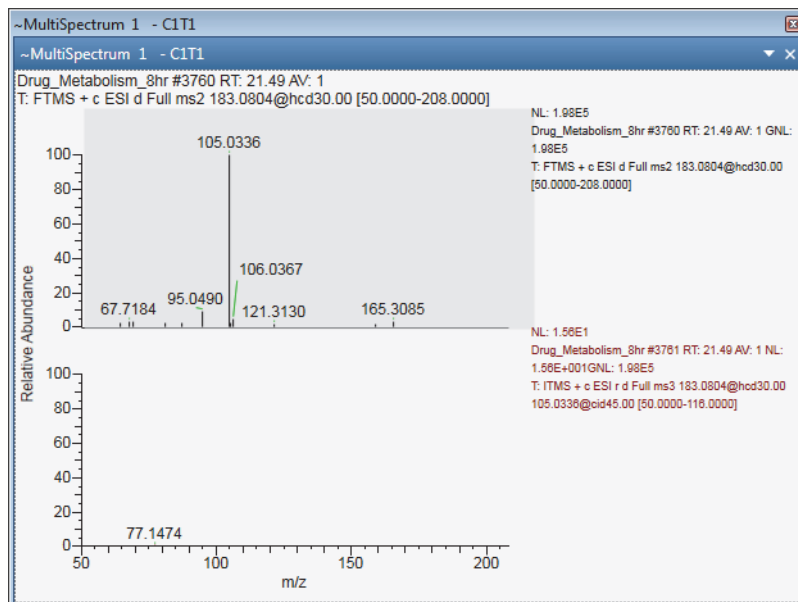
In the Normalization area of the Multi Spectrum – Display Options toolbar, click **Global**.

Figure 92 shows the effect of global normalization where both spectrum plots are normalized to the largest peak across the plots. The global normalization (GNL) is equal to 2E5, and the low intensity peak in scan# 3716 is barely visible.

4 Reviewing Spectral Data

Setting Up the Display Options for a Spectrum View

Figure 92. Global normalization



Labeling Spectrum Peaks

Use the Labels area of the Spectrum – Display Options toolbar to add or remove labels that appear above the spectrum peaks. By default, the application adds the mass-to-charge label to each mass spectrum peak.

Note The Precursor Flag and Nearby Precursor buttons do not add text labels to the spectrum peaks; these buttons add markers to the precursor scans. For instructions on how to use these markers, see [Using Precursor Markers to Open Data-Dependent Scans](#).

❖ To add or remove the labels for the mass spectrum peaks

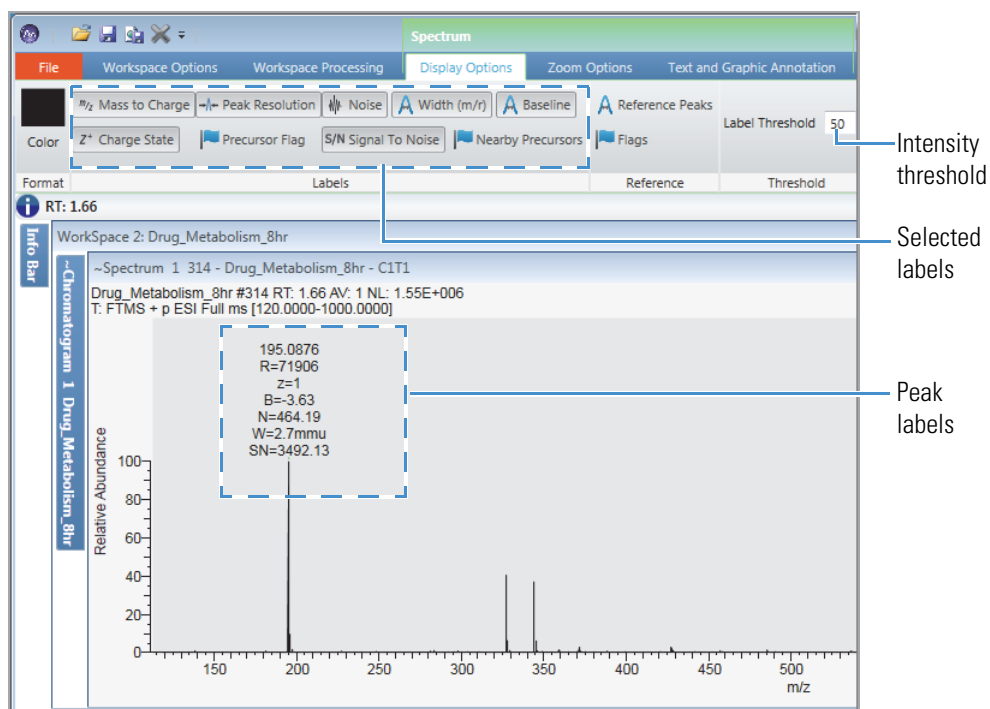
1. Open a raw data file and display a spectrum in a Spectrum view (or MultiSpectrum view).
2. Click the Spectrum (or MultiSpectrum) view of interest.
3. In the Labels area of the Display Options toolbar, click the labels that you want to display.

Note With the default value set to 0 in the Label Threshold box, the application labels all the mass peaks.

4. (Optional) To avoid labeling low-intensity mass peaks, type a higher relative intensity value in the Label Threshold box.

[Figure 93](#) shows a mass spectrum with several labels and a labeling threshold of 50% relative intensity.

Figure 93. Mass spectrum with multiple labels



Labeling Mass Spectrum Peaks with Chemical Formulas

You can add formula labels to the peaks in your mass spectra.

Note When calculating the elemental compositions of all the mass peaks in a spectrum, the application uses the settings on the Default Elemental Composition page. When you change the settings on this page, the new settings take effect after you restart the application.

Tip When determining the chemical formula for each mass peak, the application assumes that the current mass peak is an A0 peak; that is, it assumes that the peak corresponds to a monoisotopic ion, and the peaks at higher m/z values are part of the ion's isotope pattern. After determining the best matching formula for the current mass peak (by using the accurate mass and isotope pattern), the application analyzes the next mass peak, again assuming that the peak is an A0 peak.

Use the Isotope Simulation page to calculate the theoretical isotope patterns for the displayed formulas.

❖ To add formula labels to the mass peaks in a spectrum plot

1. Click a Spectrum view to select it.

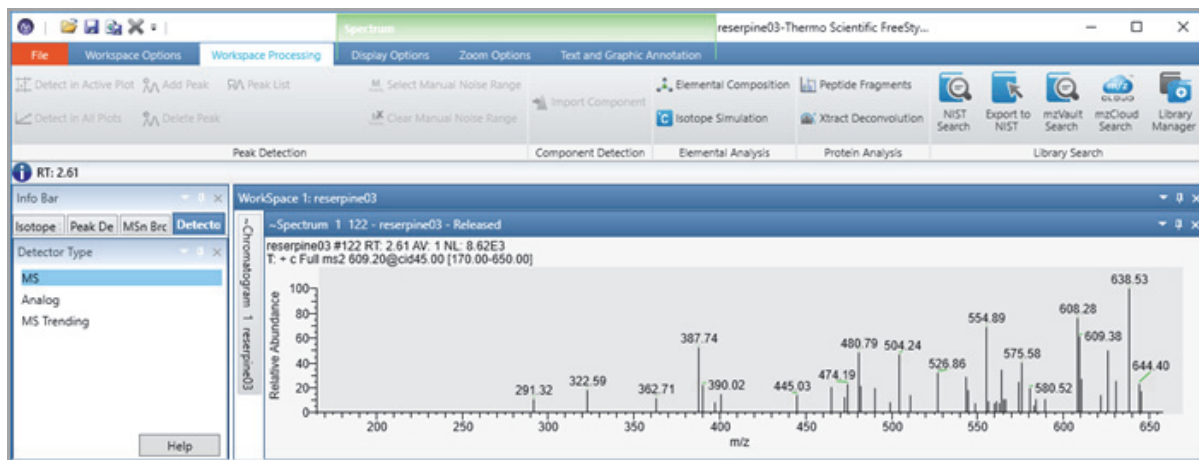
4 Reviewing Spectral Data

Setting Up the Display Options for a Spectrum View

2. Open the Spectrum – Workspace Processing toolbar.

Figure 94 shows the Spectrum – Workspace Processing toolbar in the Elemental Analysis area.

Figure 94. Workspace Processing toolbar for the Spectrum view



3. (Optional) To ignore low-intensity peaks, type a higher intensity threshold in the Label Threshold box.
4. In the Elemental Analysis area of the toolbar, click **Elemental Composition**.

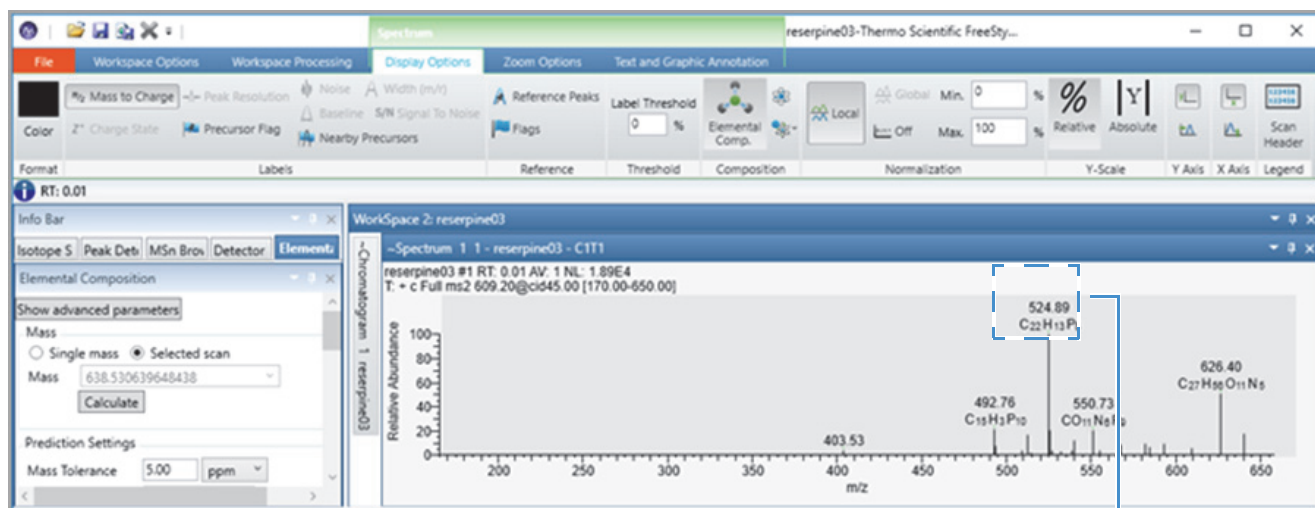
The Elemental Composition page appears in the Info Bar.

Select the **Selected Scan** check box, and then click **Calculate**.

The Theoretical Mass (Theo.) and Delta Mass buttons become available in the Composition area of the Display Options toolbar.

Figure 95 shows the parameters on the Elemental Composition page and the chemical formula labels in the Spectrum view. You can use the Elemental Composition page to find the best matching formulas for a specific mass peak.

Figure 95. Mass peaks with chemical formula labels in the Spectrum view



Experimental m/z value and calculated chemical formula for the ion

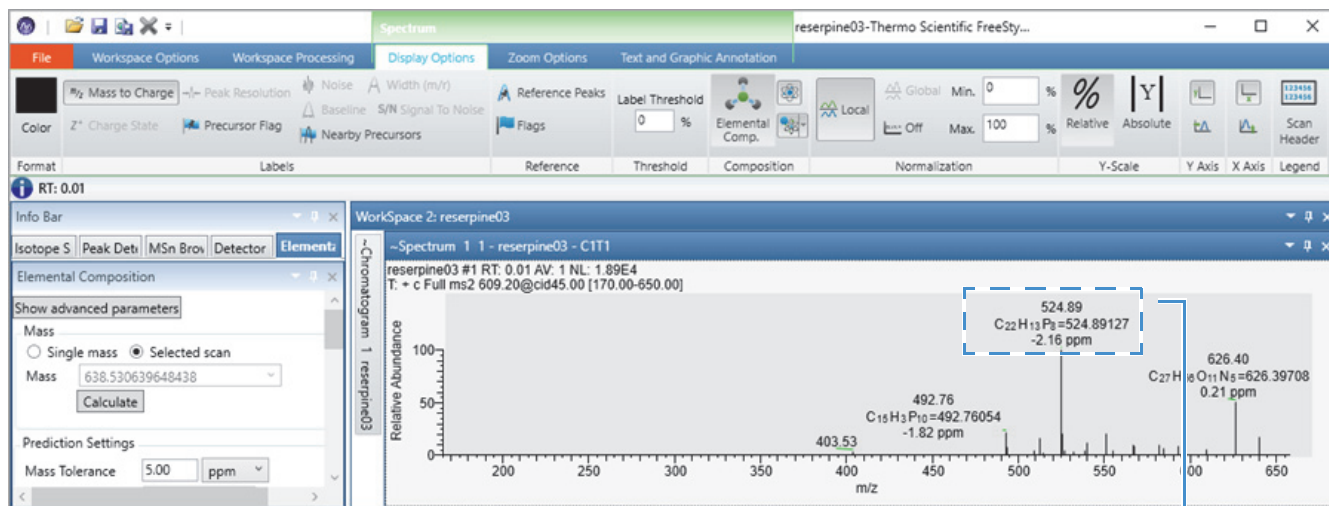
- (Optional) To label the peaks with the theoretical mass for the displayed formula and the mass difference between the experimental m/z value and the theoretical mass, click **Theo.** and **Delta Mass**, respectively.

Figure 96 shows the mass peaks labels, from top to bottom: the experimental m/z value, the formula and calculated m/z value of the ion, and the mass difference between the experimental m/z value and the calculated m/z value.

4 Reviewing Spectral Data

Setting Up the Display Options for a Spectrum View

Figure 96. Mass peaks with formula, theoretical mass, and delta mass labels



Experimental m/z value, calculated chemical formula for the ion, calculated m/z value for the formula, and mass difference between the experimental m/z value and the calculated m/z value

Changing the Y-Axis Scale of a Spectrum View

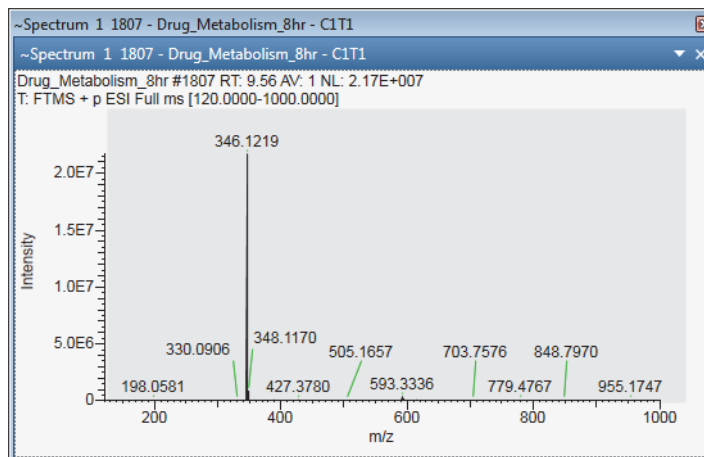
You can select the relative intensity scale or the absolute intensity scale for the y axis of a spectrum plot.

With the factory default template, a **Spectrum View** opens with the y -axis scale set to relative intensity and a y -axis label of Relative Abundance.

❖ To change the y -axis scale from relative intensity to absolute intensity

1. Click a Spectrum view to select it.
2. Click the **Display Options** toolbar tab.
3. In the Y-Scale area of the toolbar, click **10^x Absolute**.

The y -axis maximum changes from 100% to the normalization (NL) value (or slightly less than the NL value) for the scan, and the y -axis label changes from Relative Abundance to Intensity (**Figure 97**).

Figure 97. Spectrum plot of (absolute) Intensity versus m/z value

Changing the Zoom Level in a Spectrum View or MultiSpectrum View

You can zoom in on or out of a region of a spectrum by using the Spectrum – Zoom Options toolbar or the MultiSpectrum Zoom Options toolbar, the Spectrum toolbox, or the mouse pointer.


Follow these procedures:

- To incrementally zoom in and out on the x axis
- To zoom in on a specific section of the x axis
- To incrementally zoom in and out on the y axis
- To reset the zoom level of the x and y axes

❖ To incrementally zoom in and out on the x axis

1. Select the view.
2. To zoom in, do the following:
 - Open the Zoom Options toolbar and click **Zoom In X**.


–or–

- Open the Spectrum toolbox and click the **Zoom In** icon, .

Each click zooms in on a 50% smaller section of the x axis—that is, it decreases the displayed range by a factor of 2.


3. To zoom out, do the following:
 - In the Zoom Options toolbar, click **Zoom Out X**.

–or–

- Open the Spectrum toolbox and click the **Zoom Out** icon, . Each click increases the displayed m/z range by 200%.
 - ❖ **To zoom in on a specific section of the x axis**

Drag the pointer horizontally across the specific section of the x axis.
 - ❖ **To incrementally zoom in and out on the y axis**
 1. Select the view.
 2. To zoom in, open the Zoom Options toolbar and click **Zoom In Y**.

Each click zooms in on a 50% smaller section of the y axis. For example, in a Spectrum view where the y axis is set to relative abundance, clicking once zooms in on the 0 to 50% range, clicking twice zooms in on the 0 to 25% range, and so on.
 3. To incrementally undo the zoom level, click **Zoom Out Y**.

Each click zooms out to a 200% larger section of the y axis.
 - ❖ **To reset the zoom level of the x and y axes**
 - Select the view, open the Zoom Options toolbar, and click **Reset**.
 - Right-click the view and choose **Reset Scaling**.
- or–
- Select the view, open the Spectrum toolbox, and click the **Reset Zoom** icon, .

Spectra-Specific Toolbars

Use the spectra-specific toolbars as follows:

- [Spectrum – Workspace Processing Toolbar](#): For determining the elemental composition of an ion; simulating isotope distributions; Xtract deconvolution; and NIST library, mzVault library, and mzCloud.org searches.
- [Spectrum – Display Options Toolbar](#): For customizing the Spectrum view
- [MultiSpectrum – Display Options Toolbar](#): For customizing the MultiSpectrum view
- [Spectrum List – Display Options Toolbar](#): For formatting the spectrum list

These two toolbars are specific to the Spectrum, MultiSpectrum, and Chromatogram views:

- [Zoom Options Toolbar](#): For adjusting the display of the spectra (and chromatograms).
- [Text and Graphic Annotation Toolbar](#): For annotating spectra (and chromatograms) with text, lines, boxes, and symbols.

Spectrum – Workspace Processing Toolbar

Use the buttons in the Elemental Analysis and Library Search areas of the Workspace Processing toolbar to analyze spectra.

❖ To display the Spectrum Workspace Processing toolbar

1. Click a Spectrum view or a MultiSpectrum view to select it.
2. Click the **Workspace Processing** toolbar tab.

Figure 98 shows the Elemental Analysis and Library Search areas in the Workspace Processing toolbar, and Table 27 describes toolbar buttons.

Figure 98. Elemental Analysis and Library Search areas of the Workspace Processing toolbar

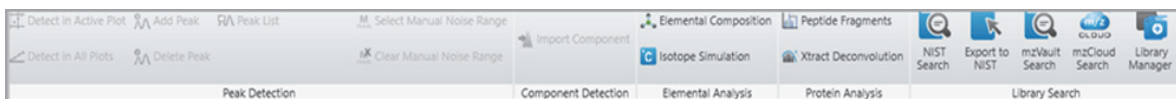


Table 27. Spectra-specific Workspace Processing buttons (Sheet 1 of 2)

Button	Description
Elemental Analysis	
Elemental Composition	Displays the Elemental Composition page, where you calculate the best matching chemical formulas for a mass in the spectrum. See Chapter 7, “Determining the Elemental Composition of Ions.”
Isotope Simulation	Displays the Isotope Simulation page of the Info Bar, where you create a simulated isotopic distribution spectrum of a chemical formula.
Protein Analysis	
Peptide Fragments	Displays the Peptide Fragments page of the Info Bar, where you set parameters to annotate the active spectrum using the CID or ETD activation types. See
Xtract	Displays the Xtract page of the Info Bar, where you set parameters for the Xtract algorithm. See Chapter 13, “Deconvolving and Deisotoping Spectra with the Xtract Algorithm.”
Library Search	
NIST Search	Opens the Library Search page in the Info Bar, where you run a NIST library search on the selected spectrum and displays the results in the NIST Search Results view. You can modify the library search parameters on the Modifying a NIST Search from the NIST Search Page of the Info Bar. See Performing a Local NIST or mzVault Library Search .

Table 27. Spectra-specific Workspace Processing buttons (Sheet 2 of 2)

Button	Description
Export to NIST	Exports a spectrum to the NIST application for a search. The NIST application opens and displays the search results (see Exporting a Mass Spectrum to the NIST MS Search Application).
mzVault Search	Opens the mzVault Search page in the Info Bar, runs an mzVault library search on the selected spectrum, and displays the results in the mzVault Search Results view.
mzCloud	Uploads a spectrum to mzCloud.org for a search. The website opens and displays the search parameters. Select the appropriate settings and click OK . The website displays the results.
Library Manager	Opens the Thermo Library Manager dialog box, where you select or create NIST libraries. See Managing Libraries .

Spectrum – Display Options Toolbar

Use the buttons in the Spectrum – Display Options toolbar to customize the Spectrum view (see [Setting Up the Display Options for a Spectrum View](#)). You can label the spectral peaks, determine the elemental composition for a peak, and specify how the application normalizes the spectrum.

❖ To display the Spectrum Display Options toolbar

1. Click a Spectrum view to select it.
2. Click the **Display Options** toolbar tab.

Figure 99 shows the Spectrum Display Options toolbar.

Figure 99. Spectrum Display Options toolbar (in two parts, left and right side)

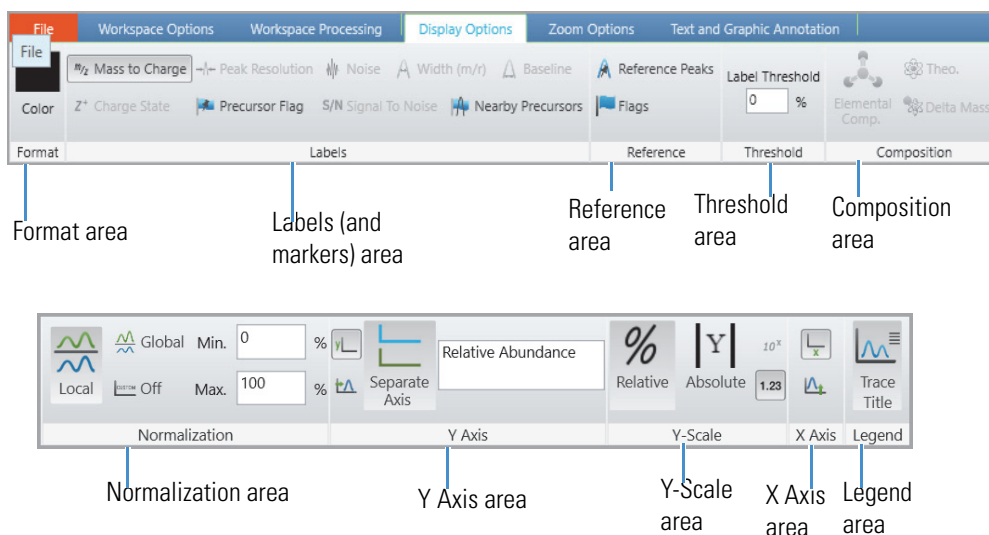


Table 28 describes the buttons in the Spectrum – Display Options toolbar.

Table 28. Spectrum – Display Options toolbar buttons (Sheet 1 of 4)

Button	Description
Color	
Color	Displays the color palette, where you select the color of the spectrum.
Labels	
Use the buttons in the Labels area to annotate the mass spectrum.	
Note The following buttons are unavailable when the raw data file does not include the requested information: Z ⁺ Charge State, Peak Resolution, Noise, Baseline, Width, and S/N Signal To Noise.	
Mass to Charge	Adds a mass-to-charge ratio label to each peak in the spectrum.
Z ⁺ Charge State	Adds a charge state label to each peak in the spectrum.
Peak Resolution	Adds a peak resolution label to each peak in the spectrum.
Precursor Flag	Adds one or more pink triangles to a peak if it triggered data-dependent scans.
Noise	Adds a noise magnitude label to each peak in the spectrum.
Baseline	Adds a baseline label to each peak in the spectrum.
Width (m/r)	Adds a width label to each peak in the spectrum. The width equals the mass, in mmu units, divided by the peak resolution.

Table 28. Spectrum – Display Options toolbar buttons (Sheet 2 of 4)

Button	Description
Signal To Noise	Adds a signal-to-noise label (intensity/noise) to each peak in the spectrum.
Nearby Precursors	Adds markers (○) to the tops of the spectrum peaks in scans that have related data-dependent scans. Depending on the configuration setting for the Nearby Precursor Plot Type parameter, clicking a peak's marker opens a Spectrum view with an average spectrum for all the related data-dependent scans or a MultiSpectrum view with a separate plot for each related data-dependent scan (within the specified RT window).
Reference	
Reference Peaks	Adds an R label to peaks that correspond to a reference compound used for an internal recalibration of a scan.
Flags	Displays one of the following letters or symbols above flagged peaks: <ul style="list-style-type: none"> • S—Saturated peaks are peaks with a signal too large to measure—that is, the signal was so high that it was outside the dynamic range of the detector, causing saturation. • R—Reference peaks are peaks from a reference compound used for an internal recalibration of a scan. • L—Lock peaks are local references used to calculate the accurate mass of nearby peaks. • E—Exception peaks are peaks from a reference compound that are not used for recalibration. These are typically small isotopes or fragments of the main references. • #—Mathematically modified peaks are peaks where the peak mass was recalculated by the instrument, usually due to a calibration process. • M—Merged peaks are peaks where the centroider combined two nearby peaks. • F—Fragmented peaks are peaks separated into multiple peaks by the centroiding activity.

Table 28. Spectrum – Display Options toolbar buttons (Sheet 3 of 4)

Button	Description
Threshold	
Label Threshold	Sets the percentage of the base peak so that the application labels only the peaks above that percentage. For example, if the base peak is 100 percent and the label threshold setting is 50 percent, the application labels only the peaks in the Spectrum view or MultiSpectrum view that are at or above 50 percent. Range: 0–100; default: 0%
Composition	
Use the buttons in the Composition area to determine the elemental composition of the ions associated with the peaks in the mass spectrum. See Chapter 7, “Determining the Elemental Composition of Ions.”	
Elemental Composition	Enables the formula labels on all the peaks in the spectrum.
Theoretical	Labels the peaks in the mass spectrum with the theoretical mass-to-charge ratios of the best matched ions.
Delta Mass	Labels the peaks in the mass spectrum with the differences between the experimentally determined mass-to-charge ratios and the theoretical mass-to-charge ratios. The unit options are amu (atomic mass units), mmu (millimass units), and ppm (parts-per-million).
Normalization	
Use the parameters in the Normalization area to specify how the application normalizes the mass spectrum.	
Local	Scales the y-axis range as a percentage.
Global	Normalizes the spectra traces so that the most intense peak of all the spectra is 100 percent. Available for a MultiSpectrum view.
Off	Scales the spectra according to the user-specified minimum and maximum absolute intensity (counts).
Min.	Displays the minimum of the y axis. Enter a value in the box to change the minimum. For local normalization, the value is a percentage. For normalization set to Off, the value is an intensity.
Max.	Displays the maximum of the y axis. Enter a value in the box to change the maximum. For local normalization, the value is a percentage. For normalization set to Off, the value is an intensity.

Table 28. Spectrum – Display Options toolbar buttons (Sheet 4 of 4)

Button	Description
Y-Scale	
Use the parameters in the Y-Scale area to specify how the application scales the <i>y</i> axis.	
Absolute	Plots the spectrum with the intensity expressed in counts.
Relative	Scales the spectrum so that the intensity of the most intense peak in the spectrum is 100 percent.
Y Axis	
Use the parameters in the Y Axis area to specify how the application labels the <i>y</i> axis.	
Show Labels	Shows or hides the <i>y</i> -axis label.
Offset Axis	Sets the location of the displayed plot a specified distance from the <i>y</i> axis. The <i>y</i> -axis offset moves the <i>x</i> axis slightly to the right of the <i>y</i> axis so that you can see plot details at low <i>x</i> -axis values.
X Axis	
Use the parameters in the X Axis area to specify how the application labels the <i>x</i> axis.	
Show Labels	Shows or hides the <i>x</i> -axis label.
Offset Axis	Sets the location for the displayed plot a specified distance from the <i>x</i> axis. The <i>x</i> -axis offset moves the <i>y</i> axis up slightly so that you can see plot details at low <i>y</i> -axis values.
Legend	
Scan Header	Opens the Scan Header dialog box, where you specify what scan header information the Spectrum view displays. For additional information, see Scan Headers and Scan Header Abbreviations .

MultiSpectrum – Display Options Toolbar

Use the buttons in the Spectrum – Display Options toolbar to customize the MultiSpectrum view.

Figure 100. MultiSpectrum – Display Options toolbar (left and right)

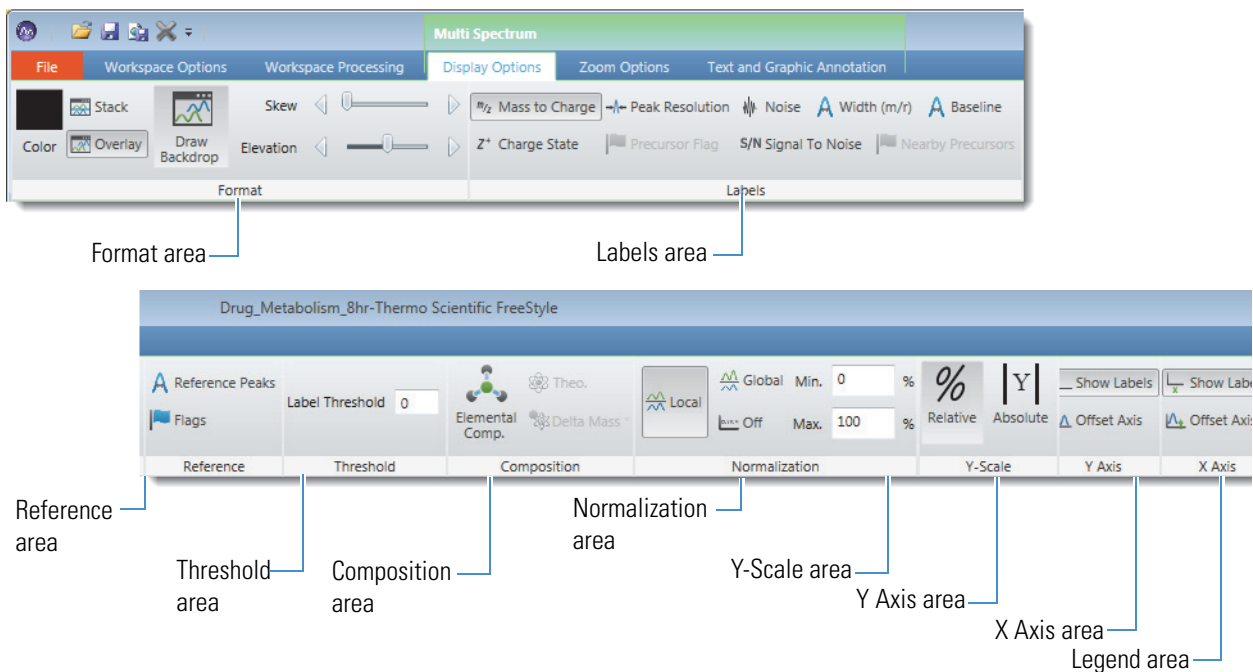


Table 29 describes the additional formatting options for the in the MultiSpectrum – Display Options toolbar. For information about the other toolbar areas, see [Spectrum – Display Options Toolbar](#).

Table 29. Spectrum Display Options toolbar buttons in the Format area (Sheet 1 of 2)

Button	Description
Format	
Color	Displays the color palette, where you select the color of the spectrum.
Stack	Stacks the spectrum traces vertically.
Overlay	Overlays the spectrum traces vertically with an optional horizontal skew (time offset).
Draw Backdrop	Draws a backdrop for overlaid plots.

Table 29. Spectrum Display Options toolbar buttons in the Format area (Sheet 2 of 2)

Button	Description
Skew	Sets the skew angle (time offset) to a value from 0–45 degrees for an overlay arrangement of spectrum traces. To set the skew, drag the Skew slider.
Elevation	Sets the vertical spacing for an overlay arrangement of spectrum traces. To set the vertical spacing, drag the Elevation slider. Move the Elevation slider to the farthest left to overlay the plots on top of each other.

Spectra-Specific Views

The FreeStyle window has these spectra-specific views:

- **Spectrum view:** Displays a spectrum corresponding to the retention time or scan number that you select in the Chromatogram view.
- **MultiSpectrum view:** Displays multiple spectra, but only the active spectrum updates corresponding to the retention time or scan number that you select in the Chromatogram view.
- **Spectrum List view:** Lists in tabular form the positions, intensities, and relative intensities of the peaks in the Spectrum view.

For information about the shortcut menu for the Spectrum and MultiSpectrum views, see [Table 31](#).

For more information, see these topics:

- [Spectrum View](#)
- [MultiSpectrum View](#)
- [Spectrum List View](#) (in the Viewing Experiment and Instrument Information chapter)

Spectrum View

The Spectrum view ([Figure 101](#)) displays a spectrum for the retention time that you either select in a chromatogram trace or specify in the Scan Ranges dialog box.

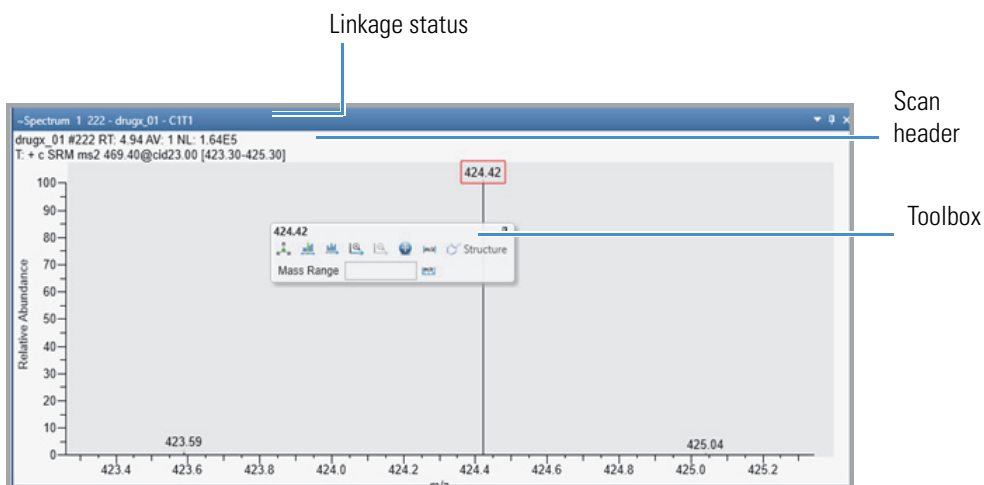
For information about displaying and reviewing the spectral data in a raw data file, see the list of topics at the beginning of this chapter, [“Reviewing Spectral Data.”](#)

❖ **To add a Spectrum view to the Workspace**

1. Click the **Workspace Options** toolbar tab.
2. Click **Spectrum**.

Figure 101 shows an example of a selected Spectrum view and the Spectrum toolbox. The view's title bar displays the scan number of the displayed spectrum, the file name of the raw data file, and the view's link status. CIT1 specifies that the view is linked to trace 1 in Chromatogram view 1.

Figure 101. Spectrum view



Note You can set a minimum trace height value in centimeters on the Default Workspace Options page. When you adjust the height of the Spectrum view, if its height becomes smaller than the set minimum value, a scrollbar automatically appears to the right of the view.

The scan header describes the experiment. [Table 30](#) describes the default scan header information. If the scan header is not visible, resize the Spectrum view.

Table 30. Default header information symbols

Symbol	Description
#S	Scan number
RT	Retention time
AV	Averaged (followed by the number of averaged scans)
SB	Subtracted (followed by the number of subtracted scans and the scan range)
NL	Normalization level
T	Scan filter string

Right-clicking a Spectrum or Multi Spectrum view displays a shortcut menu ([Figure 102](#)) with the commands described in [Table 31](#).

Figure 102. Shortcut menu for the Spectrum view (left) or the MultiSpectrum view (right)

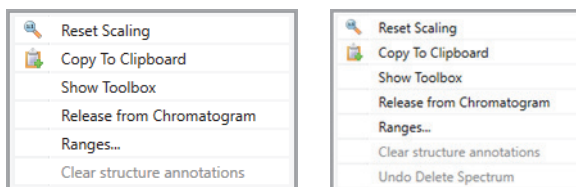


Table 31. Spectrum (or MultiSpectrum) view shortcut menu commands

Command	Description
Reset Scaling	Resets the scaling of the plot in the Spectrum view or all the plots in the MultiSpectrum view.
Copy To Clipboard	Copies an image of the Spectrum view to the Clipboard.
Show Toolbox	Displays the Spectrum toolbox. To close the toolbox, click its pin icon.
Release from Chromatogram	Releases the selected spectrum from its chromatogram.
Ranges	Opens the Spectrum Ranges dialog box for defining the spectrum.
Clear Structure Annotations	Clears the structure annotations from the view and the structure list from the Structures Annotation dialog box.
Undo Delete Spectrum (MultiSpectrum view only)	Undoes the action of deleting a spectrum in a MultiSpectrum view.

These toolbars are available when a Spectrum view or Multi Spectrum view is selected:

- [Spectrum – Workspace Processing Toolbar](#)
- [Spectrum – Display Options Toolbar](#)
- [Zoom Options Toolbar](#)
- [Text and Graphic Annotation Toolbar](#)

MultiSpectrum View

Use a MultiSpectrum view to display multiple spectrum traces when you want to compare the traces using the same normalization level or you want to simultaneously zoom in on the same mass range.

Note In the Normalization area of the Spectrum – Display Options toolbar, the Global button is unavailable when the spectra being compared are in two separate views.

Instead of zooming in on each spectrum individually in a Spectrum view, use a Multi Spectrum view to zoom in and out on all of the spectra at once. In a Multi Spectrum view, the selected spectrum (the spectrum with the darker gray background) is the only one that updates when you click the chromatogram.

You can perform background subtraction, Xtract deconvolving and deisotoping, isotope simulation, or library searches on any selected spectrum in this view. You can also use all the functions in the Spectrum toolbox.

The MultiSpectrum view's shortcut menu contains the same commands as that for the Spectrum view with the exception of one additional command—Undo Delete Spectrum.

To create and work with a MultiSpectrum view, see [Creating a MultiSpectrum View and Changing Its Spectrum Plots](#).

For more information about the MultiSpectrum view, its shortcut menu, and its toolbox, see [Spectrum View](#) and [Spectrum Toolbox](#).

Spectrum Ranges Dialog Box

Use the Spectrum Ranges dialog box to select the spectra to display, create an average spectrum, and perform background subtraction.

[Table 32](#) describes the columns in the Spectrum Ranges dialog box.

Table 32. Spectrum Ranges dialog box

Column	Description
Display	Specifies whether to add the defined spectrum to the Spectrum or MultiSpectrum view.
Release	Specifies the linkage status of the defined spectrum. Selected: Releases the spectrum from the chromatogram. Clear: Links the spectrum to the chromatogram.
File Name	Specifies the location and file name of the raw data file with the spectral data of interest. If the spectrum is not linked to a chromatogram, a browse button is available for browsing to and selecting a different raw data file.
Detector Type	Specifies the detector used to acquire the data. Unavailable when the spectrum is linked to a chromatogram.
Filter	Specifies the scan filter (same selections as the Scan Filters page of the Info bar). Unavailable when the spectrum is linked to a chromatogram.
Retention Time	Specifies the retention time for a single spectrum or the retention time range for an average spectrum. The Retention Time and Scan Number columns are interactive; that is, changing the retention time range updates the scan number range.
Scan Number	Specifies the scan number for a single spectrum or the scan number range for an average spectrum.
Subtract Background	Specifies whether to perform background subtraction or whether to subtract the scans from one or two time ranges.
Background Range 1	Specifies the time range for one set of background scans. Format: Time _{Start} -Time _{End} , where the start time point is earlier than the end time point
Background Range 2	Specifies the time range for a second set of background scans.

Spectrum Toolbox

In a Spectrum view or a Multi Spectrum view, use the Spectrum toolbox (Figure 103) to run an elemental composition analysis, add an EIC plot to the Chromatogram view, set up relative mass labels, or zoom in or out of the view. In a Deconvoluted Spectrum view, use the toolbox to set up relative mass labels or zoom in or out of the view.

How you open the floating toolbox depends on your workspace setting (see [Default Workspace Options Page](#)). With the factory default setting, you can open the toolbox in the following way:


Right-click an m/z value in the view and choose **Show Toolbox**. The toolbox opens with its pin icon in the pinned (vertical) position, .

Figure 103. Spectrum toolbox

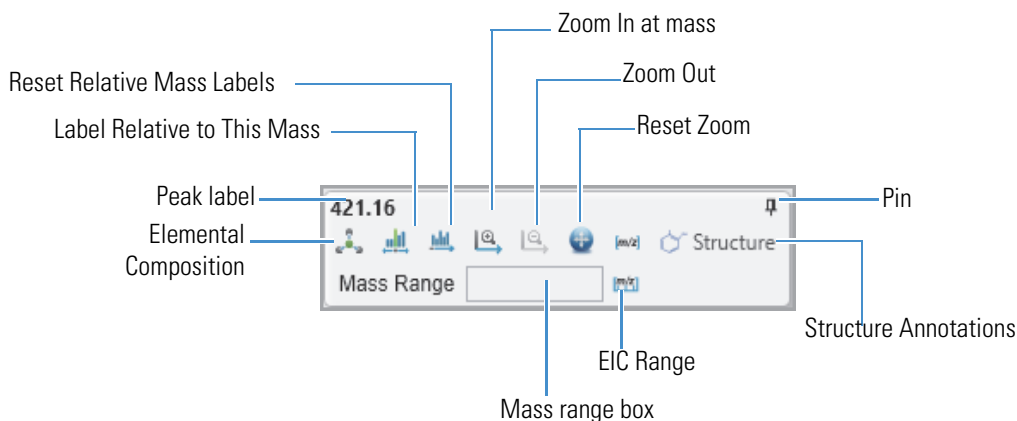


Table 33 describes the icons and the Mass Range box in the Spectrum toolbox.

Table 33. Spectrum toolbox icons (Sheet 1 of 2)

Icon/field	Description
Elemental Composition (Not available for the Deconvoluted Spectrum view)	Finds the best matching calculated chemical formulas for the selected peak in the spectrum. See Starting an Elemental Composition Analysis from the Toolbox .
Label Relative to This Mass	Labels each peak in the spectrum with the difference between its mass-to-charge ratio and the mass-to-charge ratio of the selected peak (in m/z).
Reset Relative Mass Labels	Labels each peak in the spectrum with its mass-to-charge ratio (in m/z).
Zoom In	Moves in on the spectrum.
Zoom Out	Moves out from the spectrum.
Reset Zoom	Resets the zoom to display the entire mass range.

Table 33. Spectrum toolbox icons (Sheet 2 of 2)

Icon/field	Description
EIC Mass (Not available for the Deconvoluted Spectrum view)	Uses the entered individual mass-to-charge ratio values to plot an extracted ion chromatogram (EIC).
EIC Range (Not available for the Deconvoluted Spectrum view)	Uses the entered mass-to-charge ratio range to plot an extracted ion current (EIC) chromatogram.
<p>Note For information about using the Spectrum toolbox to display an EIC chromatogram, see Displaying an EIC Trace by Using the Spectrum Toolbox.</p>	
Mass Range box (Not available for the Deconvoluted Spectrum view)	Enter a mass range within the spectrum m/z limits and then press ENTER. The view rescales to display the spectrum within the entered range.

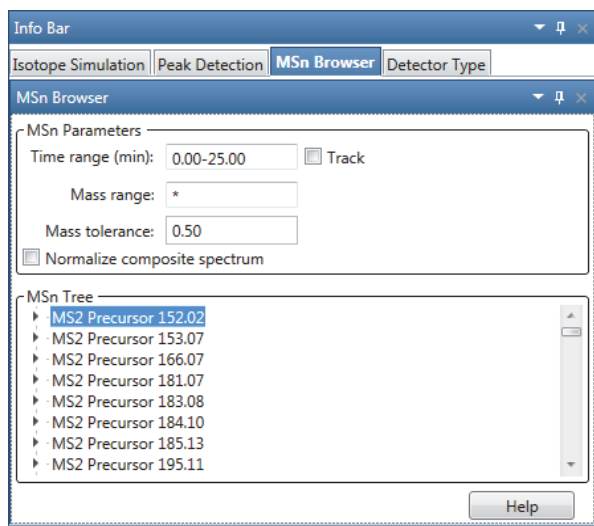
MSn Browser Page

Use the MSn Browser page in the Info Bar to select spectra of interest from an MS tree view of the spectra in a raw data file.

For information about displaying spectra, see [Selecting Spectra from an MSn Tree](#).

[Figure 104](#) shows the MSn Browser page, and [Table 34](#) describes the parameters on the this page.

Figure 104. Info Bar – MSⁿ Browser page (with a collapsed MSn tree)



Note The MSⁿ browser information page is not available for all Thermo Scientific mass spectrometers.

Table 34. MSn Browser page parameters (Sheet 1 of 3)

Parameter	Description
MSn Parameters	
Time Range (min)	<p>Specifies the time range of the MSn Tree (see To filter the MSn tree by a selected time range).</p> <p>Limits the spectra available for display in the MSn tree to MS2 or higher level scans acquired during the specified time range.</p> <p>There are two ways to change the Time Range:</p> <ul style="list-style-type: none"> Type the time range in minutes in the Time Range box. The format is <i>From-To</i>. Select the Track check box, and then drag the pointer horizontally across the Chromatogram view from the minimum time to the maximum time of interest. <p>Default: 0.00-<i>acquisition time</i> (in minutes)</p>
Track	<p>Activates the selection of a time range by using the mouse pointer. To reset the time range, right-click the Chromatogram view and choose Reset Scaling.</p> <p>Default: Clear</p>
Mass Range	<p>Specifies the mass range of the MSn Tree (see To filter the MSn tree by a selected mass range).</p> <p>The format is <i>From-To</i>.</p> <p>Default: * (Mass range from the instrument method)</p>
Mass Tolerance	<p>Specifies the mass tolerance for grouping the precursor nodes and their associated scans. Decreasing the mass tolerance to 0.00 creates an MSn Tree where most, if not all, the individual scans are associated with a separate MS2 precursor node (see To modify the grouping of the scan data in the MSn tree).</p> <p>Default: 0.05 <i>m/z</i>; Range: 0.00–10.00 <i>m/z</i></p>

Table 34. MSn Browser page parameters (Sheet 2 of 3)

Parameter	Description
Normalize Composite Spectrum	<p>Normalizes the spectral peaks in a composite spectrum so that you can view low-intensity peaks from the higher-order fragmentation spectra (see To normalize a composite spectrum).</p> <p>Each MSⁿ spectrum is individually normalized (NL) so that its highest peak is displayed at a Relative Abundance of 100%; therefore, the relative peak heights of this display are not meaningful. For example, a composite spectrum for an MS³ experiment displays both the MS² base peak and the MS³ base peak at a Relative Abundance of 100% (unless the base peaks in the MS2 and MS3 are within the specified mass tolerance) and maintains all other relative abundances of the other ions in each spectrum. If the base peaks in the MS2 and MS3 scans are within the specified mass tolerance, the application averages these peaks, causing the Relative Abundance of the MS3 base peak to be less than 100%.</p>
MSn Tree	
MS2 Precursor nodes	Display the precursor masses that triggered the MS2 level scans. To display the spectrum tree for each node, click the expand icon to the left of the node, or right-click the MSn Tree view and choose Expand List .
MSn Precursor nodes	Display the precursor masses that triggered the MSn level scans.
Average Spectrum	Double-click to display a Spectrum view with the averaged spectrum for the selected MSn level. The MSn tree includes an average spectrum when the data file includes scans for the selected scan filter.
Composite Spectrum	<p>Double-click to display a Spectrum view with the composite spectrum for all the MSn levels down to the selected MSn level.</p> <p>A composite spectrum for an MSn level is a summed spectrum of all the fragmentation scans derived from the same original precursor mass down to the current level.</p>
Single Spectrum or Scan Number at Time	Double-click to display a Spectrum view for the specified scan number.
Shortcut menu	
Include Individual Scans	Displays the individual scans (see To display all the individual scan items).
Normalize Composite Spectrum	Normalizes the composite spectra (see To normalize a composite spectrum).

Table 34. MSn Browser page parameters (Sheet 3 of 3)

Parameter	Description
Expand List	Expands all the MSn Precursor nodes (see To expand the entire MSn tree). Available when the MSn tree is collapsed.
Collapse List	Collapses all the MS2 Precursor nodes. Available when the MSn tree is expanded.
Export	Copies the MSn Tree to the Clipboard.
Print	Opens the Print dialog box where you can set up your print preferences and print the contents of the MSn Tree.

Reviewing Map Data

Use the Map View-specific toolbars and display options of the FreeStyle application to review the map view data from mass spectrometer.

These procedures describe how to set the ranges and display options for a map view.

Contents

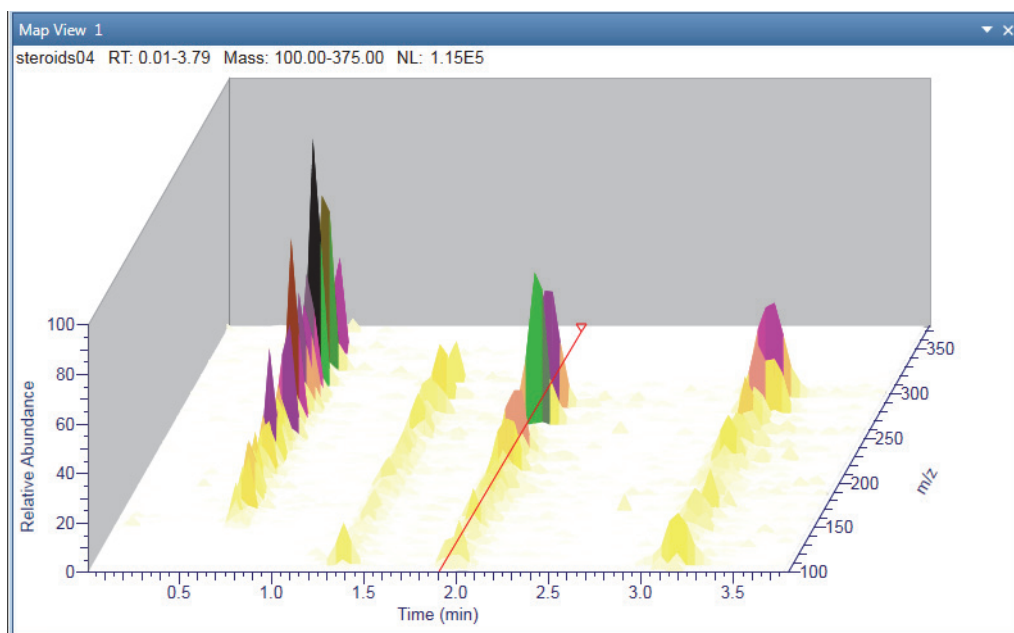
- [Working with the Map View](#)
- [Adding Map Views to the Workspace](#)
- [Changing the Zoom level of a Map View](#)
- [Linking and Releasing Map to and from a Chromatogram](#)
- [Map View - Specific Toolbars](#)
- [Map Ranges Dialog Box](#)
- [Setting the Map Display Options](#)
- [Setting the Map Format Options](#)

Working with the Map View

A map is a 2D or 3D representation of an analysis showing the mass scans acquired during an analysis. Use the Freestyle application to open a raw data file, create map views, and customize using the display options.

The Map view consists of a time point (x axis) versus an m/z value (z axis) versus a relative abundance value (y axis) map plot. The scan header contains File name, RT range, Mass range, NL values, and filter details. These values change based on the changes in the Map view.

Figure 105. Map view example



Adding Map Views to the Workspace

You can add multiple Map views to the workspace.

❖ To add a Map view

1. Open a raw data file (see [Opening Raw Data Files or Sequence Files](#)).
2. With the factory default layout, the Chromatogram view appears at the top and the Spectrum view appears at the bottom.
3. In the Workspace Options toolbar, click **Map View**.

A new Map view opens for each selected chromatogram with RT on the x axis and m/z on the z axis.

Note If the Workspace includes one or more Chromatogram views, a map view of the selected or active Chromatogram view appears.


If the Workspace does not include a Chromatogram view, a map view of the chromatogram associated with the active spectrum view appears.

You can create multiple map views, but a single map view cannot have multiple traces. To increase the intensity, use the **Up** and **Down** arrow keys on your computer.

Displaying the Scan for a Time Point in Map View

❖ To display the scan corresponding to a specific time point

1. Select the Map view.
2. Click any time point.

In the Map view, a red vertical marker, , indicates the selected data point. The Spectrum view displays the spectrum for that retention time or scan number. The Retention Time marker position in the chromatogram also is updated correspondingly to sync with the Map view.

Linking and Releasing Map to and from a Chromatogram

When you create a Map view in Workspace, the application automatically links the map view to the active chromatogram trace. The displayed map view changes when you click another trace in the linked chromatogram view. If you do not want to change the map view, you can release the map view from the trace and link to another chromatogram trace.

See these topics:

- [Linking Status of a Map View](#)
- [Releasing a Linked Map View from the Chromatogram](#)
- [Linking a Map View to a Chromatogram](#)

Linking Status of a Map View

When you link or release the map view, the application appends the linking status to the view's name.

Figure 106. Map View Linking Status

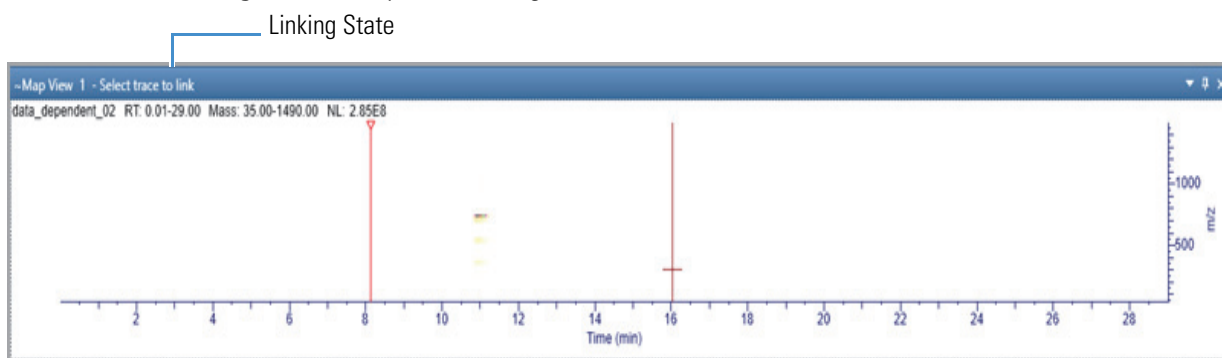


Table 35 describes the different linking status for a Map view.

Table 35. Map View Linking Status

Linking status	Description
Link to chromatogram	<p>When a Map view is linked to a chromatogram, the title bar of the Map view displays the linked chromatogram by its view number and trace number in the following format:</p> $CxTy$ <p>where</p> <p>x = the number of the Chromatogram view</p> <p>y = the number of the trace in the Chromatogram view</p>
Available for linking	<p>The following text is appended to the title bar of the Map view—Select Trace to Link.</p> <p>When a map view is available for linking, clicking any chromatogram in the Workspace links the view to the selected chromatogram, updates the RT marker in Map View and displays the spectrum based on the selected time point in the Chromatogram trace, and changes the appended text in the view's title bar to $CxTy$.</p>
Released	<p>When you release a Map view from its chromatogram, the text appended to the Map view changes to Released. In the Released state, the Map view is not affected by any actions in the Chromatogram view. The Spectrum view also does not change as you navigate to different retention times. To update a released Map view, you must use the Map Ranges dialog box.</p>

Releasing a Linked Map View from the Chromatogram

In Map view, the title bar of the view's name indicates the linking status of the Map view,

❖ To release a Map view from its linked chromatogram

Right-click the Map view and choose **Release from Chromatogram**.

The application releases the link to the chromatogram trace and appends the following text to the view's name: **Released**.

Note When you close the chromatogram trace linked to a Map view, the application automatically releases the linking status. The Map view displays the RT marker, which is not synced with a chromatogram or spectra.

Linking a Map View to a Chromatogram

❖ To link a map to a chromatogram

1. Right-click the map and choose **Link to Chromatogram**.

The application appends the following text to the view's name: **Select trace to link**.

2. Click the chromatogram trace to link.

The application links to the chromatogram trace and appends the following text to view's name:

$CxTy$

where x indicates the Chromatogram view's number and y indicates the trace number.

Changing the Zoom level of a Map View

You can reset, zoom in, or out of a Map View using the mouse pointer, Zoom options toolbar, or right-click menu options.

To understand more about zooming a map view, follow these procedures:

- [Using the Mouse Pointer](#)
- [Using the Zoom Options Toolbar](#)

Using the Mouse Pointer

❖ To zoom in and out using the mouse pointer

1. Select the view.
2. To zoom in the retention time (x scale), position the mouse pointer and drag horizontally across the specific section. Release the mouse to zoom.
3. To zoom in the m/z scale, position the mouse pointer and drag vertically across the specific section. Release the mouse to zoom.
4. To zoom in both x and m/z scales, click and drag the mouse within the map to form a box containing the region to zoom. Release the mouse to zoom.
5. To return to the previous scale, right-click and select **Undo Zoom**.

Using the Zoom Options Toolbar

❖ To zoom in and out on the x axis

1. Select the view.
2. To zoom in, from the Zoom Options toolbar, click **Zoom In X**.

Each click zooms in the x axis larger by a factor of two (2) to show more detail. For example, you can change the x -axis range from 0–20 to 5–15.

3. To zoom out, click **Zoom Out X**.

Each click increases the Time (min) range from the center. For example, you can change the x -axis range from 7.5–12.5 to 5–15.

❖ **To zoom in and out on the y axis**

1. Select the view.
2. To zoom in, from the Zoom Options toolbar, click **Zoom In Y**.

Each click zooms in on the y axis. by a factor of two from the current baseline.

3. To zoom out, click **Zoom Out Y**.

Each click zooms out on the y axis by a factor of two. For example, you can change the y -axis range from 0–25 to 0–50.

❖ **To reset, undo and redo zoom of the x and y axes**

1. To reset scaling, do the following:
 - Select the view and from the Zoom Options toolbar, click **Reset**.–or–
 - Right-click the view and choose **Reset Scaling**.
2. To undo or redo zooming, right-click the view and choose **Undo Zoom** or **Redo Zoom** respectively.

Map View - Specific Toolbars

Use the following map view-specific toolbars to create, format, and set map ranges and display options of map views.

- [Map View - Workspace Options Toolbar](#)
- [Map View - Display Options Toolbar](#)

Map View - Workspace Options Toolbar

Use the buttons in the Workspace area of the Workspace Options toolbar to create map views and set the map ranges.

❖ **To display the Map view specific Workspace Options toolbar**

1. Click the chromatogram or spectrum to select it.

- In the Workspace Options toolbar, click **Map View** to create the map view.

The Map view specific workspace options appear.

Figure 107. Map View – Workspace Options toolbar

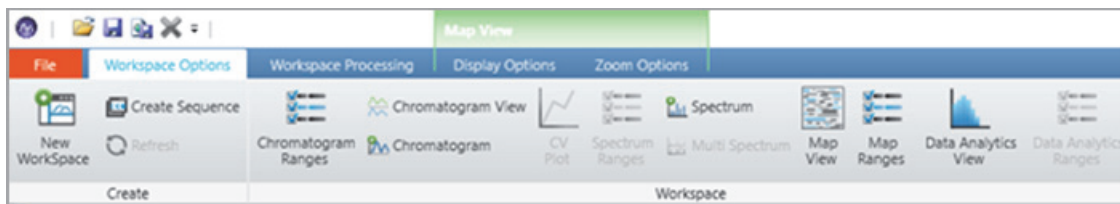


Table 36. Workspace Options–Workspace area buttons for the Map view

Button	Description
Workspace	
Use the buttons in the Workspace area to create a Map view and set the ranges.	
Map View	Creates a map view of the selected chromatogram or spectrum.
Map Ranges	Sets the filename and scan filter for the map.

Map View - Display Options Toolbar

Use the buttons in the Display Options toolbar to customize the Map view. You can format the 3D or 2D representation of the map view, define the colors based on intensity, label the axes, and specify the band width.

❖ To display the Map View – Display Options toolbar

- Click a Map view to select it.
- Click the **Display Options** toolbar tab.

The Map View specific display options appear.

Figure 108. Map View – Display Options toolbar

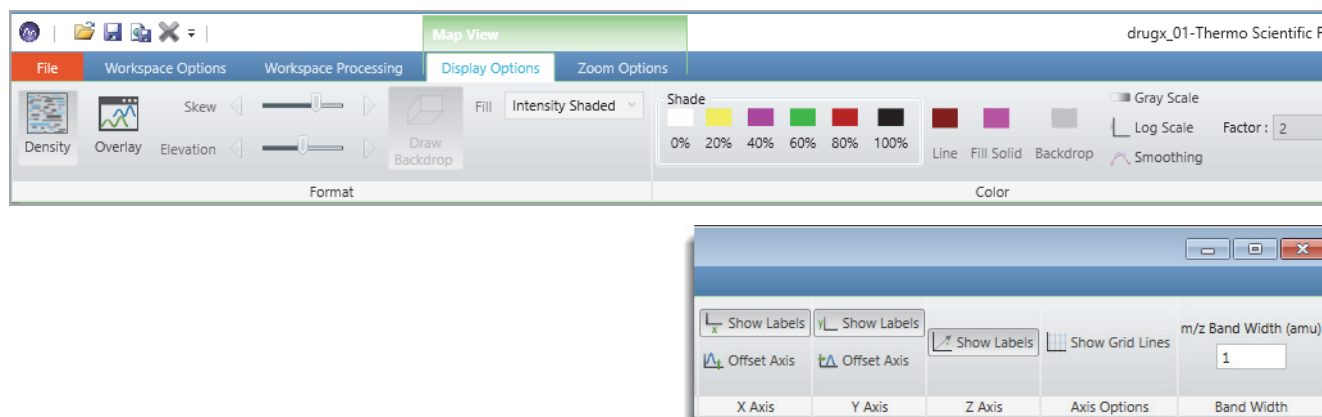


Table 37. Map View – Display Options toolbar buttons (Sheet 1 of 2)

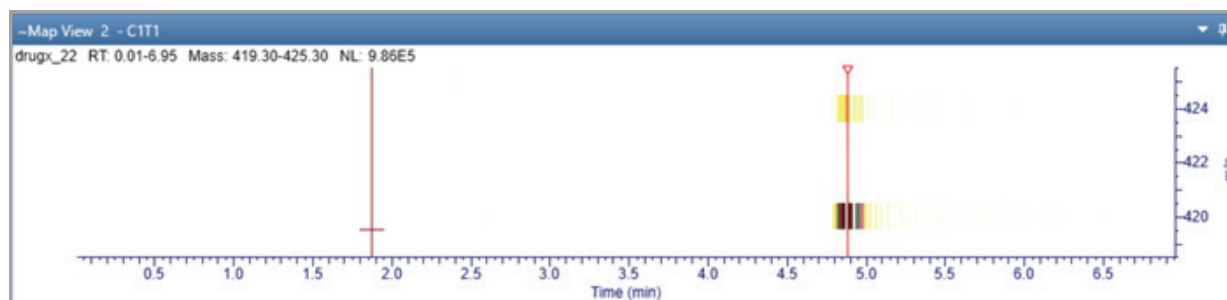
Button	Description
Format	
Use the parameters in the Format area to represent the Map view in 3D or 2D.	
Density	Displays different shades in the 2D representation of the map view, where colors represent the <i>m/z</i> bands categorized based on the intensity percentage.
Overlay	Displays the 3D representation of the map view.
Skew	Sets the skew angle (time offset) to a value from 0–45 degrees for an overlay arrangement of Map View.
Elevation	Sets the vertical spacing for an overlay arrangement of Map View.
Draw Backdrop	Adds a backdrop to overlaid views.
Fill	Sets the filling style for map view. This is applicable only for 3D representation of the map view (when Overlay is selected).
Color	
Displays the color palette, where you select the colors of the traces in the Map View.	
Line	Sets the color of the framing lines
Fill Solid	Sets the color of the solid fill.
Backdrop	Changes the color of the backdrop
Gray Scale	Plots the map in gray scale
Log Scale	Displays and clears the color of the map in a logarithmic scale.
Smoothing	Displays a smoothed density plot for the color bands (bins of <i>m/z</i> values) in the vertical direction. This is applicable only for 2D representation of the map view.
X Axis	
Use the parameters in the X Axis area to specify how the application labels and displays the <i>x</i> axis.	
Show Labels	Shows or hides the <i>x</i> -axis label.
Offset Axis	Sets the location of the displayed plot at a specified distance from the <i>x</i> axis.
Y Axis	
Use the parameters in the Y Axis area to specify how the application labels the <i>y</i> axis.	
Show Labels	Shows or hides the <i>y</i> -axis label.

Table 37. Map View – Display Options toolbar buttons (Sheet 2 of 2)

Button	Description
Offset Axis	Sets the location of the displayed plot a specified distance from the <i>y</i> axis.
Z Axis	
Use the parameter in the Z Axis area to specify how the application labels the <i>z</i> axis.	
Show Labels	Shows or hides the relative abundance label.
Axis Options	
Use the parameter in the Axis Option area to create grid lines in the map view.	
Show Grid Lines	Displays or clears the grid lines.
Band Width	
Use the parameter to specify the band width in amu units.	
<i>m/z</i> Band Width (amu)	Specifies a band width value from 0.001 to 50.

Map-Specific Views

The Map view displays a 2D or 3D representation of the map view of the selected chromatogram or spectrum. The title bar of the view displays the file name of the raw data file, RT range, Mass range, NL values, and the view's link status. C1T1 specifies that the view is linked to Trace 1 in the Chromatogram1 view.

Figure 109. Map View Title bar

Map Ranges Dialog Box

Use the Map Ranges dialog box to select the raw data file to display in the Map view, detector type used to acquire the data, scan filter, and time and mass ranges.

❖ To open the Map Ranges Dialog box

In the Workspace Options toolbar, click **Map Ranges**.

Figure 110. Map Ranges Box

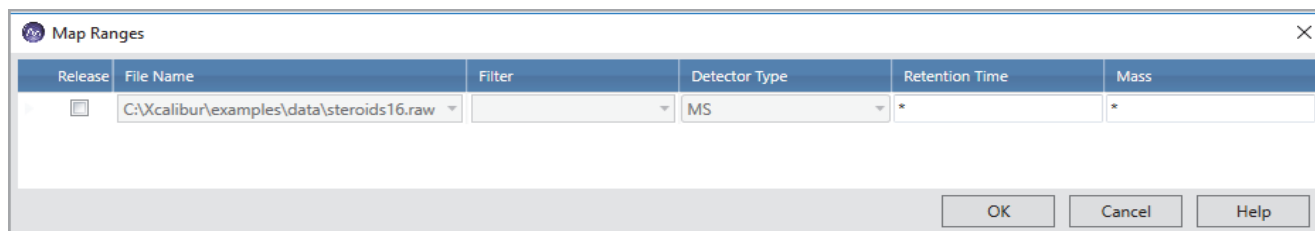


Table 38 describes the columns in the Map Ranges dialog box

Table 38. Map Ranges dialog box

Column	Description
Release	<p>Specifies the linking status of the Map view.</p> <p>Selected: Releases the map view from the chromatogram. Lets you to change the Filter name and raw file.</p> <p>Clear: Links the map view to the chromatogram. You cannot change the raw file and filter name.</p>
File Name	<p>Specifies the location and file name of the raw data file.</p> <p>If the map view is not linked to a chromatogram, a browse button is available to select a different raw data file.</p>
Detector Type	<p>Specifies the detector used to acquire the data. By default, the detector type is set to MS and is non-editable.</p> <p>Unavailable when the map view is linked to a chromatogram.</p>
Time	<p>Specifies a valid time range based on the raw file settings. Default: '*'. <ul style="list-style-type: none"> where '*' indicates All. </p>
Mass	<p>Specifies a valid mass range based on the raw file settings. Default: '*'. <ul style="list-style-type: none"> where '*' indicates All. </p>

Setting Map Ranges

❖ To set the file and filter options

1. Click a map view to make it active.
2. From the Workspace Options toolbar, click **Map Ranges**. By default, the map view is linked to the chromatogram.

3. To change the raw file and filter options, in the Map Ranges dialog box, select the release checkbox.
4. From the File Name list, click the dropdown arrow to select a file name or click the browse button to select a different raw data file.
5. In the Filter box, type or select the scan filter from the dropdown list, which displays filter options stored in the raw data file.
6. In the Retention Time and Mass boxes, enter a valid retention time and mass range values.
7. To save the settings and close the Map Ranges dialog box, click **OK**.

Defining the Retention Time and Mass Ranges

You can define the time and mass ranges in the Retention Time and Mass boxes as follows:

- Enter valid start and end numeric values for the RT and mass ranges.

-or-

- Set the start value as '*' and enter a valid end value. The application uses the default start value from the raw file as the start value.

-or-

- Defines a valid start value and enter the end value as '*'. The application takes the default end value from the raw file as the end value.

Saving a Map View

You can save customized map views in the layout in Freestyle.

❖ To save a customized Map View in the layout

1. Select the Map view you want to save.
2. Click the **Workspace Options** toolbar tab.
3. From the Layouts area, choose the option to save the map view.

The Map view is saved in the layout.

When you apply the saved map view in another view, it displays details, such as title bar number, state(s) (Dockable, Floating, Auto Hide), and the position of the map as it has been saved in the layout. It also displays the map ranges parameters and the status of the Map view (Linked or Released).

Setting the Map Display Options

Use the Display Options toolbar to set up the display options for a map view.

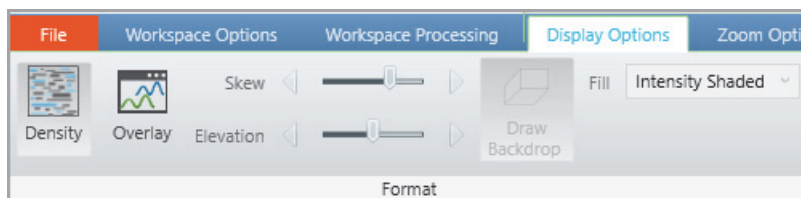
❖ To open the Display Options dialog box for a map view

1. Click the map view to make it active.
2. From the ribbon menu, click **Display Options**.
3. Follow one or more of these procedures to set up the map display options:
 - [Setting the Map Format Options](#)
 - [Setting the Map Axis Options](#)
 - [Setting the Map Color Options](#)
 - [Setting the Band Width](#)

Setting the Map Format Options

Use the Display Options toolbar to format the Map View.

Figure 111. Map Format options



❖ To set the map format options

1. Click a map view to make it active.
2. From the menu bar, click **Display Options**.
3. To specify the arrangement style, select one of these options:
 - To display a density map showing different shades for each intensity, select the **Density map** option. The default display option is Density.
 - To overlay plots vertically with optional horizontal skew (time offset) for the active map, select the **Overlay (3D)** option. The 3D view of the map appears.
4. To specify style options for overlaid (3D) plots, do the following in the Format area:
 - To set the elevation angle (from 0 to 60 degrees), drag the Elevation slider or click the left or right arrow on the Elevation slider until you reach the desired angle.
 - To set the skew angle (from 0 to 45 degrees), drag the **Skew** slider or click the left or right arrow on the Skew slider until you reach the desired angle.

The default values for the Skew and Elevation sliders are set at 30.

- To select a different filling style for the map view, select a fill option from the Fill list.

The default fill pattern is intensity shaded.

The styles control the display of filling and outlining data within the map. Outlining (wireframing) produces lines between the scans (vertical or diagonal based on the skew) and the mass bands (horizontal). Data with each scan and mass band is displayed as either unfilled, filled with a solid color, or filled with a shade based on the intensity of the data (using the shade colors as represented in the 2D density map). Some modes draw a vertical line, which represents the intensity of each mass band at each scan.

Table 39 explains the available styles, and how they control filling and outlining (wireframe).

Tip In the table, nonzero data refers to data, which when scaled to a color, would not provide a background color. If you want to view the smaller values of the plot, press the **UP** arrow key on the computer to increase the intensity scaling or set the log scale factor.

- To add a backdrop to 3D plots, click **Draw Backdrop**. To remove a backdrop, click the **Draw Backdrop** button again.

Table 39. Style options for overlaid (3D) plots (Sheet 1 of 2)

Modes	Fill	Frame	Vertical Line
Plain Lines	None	None	A vertical line at each scan and mass/band having non-zero data. All the lines are displayed of same color. To differentiate the lines, the bottom of each line is highlighted in tinted green.
Colored Lines	None	None	A vertical line at each scan and mass/band having non-zero data. The line is colored based on the intensity of the shade selected from the color palette
None	None	Creates a wireframe plot, showing small amounts of data that might otherwise hidden by large data sets	None
Solid Color	All non-zero bands are filled using the same color	Creates a wireframe that helps to separate the bands	None

Table 39. Style options for overlaid (3D) plots (Sheet 2 of 2)

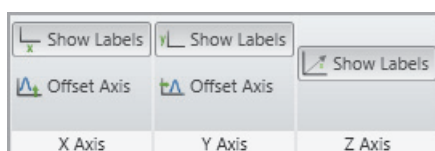
Modes	Fill	Frame	Vertical Line
Intensity shaded	All non-zero bands are filled and colored based on the shade selected from the color palette.	None	None
Shaded with Frame	All non-zero bands are filled and colored based on the shade selected from the color palette.	Creates a wireframe that helps to separate the bands	None

Setting the Map Axis Options

The Map axis display options are based on the selected map view. For Density view, the default x -axis is Time (min) and z -axis is m/z . For the overlay 3D view, there are three axes as follows:

- x axis - Time (min)
- y axis - Relative abundance
- z axis - m/z

Figure 112. Display Options – Map Axis



❖ To set the map axis options

1. Click a map view to make it active.
2. From the menu bar, click **Display Options**.
3. To show or hide the labels of x , y , or z axis, click the corresponding **Show Labels** buttons.
4. To display the grid lines, click **Show Grid Lines**.

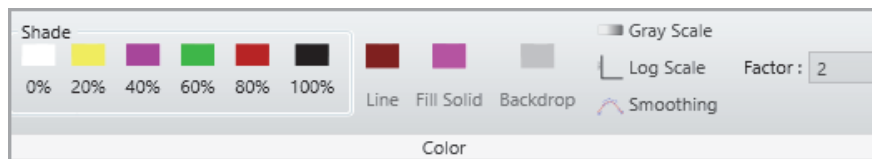
Setting the Map Color Options

Use the color options in the Display Options toolbar to set the color of the traces.

❖ To set the map color options

1. Click a map view to make it active.
2. From the ribbon menu, click **Display Options**.

Figure 113. Map Color options



3. To select the color of the framing lines, from the Fill dropdown list, select **Plain Lines**.
4. Click the Line color box above **Line**.
The Color dialog box opens with a color palette that enables you to select a preset color or customize a color.
5. To select the color of the solid fill for the active map, from the Fill dropdown list, select **Solid Color**.
6. Click the color box above **Fill Solid**.
The Color dialog box opens with a color palette that enables you to select a preset color or customize a color.
7. To select the color of the backdrop (background), click the Backdrop Color box above **Backdrop**.
The Color dialog box opens with a color palette that enables you to select a preset color or customize a color. Backdrop is active only when you select the Overlay 3D style.
8. To select Gray Scale or color, do one of the following:
 - To plot the map in gray scale, click the **Gray scale** button.
 - To plot the map in color, click the **Gray scale** button again.
9. Use the shade buttons (0%, 20%, 40%, 60%, 80%, and 100%) to change the map's color at 0%, 20%, 40%, 60%, 80%, and 100% relative abundance.
10. To select log scale or linear scale, do the following:
 - To display the color of the map in a logarithmic scale, click the **Log Scale** button.
The factor width that you set in the Factor box determines the scaling between color bands.
 - To display the map in a linear scale, click the **Log Scale** button again.
11. To select smoothing, click **Density** view and then click **Smoothing**.
The solid color bands of the map view are converted to gradient color bands and the most intense band appears at the middle and gradually fades out to the upward and downward directions.

Setting the Band Width

Use the Band Width option to specify the bandwidth in amu units. The m/z axis of the map view plot is directly related to the band width value entered in the m/z Band Width box. The default band width is set at 1.0 amu.

The Map View does not display all the peaks in scans. It merges data into mass bands displaying the most intense value in each scan. For example, when you scan a file from m/z 110 to 120, the map creates bands centered on each nominal mass, such as 109.5–110.5, 100.2–101.5, and so on. The view displays the most intense peak in each band. To increase the mass resolution, reduce the band size.

❖ To define the band width in a map view

1. Open a raw data file and make a map view the active view.
2. From the menu bar, click **Display Options**.
3. On the Band Width area, in the m/z Band Width (amu) box, type a value from **0.001** to **50**.

Figure 114. Display Options – Band Width



4. To save the setting and close the dialog box, press the **Enter** key on your computer keyboard.

Analyzing MS Data using Data Analytics View

Use the Data Analytics View display options in FreeStyle application to analyze the MS trending information from mass spectrometers. You can use the view to plot the chosen trace type properties of an MS trending detector type. You can also analyze different aspects such as frequency of occurrence, retention times, the quantity being plotted, categories, and so on. The analysis is available as a Histogram or Trend report.

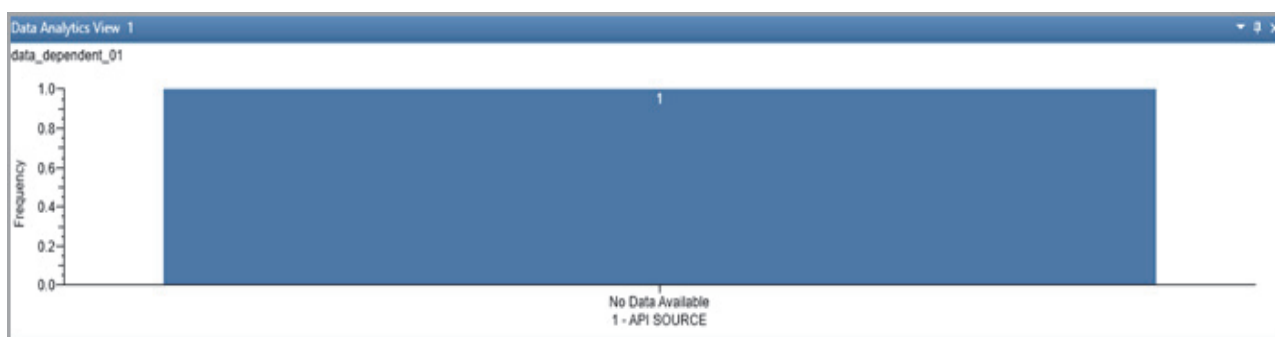
Contents

- [Adding Data Analytics View to the Workspace](#)
- [Plotting the Data as a Histogram](#)
- [Data Analytics View - Specific Toolbars](#)
- [Saving a Data Analytics Plot](#)
- [Data Analytics Ranges](#)

Adding Data Analytics View to the Workspace

You can add multiple data analytics views to the workspace.

Figure 115. Data Analytics View



❖ To add a Data Analytics View

1. Open a raw data file (see [Opening Raw Data Files or Sequence Files](#)).

6 Analyzing MS Data using Data Analytics View

Plotting the Data as a Histogram

2. With the factory default layout, the Chromatogram view appears at the top and the Spectrum view appears at the bottom.

3. In the Workspace Options toolbar, click **Data Analytics View**.

A new Data Analytics view opens for the active chromatogram. A Histogram plot displays the instrument status reading on the x axis and frequency on the y axis; and a Trend plot displays the time on the x axis and readback parameter on the y axis.

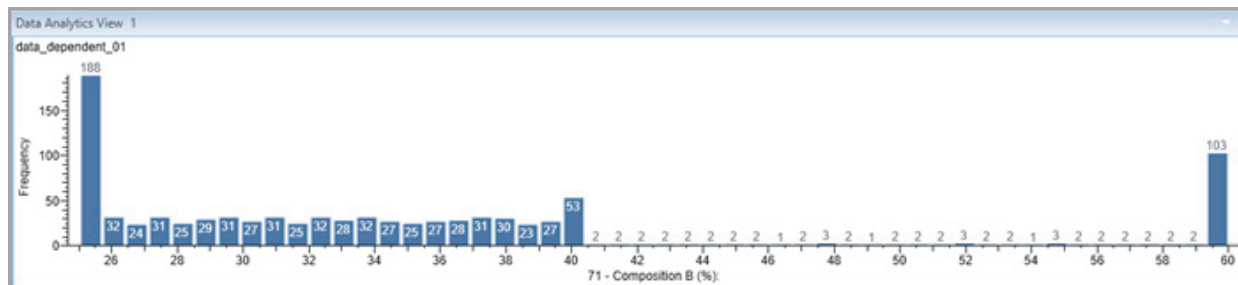
If you have selected the MS Trending detector type, the Data Analytics View process the information based on the selected trace type.; otherwise, the first trace type is used to process the information.

Note You can add multiple Data Analytics views in the workspace. If the Workspace includes one or more Chromatogram views, a Data Analytics view of the selected or active trace of the Chromatogram view appears.

Plotting the Data as a Histogram

When you create a new Data Analytics view with MS trending selected as the preferred trace type, a Histogram plot displays the instrument status reading plotted against the frequency. The data groups are represented as bins in the plot.

Figure 116. Histogram



The MS Trending detector type data is displayed as follows in the view:

- The vertical axis of the plot is Frequency (or Number of Occurrences). The title of the vertical axis is Frequency.
- The horizontal axis of the plot is a binned representation of the data groups. The title of the horizontal axis is the name of the readback parameter.
- If representing categorical data, the categories are plotted as bars on the horizontal axis and number of occurrences are plotted on the vertical axis. In this representation, bins are not created.

Note If the raw data file is a non-MS trending file, the application does not display a histogram plot.

Plotting the Data in a Trend Chart

When you plot the data in a Trend chart, the plot displays the trace type readback parameter plotted against the time.

The MS Trending detector type data is displayed as follows in the view:

- The vertical axis of the plot is the readback parameter selected as the trace type. The title of the vertical axis is the name of the selected readback parameter.
- The horizontal axis of the plot displays the readback trace during specific time period.

Data Analytics View - Specific Toolbars

Use the following Data Analytics view-specific toolbars to create, format, and set data analytics ranges and display options of the view.

- [Data Analytics View - Workspace Options Toolbar](#)
- [Data Analytics View - Display Options Toolbar](#)

Data Analytics View - Workspace Options Toolbar

Use the buttons in the Workspace area of the Workspace Options toolbar to create data analytics views and set the data analytics ranges.

❖ To display the Data Analytics view-specific Workspace Options toolbar

1. Click the chromatogram to select it.
2. In the Workspace Options toolbar, click **Data Analytics View** to create the view. The Data Analytics view-specific workspace options appear.

Figure 117. Data Analytics View-specific Workspace Options

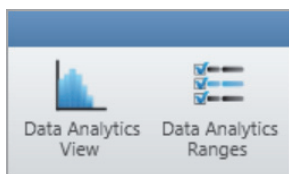


Table 40. Data Analytics View - Workspace Options toolbar buttons

Button	Description
Workspace	
Use the buttons in the Workspace area to create a Data Analytics view and set the ranges.	
Data Analytics View	Creates a data analytics view of the selected chromatogram.
Data Analytics Ranges	Sets the filename, detector type, trace type, and filter for the view.

Data Analytics View - Display Options Toolbar

Use the buttons in the Display Options toolbar to customize the Data Analytics view. You can choose the plot type, define the number of bins and bar gap, and choose the color of the plot.

❖ To display the Data Analytics View – Display Options toolbar

1. Click a Data Analytics view to select it.
2. Click the **Display Options** toolbar tab.

The Data Analytics View specific display options appear.

Figure 118. Data Analytics View-Display Options

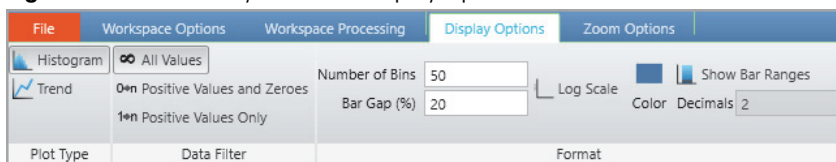


Table 41. Data Analytics View - Display Options toolbar buttons

Button	Description
Plot Type	
Use the parameters in the Plot Type area to choose the plot that the view is displayed.	
Histogram	Displays a histogram plot view in the workspace
Trend	Displays an MS trending plot view in the workspace.
Data Filter	
Use the parameters in the Data Filter area to filter the values.	
All Values	Displays all positive values and zeroes.
Positive Values and Zeroes	Displays only positive values and zeroes.
Positive Values only	Displays only positive value.
Format	
Use the parameters in the Format area to format the display of the view.	
Number of bins	Specifies the number of data groups to display. Default: 50 Ranges: 1-1000 This does not apply to the Categorical data.
Bar Gap	Specifies the width of the gap between bars. The gap is represented as a percentage (%) of the width of the bar. Default: 0-500
Log Scale	Displays the bar in a logarithmic scale. When you click the button, logarithmic changes are applied to y-axis values.
Color	Displays the color palette where you specify the color of the bars in the histogram plot.
Show Bar Ranges	Displays the ranges of data represented by bars in the histogram plot.
Decimals	Specifies the required number of decimal places for the bar range values.

Saving a Data Analytics Plot

You can modify and save the histogram or trend plots in the layout.

❖ **To save a customized Data Analytics View in the layout**

1. Select the Data Analytics view that you want to save.
2. Click the **Workspace Options** toolbar tab.
3. From the Layouts area, choose the option to save the view.

The Data Analytics view is saved in the layout.

When you apply the saved Data Analytics plot in another view, the position of the bins is displayed as it has been saved in the layout. If the imported layout has the binned distribution of quantity plotted in a range that is not present in the current view, the application displays the default distribution group.

Data Analytics Ranges

Use the Data Analytics Ranges option to create or update a Histogram or Trend plot.

Figure 119. Data Analytics Ranges dialog box

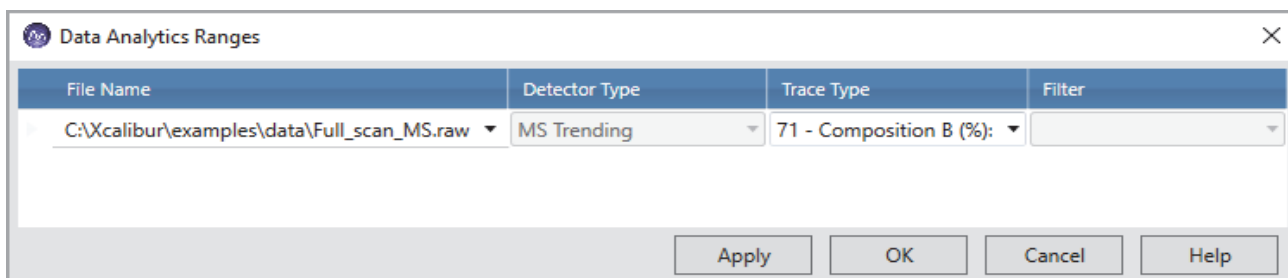


Table 42 describes the columns in the Data Analytics Ranges dialog box.

Table 42. Data Analytics Ranges dialog box

Column	Description
File Name	Specifies the location and file name of the raw data file with the MS Trending data. To import a new raw data file, use the browse button to select a different file.
Detector Type	Specifies the detector used to acquire the data. By default, the detector type is MS Trending.

Table 42. Data Analytics Ranges dialog box

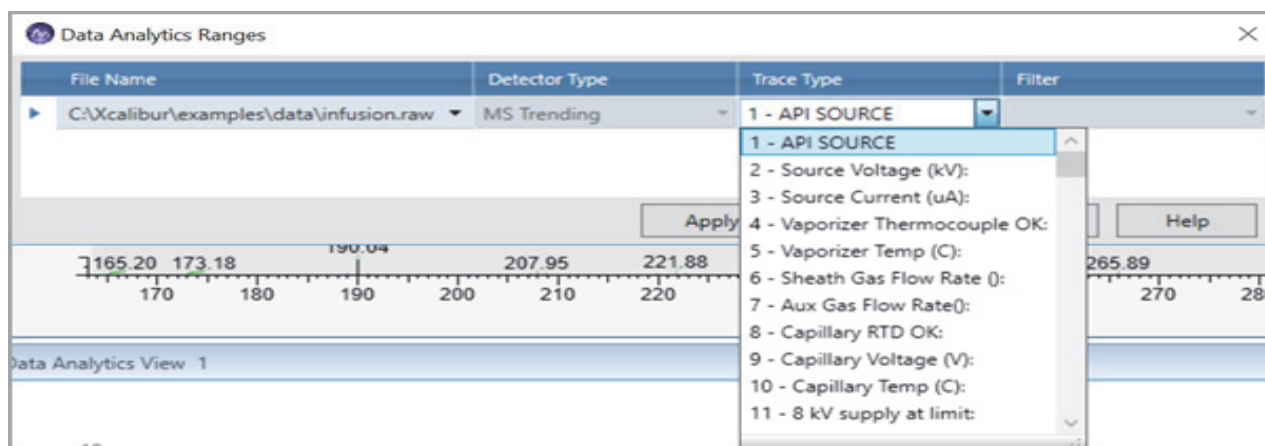
Column	Description
Trace Type	Specifies the trace type. For MS Trending, select one of the instrument status parameters, such as the API source.
Filter	Specifies the scan filter (same selections as the Scan Filters page of the Info bar). You can apply scan filters to filter the Trailer Extra information and Custom trace scan time values. The Status trace type is not associated with scans, and you cannot apply filters to the Status trace type.

Filtering Data Analytics Views by Trace Types

You can create data analytics plots based on selected trace types. The different trace type categories in FreeStyle are as follows:

- Status
- Trailer Extra
- Custom

Figure 120. Data Analytics Plot Trace Types



Using the Status Trace Type

The first group in the Trace Type dropdown list is the Status trace type that displays the MS detector status readback that was collected over time. You cannot apply filters to Status trace types. It contains information about the different status of instrument, such as Temperature, Pressure, Turbo Pump speed, and so on.

6 Analyzing MS Data using Data Analytics View

Data Analytics Ranges

For example, [Figure 121](#) shows the Trend plot created by the drugs_06 raw data file when the MS Trending detector and Ambient Temperature trace type are selected. The plot displays highest and lowest temperatures recorded between 28.4 °C and 28.55 °C due to the temperature variations in the instrument.

Figure 121. Trends plot: Ambient Temperature

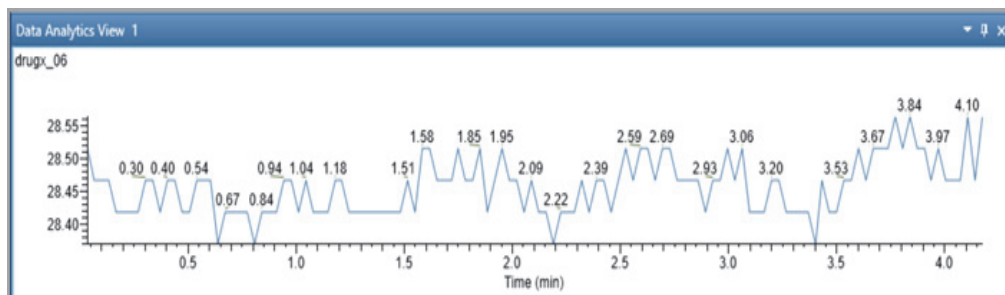
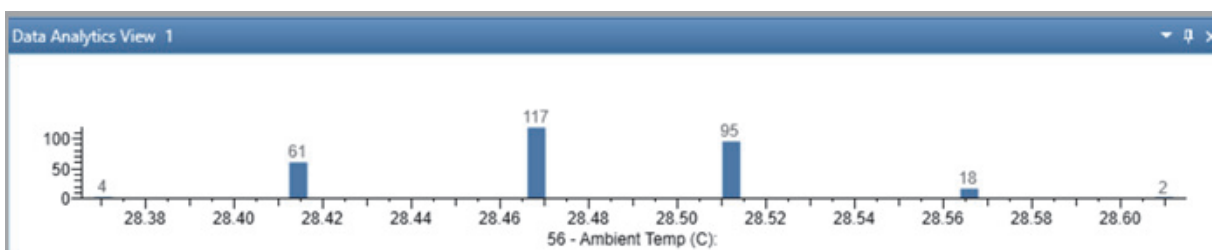
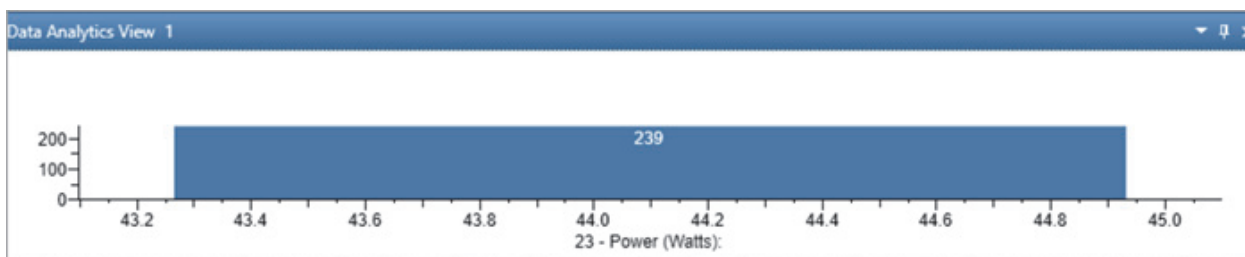


Figure 122. Histogram Plot: Ambient Temperature



Similarly, in [Figure 123](#), the raw data file displays a histogram plot with one bar when Power is selected. The *x* axis of the plot shows that the power remains static between 43.5 and 44.9.

Figure 123. Histogram Plot: Power



Using Trailer Extra

The scan header displays important information about a scan, such as retention time, and depends on the data acquisition settings for the mass spectrometer used to acquire the raw data file. The detector can record additional information about each scan called Trailer Extra. You can apply scan filters to filter the trailer extra information of each scan.

Figure 124. Data Analytics Plot: Number of Ions

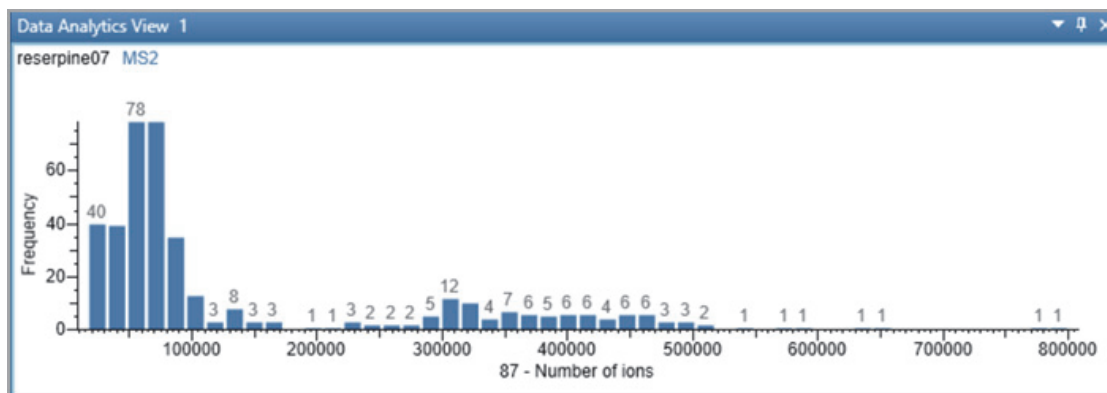


Figure 124 displays an MS2 scan filtered by number of ions.

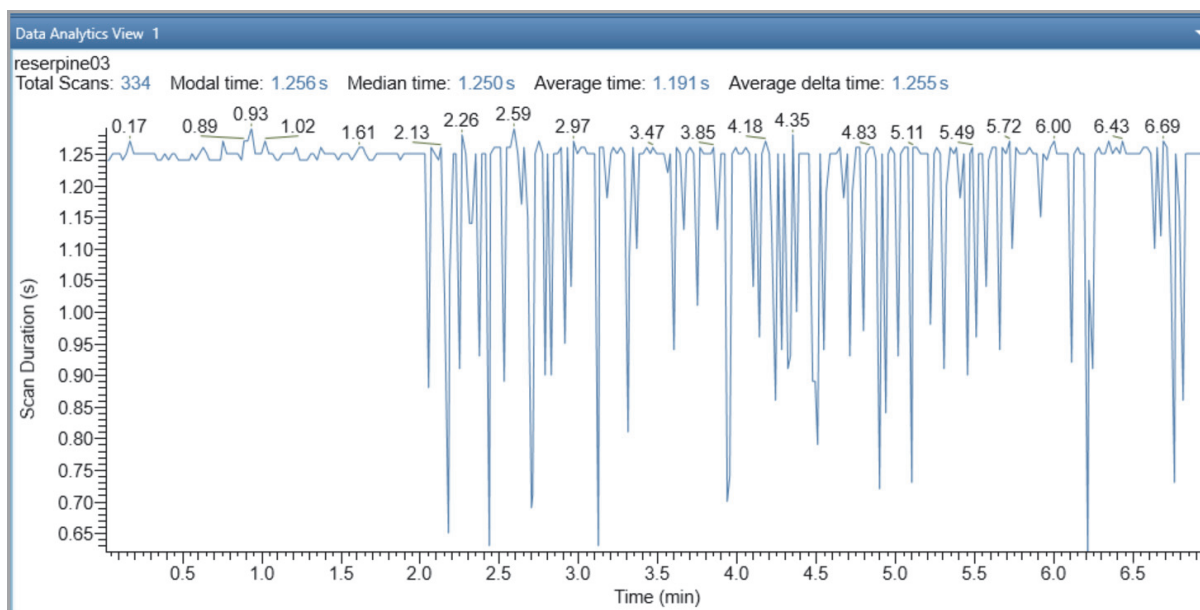
Using Custom Trace Type

The application includes the following Custom Trace types to filter an MS Trending scan:

- Scan Time
- MS Order
- Cycle Time

Scan Time: The amount of time required to accomplish one scan, from the lowest mass to the highest mass of a specified scan range.

Figure 125. Scan Time



If a Scan Time Trend plot is created, the scan header displays the total number of scans, modal time, median time, average time, and average delta time for the selected data.

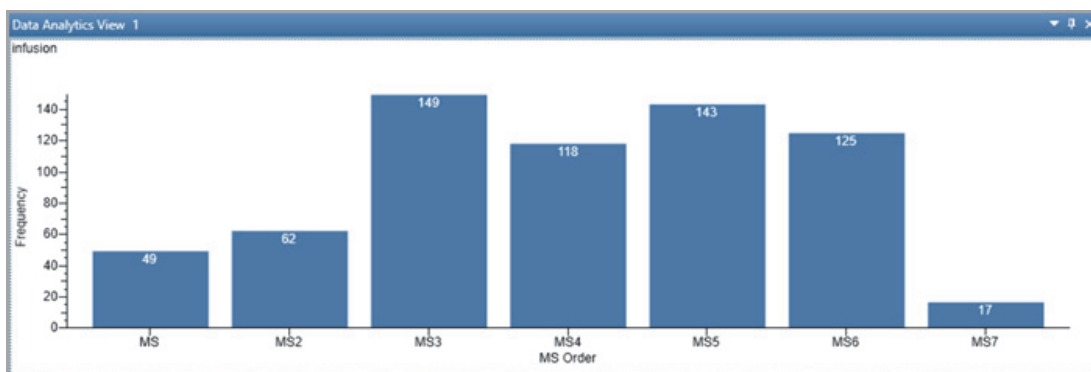
Table 43. Scan Time Trend Plot

Median time	The median is a simple measure of central tendency. Median is the middle value.
Average Time	The average time taken per scan.
Average Delta Time	Average of the time taken between two consecutive MS scans.

MS Order: The MS order of the mass spectra that makes up the chromatogram. It indicates the order of the MS scan.

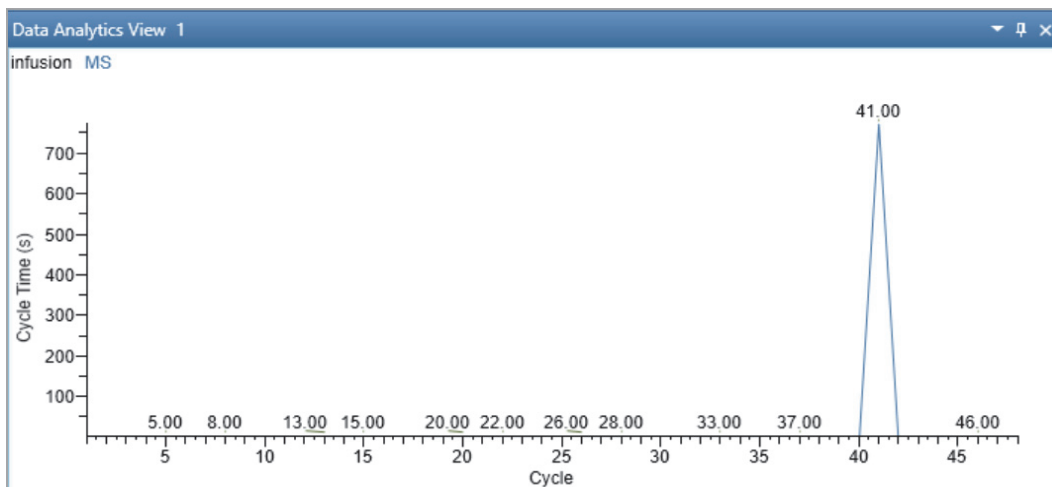
- MS1 is one MS scan.
- MS2 is an MS/MS scan.
- MS3 is an MS3 scan.

Figure 126. Data Analytics: MS Order Trace Type



Cycle Time: Cycle time is the time between two consecutive full MS scans. One cycle is the start of one full MS scan to the start of the next/successive full MS scan.

Figure 127. Data Analytics: Cycle time trace type



Creating a Data Analytics Plot

❖ To display a new plot for MS trending detector type

1. Click a **Data Analytics View** to make it active.
2. From the Workspace Options menu bar, click **Data Analytics Ranges**.

The Data Analytics Ranges dialog box opens with a list of the scans currently displayed in the selected view.

3. In the Data Analytics Ranges dialog box, from the File Name list, click the dropdown arrow to select a file name or click the browse button to select a different raw data file.

In the Detector Type list, the **MS Trending** type is selected by default. You cannot edit the detector type.

4. From the Trace Type dropdown list, select the trace type. For MS Trending, select one of the instrument status parameters, such as the Ion Injection Time (ms).
5. In the Filter box, type or select a different scan filter from the dropdown list that displays the filter options stored in the raw data file.
6. Click **Apply**, and then click **OK**.


Determining the Elemental Composition of Ions

To determine the elemental composition of ions, follow these topics.

Contents

- [Overview of an Elemental Composition Analysis](#)
- [Starting an Elemental Composition Analysis](#)
- [Using the Fragments Matching Algorithm for Confirmation](#)
- [Reviewing the Elemental Composition Results](#)
- [Modifying the Elements in the Elements in Use Table](#)
- [Using Custom Periodic Table for Analysis](#)
- [Elemental Composition Results View](#)
- [Elemental Composition Page](#)

Note In the FreeStyle application, you can add formula labels to the peaks in a mass spectrum, or you can run an elemental composition analysis for a specific mass peak.

- Clicking Elemental Composition in the Workspace Processing toolbar opens the Elemental Composition page in the Info Bar.
- Clicking Elemental Composition in the Spectrum – Display Options toolbar, after you run the elemental composition in the Info Bar, adds formula labels to the peaks in a selected spectrum plot.
- Clicking the Elemental Composition icon, , in the Spectrum view toolbox runs an elemental composition analysis for the selected mass peak, and opens the Elemental Composition page in the Info Bar.

Overview of an Elemental Composition Analysis

The FreeStyle application assigns chemical formulas to components by using an isotopic pattern-matching algorithm that accounts for isotope accurate mass and intensity ratios.

7 Determining the Elemental Composition of Ions

Overview of an Elemental Composition Analysis

The basic elemental composition algorithm uses a single mass, usually the monoisotopic mass of a measured isotope pattern, to calculate all possible elemental compositions that lie within a tolerance window. Then, the algorithm calculates a theoretical isotope pattern for each elemental composition candidate. It calculates the fit between the theoretical and measured isotope pattern, sorting the identified candidates in decreasing order of isotopic pattern score. The isotopic pattern score value is a number between 0 percent (where the patterns are completely different) and 100 percent (where the patterns are indistinguishable by using the scoring parameters specified in the processing method).

The additional fragments matching algorithm provides confirmation of the chemical formula by comparing the fragmentation scan of the selected precursor mass to the theoretical fragmentation pattern for the chemical formula.

To determine the elemental composition of a specific ion, you select the ion of interest in a spectrum of interest, and then start the elemental composition analysis from the Workspace Processing toolbar or the Spectrum view toolbox. To increase your confidence in the best matching formula, you can refine the analysis by adding the information from a fragmentation scan for the selected ion.

When an elemental composition analysis finishes, the application displays the following:

- The [Elemental Composition Results View](#), which lists the best matching chemical formulas.
- The [Elemental Composition Page](#) of the Info Bar, where you can modify the parameters that the elemental composition algorithm uses.
- A theoretical spectrum for the best matching formula in the Spectrum view or MultiSpectrum view. The peaks in the theoretical spectrum are color-coded. Peaks above the specified intensity threshold are green, and peaks below this threshold are red.

[Figure 128](#) shows the Elemental Composition page and the Spectrum view with the theoretical isotope pattern for the calculated elemental composition.

Figure 128. Elemental Composition page, Spectrum view, and Elemental Composition Results view



Starting an Elemental Composition Analysis

After you select a spectrum of interest, you can start an elemental composition analysis from the Workspace Processing toolbar or the Spectrum view toolbox.

- When you start the analysis from the toolbar, you can set up the analysis before the application processes the data.
- When you start the analysis from the toolbox, the application immediately runs the analysis on a selected ion by using the default parameter settings.

To start an elemental composition analysis, follow one of these topics:

- [Starting an Elemental Composition Analysis from the Toolbar](#)
- [Starting an Elemental Composition Analysis from the Toolbox](#)

Starting an Elemental Composition Analysis from the Toolbar

❖ To start an elemental composition analysis from the Workspace Processing toolbar

1. In the Workspace, click a Spectrum view to select it, or select the spectrum of interest in a MultiSpectrum view.

For information about selecting a spectrum of interest, see [Chapter 4, “Reviewing Spectral Data.”](#)

7 Determining the Elemental Composition of Ions

Starting an Elemental Composition Analysis

2. Click the **Workspace Processing** toolbar tab. Then, in the Elemental Analysis area, click **Elemental Composition**.

The Elemental Composition page opens in the Info Bar. The Mass box at the top of the page is populated with the m/z value for the selected spectrum's base peak.

3. (Optional) To select another peak, double-click its m/z label, and then confirm the value in the Mass box.
4. (Optional) To modify the settings for the advanced parameters, click **Show Advanced Parameters**.
5. (Optional) To use a custom periodic table, select the **Use Custom Periodic Table** check box, and then click **Custom Table**.
6. (Optional) Add the fragments matching algorithm to the analysis (see [Using the Fragments Matching Algorithm for Confirmation](#)).
7. Click **Calculate**.

The application runs the elemental composition algorithm on the selected m/z value.

Starting an Elemental Composition Analysis from the Toolbox

❖ To start an Elemental Composition Analysis from the Spectrum view toolbox

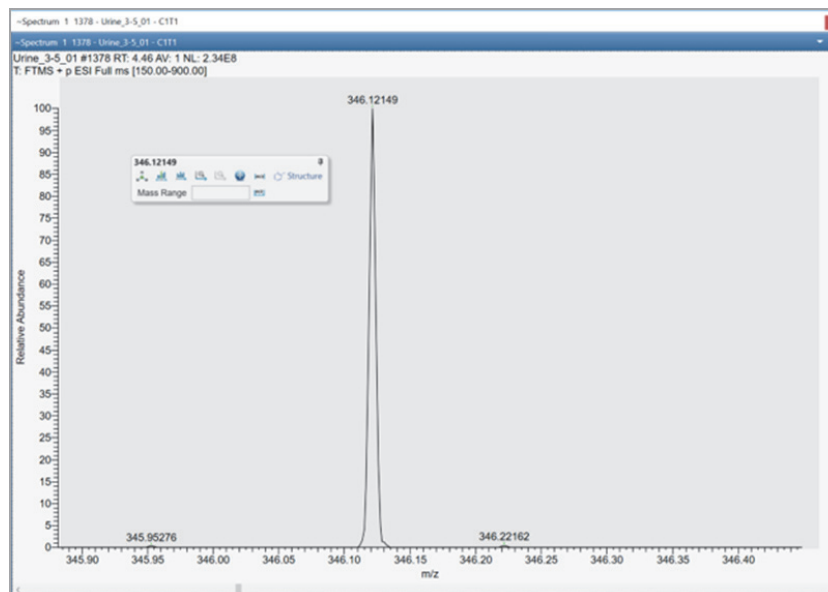
1. In the Workspace, click a Spectrum view to select it, or select the spectrum of interest in a MultiSpectrum view.

For information about selecting a spectrum of interest, see [Chapter 4, "Reviewing Spectral Data."](#)

2. Open the Spectrum toolbox (see [Spectrum Toolbox](#)).
3. In the Spectrum view, place the pointer over the m/z value of interest.

A red rectangle highlights the selected m/z value ([Figure 129](#)).

Figure 129. Spectrum view with a selected m/z value



4. Click the **Elemental Composition** icon, .

The application opens the Elemental Composition page in the Info Bar, populates the Mass box with the m/z value of the selected ion, runs the elemental composition algorithm, and displays the results in the [Elemental Composition Results View](#).

5. (Optional) Restart the analysis by doing one of the following:
- Select another m/z value in the experimental spectrum, and then click the **Elemental Composition** icon.

—or—

- Change the parameter settings on the Elemental Composition page, and then click **Apply**.

For example, you can add the fragments matching algorithm to the analysis (see [Using the Fragments Matching Algorithm for Confirmation](#)).

Using the Fragments Matching Algorithm for Confirmation

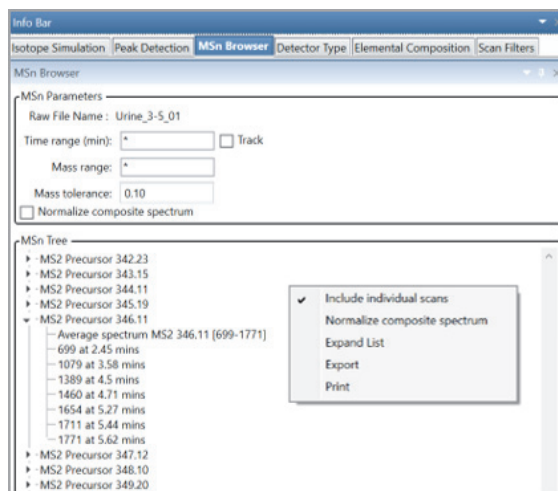
Use the Fragments Matching area of the Elemental Composition page to specify the fragmentation scan number for the fragments matching algorithm. From the Workspace Processing toolbar, you can specify the fragmentation scan before you start the analysis. From the Spectrum view toolbox, you can apply the fragments matching algorithm after the initial analysis finishes.

❖ To use the fragments matching algorithm

1. Find an appropriate fragmentation scan by doing one of the following:
 - a. In the Info Bar, click the **MSn Browser** tab.
 - b. Expand the MS2 node for the precursor mass that you selected on the Elemental Composition page.
 - c. Right-click the MSn Browser page, and choose **Include Individual Scans**.

Figure 130 shows the scan numbers of the individual scans for the MS2 precursor ion at m/z 346.12.

Figure 130. MSn Browser page showing the expanded list for m/z 346.11



—OR—

- a. In the Chromatogram view, insert an MS2 filtered chromatogram for the selected ion.
 - b. Select the MS2 filtered chromatogram and use the left and right arrows to display a data-dependent MS2 scan for the precursor ion in the Spectrum view.
2. On the Elemental Composition page, in the Mass box, check the m/z value of the selected ion.
 3. In the Fragments Matching area, select the **Use Fragments Matching** check box (Figure 131).

Figure 131. Fragments Matching area of the Elemental Composition page

Fragments Matching	
Use Fragments Matching	<input checked="" type="checkbox"/>
Mass Tolerance	5.00 ppm
MSMS Scan No	0
S/N Threshold	3.00
Min. Spectral Fit [%]	10.0

4. In the MSMS Scan No. box, type the scan number that you found in [step 1](#).
5. Click **Apply**.

Review the results in the MSMS Coverage and MSMS Matched Peaks columns of the Elemental Composition Results view.

Reviewing the Elemental Composition Results

After you run an elemental composition analysis, you can review the results in the Elemental Composition Results view.

See these topics:

- [Reviewing the Best-Matching Formulas](#)
- [Reviewing the MS/MS Coverage Score for Matching Fragments](#)
- [Reviewing MS/MS Annotation Results](#)

Reviewing the Best-Matching Formulas

Review the results of an elemental composition analysis in the [Elemental Composition Results View](#). The application displays up to the maximum number of candidates specified on the Elemental Composition page or the Default Elemental Composition page and orders the best matching formulas by rank.

❖ To review the best-matching formulas

1. For each row in the Elemental Composition Results view, review the formula, number of matched isotopes, and MS coverage score.
2. Compare the theoretical isotope pattern for the proposed chemical formula, and then zoom in on the mass peak of interest in the Spectrum view.

In the theoretical isotope pattern, a green marker represents the matched mass peaks that fall inside the spectral distance and a red marker represents the missing mass peaks that fall outside the spectral distance (mass tolerance plus intensity tolerance) at the experimental resolution. An Orange marker represents the matched mass peaks that fall outside the spectral distance in the Spectrum view.

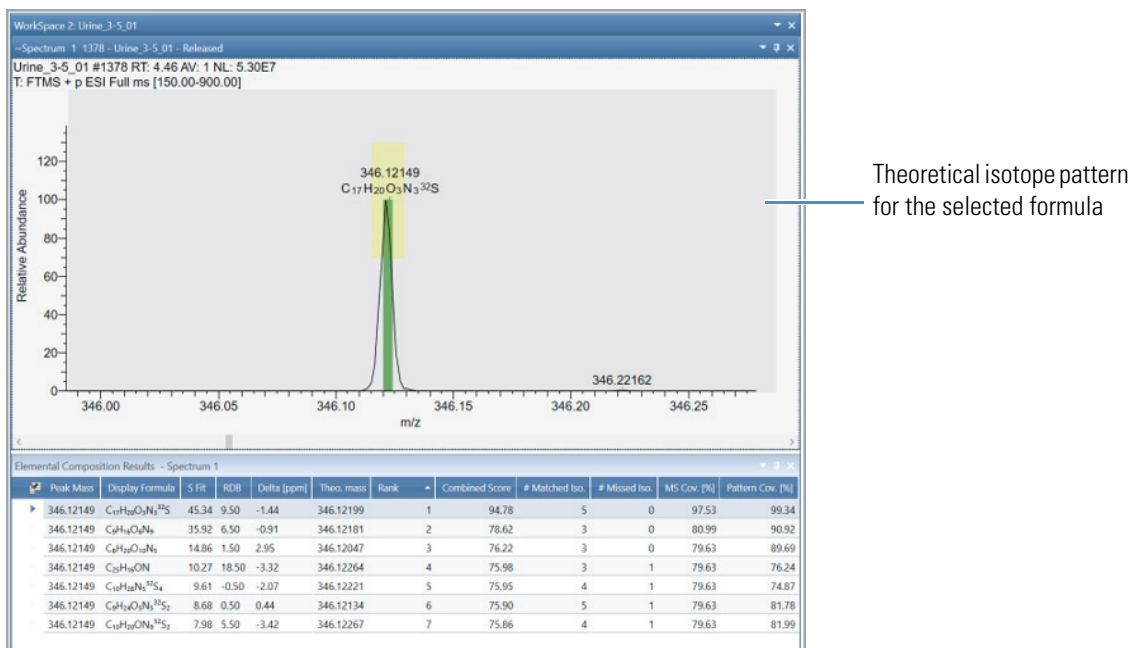
7 Determining the Elemental Composition of Ions

Reviewing the Elemental Composition Results

The isotope pattern uses the resolution provided with the scan data to perform the isotope pattern and calculate the spectral distance.

Figure 132 shows the elemental composition results for a mass peak at m/z 346.12.

Figure 132. Theoretical isotope pattern with the markers



You can view the chemical formula of each mass spectrum peak by clicking Elemental Comp. button in the Spectrum Display Options toolbar. You can also view the theoretical mass and delta mass units of the simulated isotope profile when you point to the isotope markers.

Note The data that displayed when you point to the isotope markers is different from the Elemental Composition Results table. The results table displays the theoretical and delta mass units for the monoisotopic peaks; pointing to the markers display data for the simulated profile.

Reviewing the MS/MS Coverage Score for Matching Fragments

After you run the elemental composition algorithm with the additional fragments matching algorithm (see [Using the Fragments Matching Algorithm for Confirmation](#)), review the results in the MS/MS Coverage [%] and MS/MS Matched Peaks columns of the Elemental Composition Results table.

❖ To compare the number of fragmentation peaks to the value in the MS/MS Matched Peaks column

1. In a Spectrum view, display the fragmentation scan that you entered on the Elemental Composition page, in the Fragments Matching area, in the MS/MS Scan No box.

2. Compare the values for MS/MS Coverage [%] and MS/MS Matched Peaks to the number of mass spectrum peaks in the fragmentation scan.

Reviewing MS/MS Annotation Results

After you perform the elemental composition, you can review the results of an MS/MS scan in the MSMS Annotation result window. The application displays the result when you type the spectral scan corresponding to the MS/MS scan number in the info bar.

❖ To review the MS/MS annotation results

1. The spectral scan corresponding to MS/MS scan number opens only when you enable the **Use fragment matching** option and type a valid MS/MS scan number in the info bar.
2. When you change the MS/MS scan number and re-perform the elemental composition analysis, the result of the MS/MS scan is automatically replaced in the MS/MS Annotation result window.
3. After completing the Elemental composition analysis, the MS/MS Annotation result window is annotated with the elemental composition of the fragment masses.

Note Zoom in on the MS/MS Annotation Result page to view the annotation of the lower intensity peaks.

Modifying the Elements in the Elements in Use Table

The elemental composition algorithm uses only the minimum and maximum number of atoms for each element in the Elements in Use table to determine the elemental composition of ions. You can edit the Elements in Use table in two locations: the Elemental Composition page (in the Info Bar) or the Default Elemental Composition page (in the Default Options Configuration dialog box).

To add elements to or remove elements from the table, follow the appropriate topic:

- [Adding Elements to the Elements in Use Table](#)
- [Elemental Composition Results View](#)

Adding Elements to the Elements in Use Table

Use the Periodic Table of Elements dialog box to add isotopes to the Elements in Use table on the Elemental Composition page or Default Elemental Composition page.

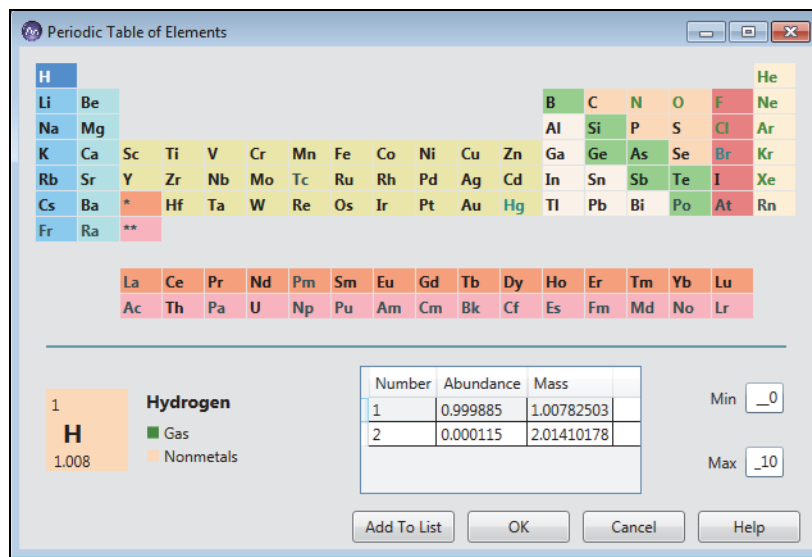
❖ To add an element to the Elements In Use table

1. To display the dialog box of the periodic table of elements, on the Elemental Composition page or the Default Elemental Composition page, click **Add** below the Elements in Use table.

7 Determining the Elemental Composition of Ions

Modifying the Elements in the Elements in Use Table

Figure 133. Periodic table of elements



2. Do the following for each element that you add:
 - a. Click the element's symbol in the table.

The element's isotopes appear below the periodic table.
 - b. If the selected element has several isotopes, select the isotope with the highest abundance for non-labeled analytes or the appropriate isotope for labeled analytes.
 - c. Specify the limits for the element as follows:
 - In the Min box, type an integer to specify the minimum number of atoms of this element that must be in the calculated elemental composition.
 - In the Max box, type an integer to specify the maximum number of atoms of this element that can be in the calculated elemental composition.
 - d. Click **Add to List**.
3. Click **OK** to add the selected elements to the table.

Removing Elements from the Elements In Use Table

❖ To remove an element from the Elements In Use Table

1. On the Elemental Composition page of the Info Bar or the Default Elemental Composition page of the Default Configuration Options dialog box, below the Elements in Use table, click to the left of the element to highlight its row (Figure 134).

Figure 134. Elements in Use area on the Elemental composition page of the Info Bar

Elements in Use				
Isotope	Mass	Min	Max	
14 N	14.0031	0	10	
16 O	15.9949	0	15	
12 C	12.0000	0	30	
1 H	1.0078	0	60	
32 S	31.9721	0	10	
35 Cl	34.9689	0	4	
31 P	30.9738	0	10	Selected row

Remove Add

2. Click **Remove**, and in the confirmation box, click **OK**.

Using Custom Periodic Table for Analysis

Use the Custom Periodic table for the elemental composition calculation by adding custom isotope abundance to the elements. The table provides a list of all known isotopes, including those having zero abundance. When you select an element from the list, the elemental composition details of the selected element display separately. You can edit the abundances of each isotope and change the total abundance of the element. The abundance of each isotope must be greater than 0 and the total abundance must be equal to 1. You can save the changes and also load an existing custom periodic table.

Figure 135. Custom Periodic Table

Custom Periodic Table of Elements

C:\Users\julie.erinjery\Desktop\JulieE_custom.xml

Number	Abundance	Mass
<input checked="" type="checkbox"/> 14	0.99632	14.003074
<input checked="" type="checkbox"/> 15	0.00368	15.0001089

Revert to Default Total Abundance: 1

Save Save As Load Cancel Help

❖ To add custom isotope abundance

1. On the Elemental Composition Info Bar, below Custom Periodic Table, select the **Use Custom Periodic Table** check box.
2. Click **Custom Table**.
3. Click the element's symbol in the table.

7 Determining the Elemental Composition of Ions

Using Custom Periodic Table for Analysis

The element's isotopes appear below the periodic table.

4. Select the isotope and edit the abundance.
5. Click **Save**.

The Elemental Composition Info page displays the name and path of the loaded custom periodic table.

Note The application highlights the Total Abundance value column in a red rectangle if the abundance is greater or less than 1.

6. Click **Revert to Default** to revert and use the default periodic table.
7. Click **Load** to use an existing custom periodic table saved in your computer.

In the following example, when N is selected from the Element list, the isotopes 14 and 15 is displayed below with an abundance of 0.99632 and 0.00368 respectively. The Total Abundance of the isotopes is 1.

Figure 136. Total Abundance equals to 1

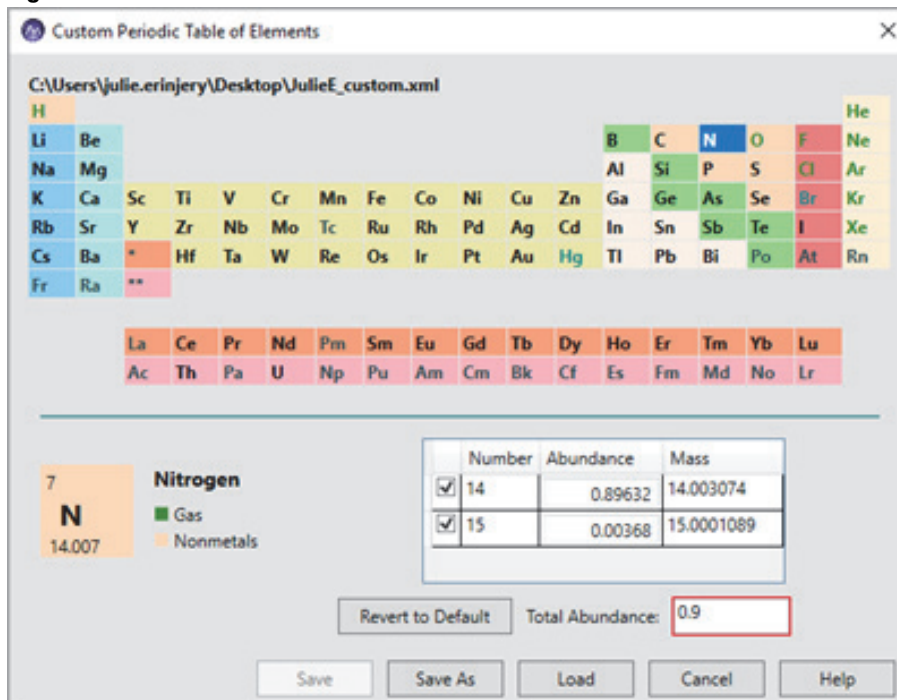
The screenshot shows a window titled "Custom Periodic Table of Elements" with the file path "C:\Users\julie.erinjery\Desktop\JulieE_custom.xml". The periodic table is displayed with Nitrogen (N) highlighted in blue. Below the table, the element information for Nitrogen is shown: "7 Nitrogen", "N", "14.007", "Gas", and "Nonmetals". A table of isotopes is displayed:

Number	Abundance	Mass
<input checked="" type="checkbox"/> 14	0.99632	14.003074
<input checked="" type="checkbox"/> 15	0.00368	15.0001089

Below the table, there is a "Revert to Default" button and a "Total Abundance:" field with the value "1". At the bottom, there are buttons for "Save", "Save As", "Load", "Cancel", and "Help".

When the abundance of isotope 14 is changed to 0.89632, the Total Abundance becomes 0.9 and the value is highlighted in a red rectangle as the abundance is less than 1.

Figure 137. Total Abundance less than 1



When you run an elemental composition analysis using the custom periodic table, the application displays the matching chemical formula for the selected peak in the spectrum based on the modified abundance. The application uses default periodic table if the custom periodic table is not selected for analysis.

Elemental Composition Results View

After you run an elemental composition analysis, the Elemental Composition Results view appears with the best matching calculated chemical formulas for the selected peak in the spectrum.

Table 44 describes the columns in the Elemental Composition Results view. By default, the S Fit, RDB, Combined Score, MSMS Cov. [%], MSMS Shift Measure, and MSMS Matched Peaks columns are hidden. To display these columns, use the Field Chooser dialog box.

Table 44. Elemental Composition Results view parameters (Sheet 1 of 3)

Parameter/icon	Description
Field Chooser icon (☰)	Opens the Field Chooser dialog box for selecting the result columns to display.
Peak Mass	Displays the mass-to-charge ratio of the selected isotopic peak.

Table 44. Elemental Composition Results view parameters (Sheet 2 of 3)

Parameter/icon	Description
Display Formula	Displays the chemical formula of the match by using the elements that you specified on the Elemental Composition Page . Also displays isotope labels for elements listed in the Elements in Use table.
S Fit	Displays the spectral similarity score between the measured and theoretical isotope patterns. Fragments matching does not affect this value.
RDB	Displays the ring and double-bond (RDB) equivalents that the algorithm calculated for the proposed match.
Delta (ppm)	Displays the difference between the measured mass-to-charge ratio and the theoretical mass-to-charge ratio. The table lists only formulas whose mass is within the tolerance that you specified on the Elemental Composition Page , in the units (ppm, mmu, or amu) that you specified there.
Theoretical Mass	Displays the theoretical mass-to-charge ratio of the proposed match.
Rank	Displays the ranking of the proposed match by decreasing Combined Score values.
Combined Score	Displays a percentage that conveys how close the measured spectrum matches the theoretical spectrum. Fragments matching refines this value by including the relative number of matching fragments in the score. The Combined Score for a proposed formula increases if the fragmentation spectrum includes more matching fragments and decreases if the fragmentation spectrum includes fewer matching fragments relative to other formulas.
#Matched Isotopes	<p>Displays the number of isotopes in the measured isotope pattern that match the theoretical isotope pattern for the chemical formula.</p> <p>A matching isotope matches the delta mass from the A0 peak (the peak with the lowest m/z value in an isotope pattern) and the relative intensity of the theoretical isotope pattern within the specified tolerances.</p>
#Missed Isotopes	Displays the number of isotopes that are missing in the measured isotope pattern for the precursor ion. An isotope peak is missing if it is part of the theoretical isotope pattern and not found in the measured spectrum within the specified mass and intensity tolerances.

Table 44. Elemental Composition Results view parameters (Sheet 3 of 3)

Parameter/icon	Description
MS Coverage [%]	<p>Displays the summed intensity of matching isotope peaks in the measured pattern relative to the summed intensity of all the peaks in the measured pattern.</p> $\frac{\text{Summed intensity of the matching isotope peaks} \times 100}{\text{Summed intensity of all the peaks in the measured pattern}}$ <p>IMPORTANT Low values for all the candidates might indicate an overlapping pattern rather than a lack of good matches.</p>
Pattern Coverage [%]	<p>Displays the summed intensity of the matching isotope peaks in the measured MS1 spectrum relative to the summed intensity of the theoretical isotope pattern.</p> $\frac{\text{Summed intensity of the matching isotope peaks} \times 100}{\text{Summed intensity of the theoretical isotope pattern}}$ <p>Provides a quantitative measure of how well the measured isotope pattern matches the theoretical isotope pattern.</p> <p>Note Because the A0 peak (peak with the lowest m/z value in the pattern) is typically responsible for most of the pattern intensity, even a small decrease in the percent coverage might be important. For example, a missing peak for an isotope with two ^{13}C atoms might cause only a small decrease in the summed intensity of the measured isotope pattern.</p>
The following columns display nonzero values when you add the fragments matching algorithm to the elemental composition analysis.	
MSMS Cov. [%]	<p>Displays the summed intensity of the matched fragment peaks relative to the summed intensity of all the fragment peaks in the selected MSMS scan.</p> <p>Note Low values for all the candidates might indicate a contaminating compound within the isolation window for the fragmentation scan.</p>
MSMS Shift Measure	<p>Not used for scoring.</p> <p>The application searches for each fragment in a range from the expected mass based on the chemical formula and the expected mass plus the delta mass between the precursor ion and its displayed formula. The MSMS Shift Measure value increases when the fragments are closer to their expected masses rather than their shifted masses.</p>
MSMS Matched Peaks	Displays the number of matching fragments in the MSMS spectrum. Click the expand icon, +, to view the fragment list.

Elemental Composition Page

Use the Elemental Composition page (Figure 138) of the Info Bar for specifying the criteria to calculate the best matching chemical formula for a mass-to-charge ratio in a mass spectrum.

Figure 138. Elemental Composition page

The screenshot shows the 'Elemental Composition' panel in the software's Info Bar. It is divided into several sections for configuring the calculation of a chemical formula based on a mass-to-charge ratio.

Mass

Single mass Selected scan

Mass: 420.411071777344

Calculate

Prediction Settings

Mass Tolerance: 5.00 ppm

Max. # Candidates: 10

Charge: 1

Fragments Matching

Use Fragments Matching

Mass Tolerance: 5.00 ppm

MSMS Scan No: 0

Custom Periodic Table

Use Custom Periodic Table **Custom Table**

C:\Users\Standard\AppData\Local\Thermo Scientific\Freestyle\Templates\CustomElements.xml

Elements in Use

Isotope	Mass	Min	Max
14 N	14.0031	0	10
16 O	15.9949	0	15
12 C	12.0000	0	30
1 H	1.0078	0	60
32 S	31.9721	0	10
35 Cl	34.9689	0	4
31 P	30.9738	0	10

Remove Add

Simulated Spectrum

Profile Resolution: 1000000

Load Save As Apply Help

Save as Defaults

Table 45. Elemental Composition page parameters (Sheet 1 of 6)

Parameter	Description
Show/Hide Advanced Parameters	The advanced parameters to show or hide are as follows: <ul style="list-style-type: none"> • Under Prediction Settings: Min and Max RDBE, Nitrogen Rule, Centroid Algorithm, Normalization Mode, and Use Representative Elements • Pattern Matching: Intensity Threshold and Intensity Tolerance • Under Fragments Matching: S/N Threshold and Min. Spectral Fit
Mass	
Single Mass	Specify the mass-to-charge ratio that the algorithm uses to calculate probable chemical formulas. Range: Minimum m/z of the measured spectrum to the maximum m/z of the measured spectrum
Selected Scan	Specify the selected scan that the elemental composition calculation is performed. The spectrum plot and each row of the spectrum list displays the highest scoring result for each peak. The results view is not available for this analysis.
Calculate	Calculate formulas and display them in the Elemental Composition Results View .
Prediction Settings	
Mass Tolerance	Specify a mass tolerance to restrict the number of possible elemental compositions. The elemental composition algorithm returns results of the search only if the theoretical mass matches the submitted mass within the specified tolerance. Range: 0.00–100.00; default: 5.00 ppm
Units	Select the units that you want to associate with the mass tolerance. The options are amu (atomic mass units), mmu (millimass units), and ppm (parts-per-million). If you specify the error limits in ppm, the errors are mass-dependent and get larger at higher masses and smaller at lower masses. Default: ppm
Max #Candidates	Specify the maximum number of formulas to display. Range: 0–400; default: 10

Table 45. Elemental Composition page parameters (Sheet 2 of 6)

Parameter	Description
Charge	Select the charge state that you want to use to calculate the probable formulas. Range: -99 to 99; default: 1
Min. and Max. RDBE	Specify a range of values for double bonds and ring equivalents—a measure of the number of unsaturated bonds in a compound—that limits the calculated formulas to only those that make sense chemically. Limits range: -1000.0 to 1000.0 The value is calculated by the following formula: $D = 1 + \frac{\left[\sum_i^{i_{\max}} N_i (V_i - 2) \right]}{2}$ where: <ul style="list-style-type: none">• D is the value for the RDB equivalents• i_{\max} is the total number of different elements in the composition• N_i is the number of atoms of element i• V_i is the valence of atom i The calculation produces an integer such as 3, which indicates an odd-electron ion, or a number with a remainder of 0.5, which indicates an even-electron ion. Minimum value: -0.5, corresponding to a protonated, saturated compound

Table 45. Elemental Composition page parameters (Sheet 3 of 6)

Parameter	Description
Nitrogen Rule	<p>Select whether or how to use the Nitrogen Rule in the formula calculation:</p> <ul style="list-style-type: none"> • Do Not Use: Do not use the Nitrogen Rule. • (Default) Even Electron Ions: Select for even-electron ions, such as protonated species. • Odd Electron Ions: Select for odd-electron ions, such as radical cations. <p>Note McLafferty³ states the Nitrogen Rule as follows: “If an odd-electron ion contains no (or an even number of) nitrogen atoms, its molecular ion will be at an even mass number...[Similarly,] an odd-electron ion will be at an odd mass number if it contains an odd number of nitrogen atoms.”</p>
Centroid Algorithm	<p>Select the centroiding algorithm for profile data.</p> <p>Default: FT Orbitrap Selections: FTOrbitrap, GCQ, TSQ, or MAT</p>
Normalization Mode	<p>Select one of the following normalization modes for the spectral peak intensities:</p> <ul style="list-style-type: none"> • (Default) Base peak: The most common normalization mode. Normalizes the isotope peaks to the base peak height of 100%. Its disadvantage is the propagation of the intensity error of the base peak to all isotope intensities. • Linear: Normalizes the theoretical and measured patterns so that the sum of all isotopic pattern intensities is the same. • Quadratic: Normalizes the theoretical and measured patterns so that the error squares (intensity differences) are minimized.

Table 45. Elemental Composition page parameters (Sheet 4 of 6)

Parameter	Description
Use Representative Elements	<p>Select the check box to use the Representative Elements table for the elemental composition calculations. Clear the check box to use the Protein Elements table.</p> <p>Currently the elemental composition algorithm supports two types of element tables, differing mainly in their carbon (C) abundances.</p> <ul style="list-style-type: none"> The Representative Elements table has the following C abundances: <ul style="list-style-type: none"> 12C: 0.9893 13C: 0.0107 The Protein Elements table has the following C abundances: <ul style="list-style-type: none"> 12C: 0.989136445 13C: 0.010863555 <p>Default: Selected</p>
Pattern Matching	
Intensity Threshold [%]	<p>Specify the isotope intensity threshold, relative to the base peak of the theoretical isotope pattern that the algorithm uses for pattern simulation. The algorithm skips isotopes below the threshold—that is, if the expected intensity of an isotopic peak is below the threshold, the algorithm does not look for the peak in the experimental spectrum, and the peak is not part of the score calculation.</p> <p>Range: 0.00 to 10 000.00%; default: 0.10%</p>
Intensity Tolerance [%]	<p>Specify the relative intensity tolerance for the isotope search.</p> <p>Range: 0.00 to 10 000.00%; default: 30.00%</p>
Fragments Matching	
Use Fragments Matching	<p>Select this check box to turn on the fragment matching algorithm, which ranks the identified candidates (chemical formulas) by the number of matching peaks in the fragmentation scan for the precursor ion.</p> <p>The fragment matching algorithm requires a fragmentation scan for the selected precursor ion.</p> <p>To perform MS/MS matching, you must enter the scan number of the fragmentation scan for the precursor ion (m/z value) of interest.</p>

Table 45. Elemental Composition page parameters (Sheet 5 of 6)


Parameter	Description
Mass Tolerance	Specify the mass tolerance for the peaks in the fragmentation spectrum (expected mass versus theoretical mass).
MSMS Scan No	Specify the MS/MS scan number of the fragmentation scan for the selected precursor ion. For instructions on how to find an appropriate scan number, see Using the Fragments Matching Algorithm for Confirmation . Default: 0 (No fragment matching); range: 1 to the last scan number in the raw data file
S/N Threshold	Specify the signal-to-noise threshold for the peaks in the MS/MS spectrum. The fragments matching algorithm ignores peaks below this S/N threshold.
Min. Spectral Fit [%]	Restrict the candidate list to candidates that meet or exceed the minimum spectral fit value. Range: 0.00 to 100.00%; default: 10.00%
Custom Periodic Table	
Defines elements having customized isotope abundance that you want to consider when the algorithm calculates the elemental composition.	
Use Custom Periodic Table	Use the custom periodic table for elemental composition analysis.
Button	
Save	Saves the modified isotope abundances.
Save As	Saves the custom periodic table to your computer.
Load	Selects an existing custom periodic file that contains the modified isotope abundances to consider for the elemental composition analysis.
Elements in Use	
Lists the isotopes and the number of occurrences for each isotope to consider when the algorithm calculates possible elemental compositions for the submitted mass value. See Modifying the Elements in the Elements in Use Table .	
Field Chooser 	Displays the Field Chooser dialog box for selecting which fields appear in the Elements in Use table (see Selecting the Columns to Display in a View or Dialog Box with Tabular Data).

Table 45. Elemental Composition page parameters (Sheet 6 of 6)

Parameter	Description
Isotope	<p>Select the isotopes that you want the data system to consider when it calculates the possible elemental compositions for a monoisotopic ion with the given mass (A0 mass peak).</p> <p>To add an isotope, click Add. The dialog box for the periodic table of elements opens. You can also right-click in the grid and choose Add from the shortcut menu.</p> <p>To remove an isotope, click to the left of the element in the table to highlight its row, and then click Remove or right-click and choose Remove from the shortcut menu.</p>
Min	Specify the minimum number of occurrences of an isotope in the chemical formula.
Max	Specify the maximum number of occurrences of an isotope in the chemical formula.
Mass	<p>Displays the exact isotopic mass for each isotope in the Elements in Use list.</p> <p>You cannot edit this value.</p>
Simulated Spectrum	
Profile	<p>Select the Profile check box to add an overlay of the simulated spectrum. The Resolution box displays the value used to perform the elemental composition. You cannot edit the resolution value. The scores are calculated based on the peaks determined from the profile.</p> <p>Tip You can view the resolution values for each mass spectrum peak by clicking Peak Resolution in the Spectrum Display Options toolbar (see Spectrum – Display Options Toolbar).</p>
Button	
Load	Select a file (with an .limx file name extension) that contains a set of isotope limits.
Save As	Save a list of isotope limits to a file (with an LIMX file name extension).
Apply	Apply the elemental composition settings to the active mass spectrum.
Help	Opens the FreeStyle Help to the elemental composition topic.

^a McLafferty, F. W. *Interpretation of Mass Spectra*; University Science Books: Mill Valley, CA, 1980.

Searching Mass Spectrum Libraries

To compare a query spectrum against a set of mass spectrum libraries, follow these topics.

Contents

- [Performing a Local NIST or mzVault Library Search](#)
- [Reviewing the Results of a Local NIST or mzVault Library Search](#)
- [Modifying the Settings for a Local NIST or mzVault Library Search](#)
- [Searching the Online mzCloud Mass Spectral Database](#)
- [Exporting a Mass Spectrum to the NIST MS Search Application](#)
- [Managing Libraries](#)
- [Spectrum Workspace Processing Toolbar – Library Search Buttons](#)

Performing a Local NIST or mzVault Library Search

Search your local NIST or mzVault mass spectrum libraries to identify unknown compounds.

Follow these topics in order:

1. [Setting Up the Default Library Search Parameters](#)
2. [Selecting the Query Spectrum](#)
3. [Starting a Library Search](#)

Setting Up the Default Library Search Parameters

Before you can perform an mzVault library search, you must specify the location of at least one of your local mzVault database files. The application automatically links to the NIST libraries installed during the software installation process.

Follow these procedures:

- To open the Library Search page of the Default Options Configuration dialog box
- To specify the location of your local mzVault database files
- To specify the default parameter settings for a library search

❖ To open the Library Search page of the Default Options Configuration dialog box

1. In the Workspace Options toolbar, click **Default Options**.

The Default Options Configuration dialog box opens.

2. In the left pane, click **Library Search**.

❖ To specify the location of your local mzVault database files

1. In the Library Type area on the Library Search page, select the **mzVault** option.
2. For each database file that you want to add to the Search list, do the following:
 - a. In the Search List area, browse to the location of your mzVault database files.
 - b. Select an mzVault database file (DB).
 - c. Click **Open**.
3. Click **Save**.

Tip After you run your first mzVault search, you can also specify the location of your mzVault database files on the mzVault Search page in the Info Bar.

❖ To specify the default parameter settings for a library search

1. In the Library Type area on the Default Library Search page, select either the **NIST** or **mzVault** option.
2. Make the appropriate selections and entries.
3. Click **Save**.

Note If you plan to run both NIST and mzVault searches, make sure to select and save the settings for both search types.

Tip After you run a library search, selecting a different setting from the Display list on the mzVault Search page or the NIST Search page (in the Info Bar) automatically changes how the Spectrum view displays the query spectrum versus the matching library spectrum. You can also modify the search criteria from the Info Bar; however, to start the new search, you must click Apply.

Go to the next topic [“Selecting the Query Spectrum.”](#)

Selecting the Query Spectrum

The query spectrum is the spectrum that you want to search against a library of mass spectra.

❖ To select the query spectrum

1. Open a raw data file that contains mass spectral data.

IMPORTANT You can perform a library search on a mass spectrum, an average mass spectrum, or a composite mass spectrum. For an mzVault or a NIST search for a matching MS/MS spectrum, the query spectrum must be a data-dependent scan.

2. Select a mass spectrum by doing any of the following:

- Select a data point in a chromatogram plot. If a Spectrum or Multi Spectrum view is not open, in the Workspace Options toolbar, click **Spectrum**.

The mass spectrum for the selected time point appears in the Spectrum view or the Multi Spectrum view.

- Click the **MSn Browser** tab, and then double-click an item in the MSn Tree.

The selected spectrum opens in a new Spectrum view.

–or–

- a. Click the Spectrum (or Multi Spectrum) view of interest.
- b. In the Workspace Options toolbar, click **Spectrum Ranges**.
- c. In the Spectrum Ranges dialog box, specify the spectrum of interest.

Note You can perform a library search on a mass spectrum, an averaged mass spectrum, or a composite mass spectrum.

Go to the next topic [“Starting a Library Search.”](#)

Starting a Library Search

After you set up the default options for an mzVault or NIST library search and select a query spectrum, you are ready to start the search.

❖ To start a library search

1. Set up the default search settings (see [Setting Up the Default Library Search Parameters](#)).
2. Select a spectrum plot (see [Selecting the Query Spectrum](#)).
3. Click the **Workspace Processing** tab.

8 Searching Mass Spectrum Libraries

Reviewing the Results of a Local NIST or mzVault Library Search

4. Click **mzVault Search** or **NIST Search**.

The selected library search page appears in the Info Bar, and the application automatically searches the specified libraries.

When the search ends, the following items appear:

- The library spectrum for the best hit appears in the Spectrum view as a stacked plot, a mirror plot, or an exclusion plot.

Note From the Info Bar, you can change the display setting on the mzVault Search page or the NIST Search page without rerunning the search.

- A list of search hits appears in the search results view.
- The structure of the best hit appears in the chemical structure view.

Go to the next procedure “[Reviewing the Results of a Local NIST or mzVault Library Search.](#)”

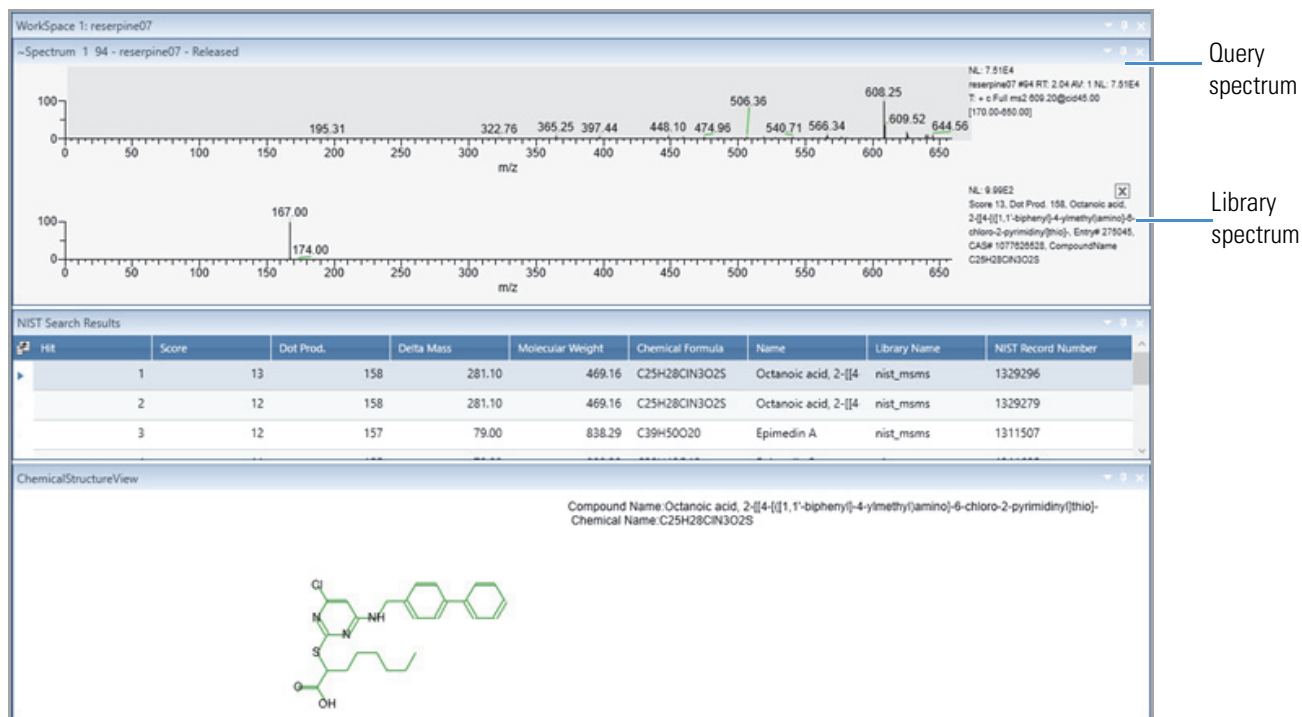
Reviewing the Results of a Local NIST or mzVault Library Search

To review the results of a library search, see the appropriate topics:

- [Reviewing the Results of a NIST Library Search](#)
- [NIST Search Results View](#)
- [Reviewing the Results of an mzVault Search](#)
- [mzVault Search Results View](#)

Reviewing the Results of a NIST Library Search

[Figure 139](#) shows the results of a NIST library search. The Spectrum view contains a mirror plot. The NIST Search Results view lists 10 hits sorted in order of probability. The Chemical Structure View displays the two-dimensional structure of the selected hit.

Figure 139. Results of a NIST library search displayed as a mirror plot

NIST Search Results View

The NIST Search Result view appears in the workspace after you submit a NIST library search. The library search algorithm returns a list of the best matches from the selected libraries. See [Managing Libraries](#).

❖ To display the NIST Search Results view

Perform a NIST library search (see [Performing a Local NIST or mzVault Library Search](#)).

[Table 46](#) describes the columns for the NIST Search Results view.

Table 46. NIST Search Result view columns

Column	Description
Hit	Relative ranking of library search matches based on decreasing SI (Search Index) values.
SI (Search Index)	Direct matching factor for the query spectrum and the library spectrum.
RSI (Reverse Search Index)	Reverse search matching factor that ignores any spectrum peaks in the query spectrum that are not in the library spectrum.
Prob	Probability factor based on the differences between adjacent hits in an SI ordered list.
Molecular Weight	Molecular weight (in daltons) of the library search match.
Chemical Formula	Chemical formula of the library search match.
Name	Name of the matched compound in the library.
Library Name	Name of the library that contains the matching compound.
NIST Record Number	Record number of the matched compound in the NIST library.
Score	Number of results returned from NIST.
Dot. Prod.	Defines the score returned from the NIST library.
Delta Mass	Delta Mass of the matched compound in the library. It is the differences between the experimentally determined mass-to-charge ratios and the theoretical mass-to-charge ratios. Displays delta mass only for the MS/MS Hybrid search type.

Three factors describe the accuracy of the match to the submitted spectrum: SI, RSI, and Prob. With the SI and RSI matching factors, a perfect match results in a value of 1000. As a general guide, 900 or greater is an excellent match; 800–900, a good match; and 700–800, a fair match. A matching factor less than 600 is a poor match. Unknown spectra with many peaks tend to yield lower match factors than similar spectra with fewer peaks.

The probability factor is a complex parameter based on the SI matching factor and the difference between adjacent matches. If a match has an SI match factor greater than 900 and the next best match has a match factor of 300, the probability of the compound being correctly identified is high. Conversely, if several matches have very similar SI matching factors, the probability of a correct assignment is low.

The Chemical Structure View appears when you perform a NIST library search. The Chemical Structure view displays the chemical structure, the compound name, and the chemical formula of the compound that you select in the NIST Search Result view.

Reviewing the Results of an mzVault Search

For information about starting an mzVault library search, see [Starting a Library Search](#).

Follow these procedures:

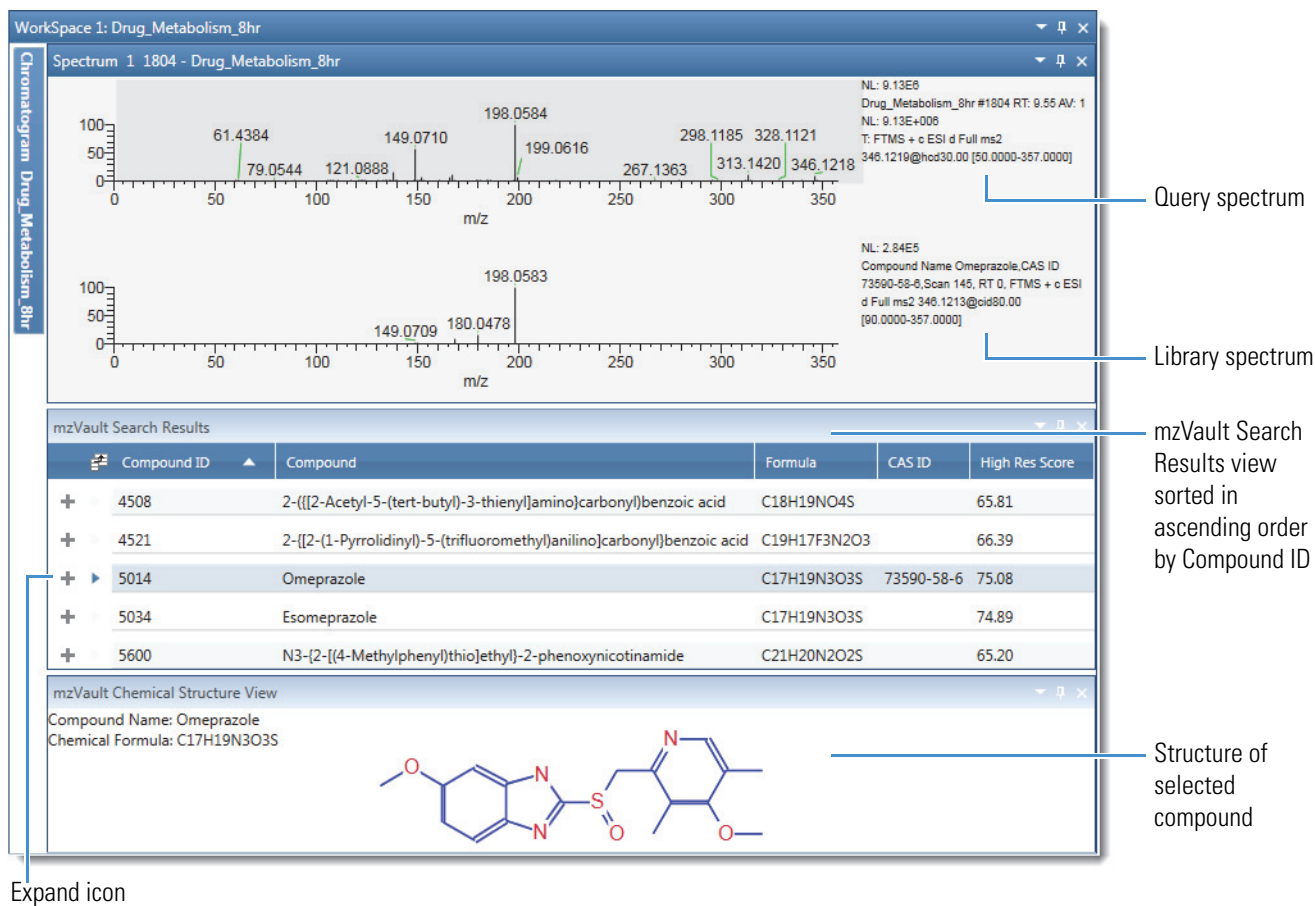
- To review the matching compounds found by an mzVault library search
 - To view all the matching scans for a selected compound
- ❖ **To review the matching compounds found by an mzVault library search**

In the mzVault Search Results view, select a compound.

The best matching spectrum for the library compound appears in the Spectrum view and its chemical structure appears in the mzVault Chemical Structure View.

Figure 140 shows the results of an mzVault library search. The query spectrum's plot is stacked above the library spectrum's plot. The mzVault Search Results view lists matching compounds in order of their compound ID numbers. The mzVault Chemical Structure view displays the two-dimensional structure of the compound with the highest score, from 0 to 100%.

Figure 140. mzVault library search results



- ❖ **To view all the matching scans for a selected compound**

1. Click the expand icon, **+**, to open the complete list of matching scans.

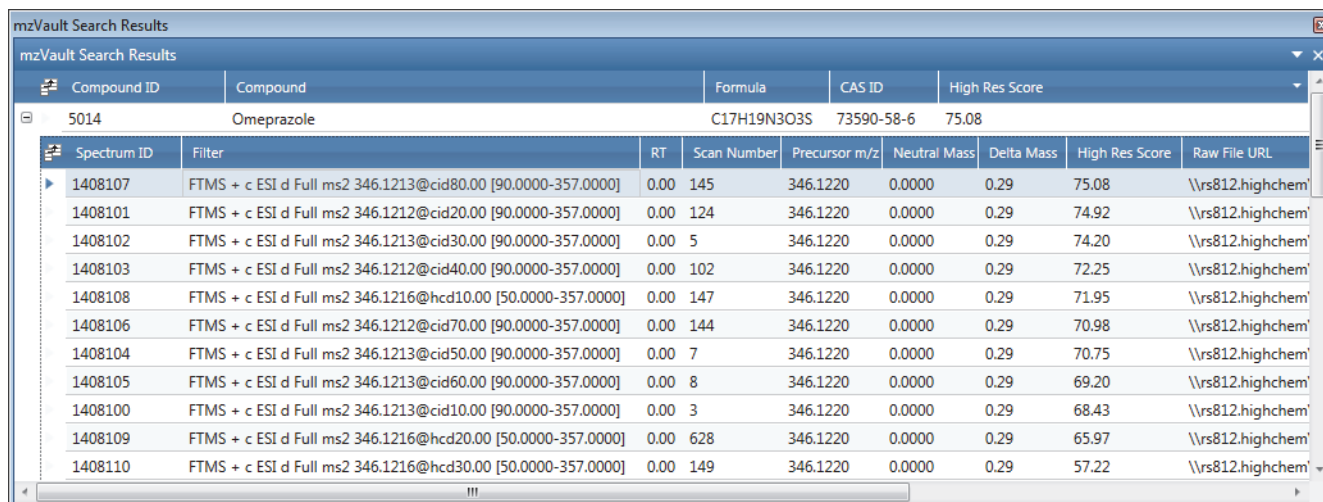
8 Searching Mass Spectrum Libraries

Reviewing the Results of a Local NIST or mzVault Library Search

- To view a scan in the Spectrum view, select it in the subtable.

Figure 141 shows the subtable of matching scans in the mzVault database file. For details about the table columns, see the next topic “mzVault Search Results View.” To modify and rerun the search, see [To reset the mzVault search parameters and query spectrum](#).

Figure 141. Subtable of matching scans in the mzVault Search Results view



Compound ID	Compound	Formula	CAS ID	High Res Score				
5014	Omeprazole	C17H19N3O3S	73590-58-6	75.08				
Spectrum ID	Filter	RT	Scan Number	Precursor m/z	Neutral Mass	Delta Mass	High Res Score	Raw File URL
1408107	FTMS + c ESI d Full ms2 346.1213@cid80.00 [90.0000-357.0000]	0.00	145	346.1220	0.0000	0.29	75.08	\\rs812.highchem
1408101	FTMS + c ESI d Full ms2 346.1212@cid20.00 [90.0000-357.0000]	0.00	124	346.1220	0.0000	0.29	74.92	\\rs812.highchem
1408102	FTMS + c ESI d Full ms2 346.1213@cid30.00 [90.0000-357.0000]	0.00	5	346.1220	0.0000	0.29	74.20	\\rs812.highchem
1408103	FTMS + c ESI d Full ms2 346.1212@cid40.00 [90.0000-357.0000]	0.00	102	346.1220	0.0000	0.29	72.25	\\rs812.highchem
1408108	FTMS + c ESI d Full ms2 346.1216@hcd10.00 [50.0000-357.0000]	0.00	147	346.1220	0.0000	0.29	71.95	\\rs812.highchem
1408106	FTMS + c ESI d Full ms2 346.1212@cid70.00 [90.0000-357.0000]	0.00	144	346.1220	0.0000	0.29	70.98	\\rs812.highchem
1408104	FTMS + c ESI d Full ms2 346.1213@cid50.00 [90.0000-357.0000]	0.00	7	346.1220	0.0000	0.29	70.75	\\rs812.highchem
1408105	FTMS + c ESI d Full ms2 346.1213@cid60.00 [90.0000-357.0000]	0.00	8	346.1220	0.0000	0.29	69.20	\\rs812.highchem
1408100	FTMS + c ESI d Full ms2 346.1213@cid10.00 [90.0000-357.0000]	0.00	3	346.1220	0.0000	0.29	68.43	\\rs812.highchem
1408109	FTMS + c ESI d Full ms2 346.1216@hcd20.00 [50.0000-357.0000]	0.00	628	346.1220	0.0000	0.29	65.97	\\rs812.highchem
1408110	FTMS + c ESI d Full ms2 346.1216@hcd30.00 [50.0000-357.0000]	0.00	149	346.1220	0.0000	0.29	57.22	\\rs812.highchem

mzVault Search Results View

The mzVault Search Results view appears in the workspace after you submit an mzVault search. The mzVault search returns a list of matching compounds.

❖ To display the mzVault Search Results view

Perform an mzVault library search (see [Performing a Local NIST or mzVault Library Search](#)).

The mzVault Search Results view consists of the top-level compound entries and the secondary-level spectrum entries for each compound hit.

Table 47 describes the columns in the compounds table in the mzVault Search Results view.

Table 47. Compounds table in the mzVault Search Results view (Sheet 1 of 2)

Column	Description
Compound ID	Displays the assigned entry number in the mzVault library. The mzVault application assigns entry numbers in sequential order, beginning with the number 1 for the first entry.
Compound	Displays the name of the compound associated with the matching spectrum in the library.
Formula	Displays the chemical formula for the library entry.

Table 47. Compounds table in the mzVault Search Results view (Sheet 2 of 2)

Column	Description
CAS ID	Displays the unique Chemical Abstracts Service™ registry number.
High Res Score	Displays a score calculated by a proprietary algorithm that indicates how well the library spectrum and the query spectrum match.

Table 48 describes the spectrum information for each compound entry.

Table 48. Spectrum information for each compound entry

Column	Description
Spectrum ID	Displays the identification number for the library spectrum.
Filter	Displays the scan filter from the raw data file.
RT	Displays the retention time from the raw data file.
Scan Number	Displays the scan number from the raw data file.
Precursor <i>m/z</i>	Displays the mass-to-charge ratio of the precursor ion from the scan filter for the library spectrum.
Neutral Mass	Displays the uncharged, neutral mass of the molecule.
Delta Mass	Displays the difference in mass between the mass-to-charge ratio of the precursor ion for the library entry and the mass-to-charge ratio of the precursor ion for the query spectrum, in parts per million.
High Res Score	Displays the score calculated by a proprietary algorithm that indicates how well the library spectrum and the query spectrum match.
Raw File URL	Displays the name and path of the raw data file where the matching spectrum in the library entry came from.

Modifying the Settings for a Local NIST or mzVault Library Search

After you run a NIST or an mzVault library search, you can modify the search settings on the respective page in the Info Bar, and then restart the search.

To modify and rerun a search, see the appropriate topic:

- [Modifying a NIST Search from the NIST Search Page](#)
- [Modifying an mzVault Search from the mzVault Search Page](#)

Modifying a NIST Search from the NIST Search Page

After running a NIST search, use the NIST Search page of the Info Bar to modify the search or the display in the NIST Search Results view.

Follow these procedures:

- To display the NIST Search page
- To modify the display in the NIST Search Results view
- To modify the NIST search

❖ To display the NIST Search page

Run a NIST library search (see [Performing a Local NIST or mzVault Library Search](#)).

When the search is complete, you can reset the search parameters.

Figure 142 shows the NIST Search page.

Figure 142. NIST Search page

Info Bar

Isotope Simulation | Peak Detection | MSn Browser | Detector Type | **NIST Search**

NIST Search

Search List

Library: mainlib, nist, msms, nist, msms2, nist, ri, replib

Hits: 10

Display: Stacked

Search Parameters

Search Type

Identity Similarity

Normal Simple

Quick Hybrid

MS/MS Neutral Loss

In-source HiRes MS/MS Hybrid

Options

Search With MW= 0

Reverse Search

Penalize Rare Compounds

Mass Defect

Enable

Defect (mmu) At Mass (amu)

0 1

Defect (mmu) At Mass (amu)

0 2

MS/MS and In-source HiRes Search Options

Search Tolerance

Precursor +- 10 ppm

Product ions +- 10 ppm

Make Default Apply Help

❖ **To modify the display in the NIST Search Results view**

In the Display list on the NIST Search page, make a different selection (see [Display](#) in [Table 49](#)).

❖ **To modify the NIST search**

Modify the parameter settings on the NIST Search page, and then click **Apply**.

[Table 49](#) describes the parameters for the NIST Search page.

Table 49. NIST Search page parameters (Sheet 1 of 5)

Parameter	Description
Search List	
Use the settings in this area to choose the library for library searches. Specify how many search results to display and how they are displayed.	
Library	Select the search libraries.
Hits	Select how many search results to display in the NIST Search Results View . Range: 1–100; default: 10
Display	Select how to display the library spectrum and measured spectrum, from these options: <ul style="list-style-type: none"> Stacked: Displays the submitted spectrum above the currently selected library spectrum in the match list. Exclusion: Subtracts the library spectrum from the submitted spectrum. Mirror: Plots the submitted spectrum that points up from the x axis and the library spectrum that points down from the x axis.
Search Type	
Use the options in this area to choose the type of library search to apply. The two main search types are Identity and Similarity. They differ primarily in the weightings of the spectrum as a function of mass.	
Identity	Applies an identity-search algorithm for the library matching of spectra.
Normal	(Default) Applies a normal identity-search algorithm for library matching of spectra. Use a normal identity search for low quality or unusual spectra. The search algorithm uses a standard prescreen search filter.

Table 49. NIST Search page parameters (Sheet 2 of 5)

Parameter	Description
Quick	<p>Applies a quick identity-search algorithm for library matching of spectra.</p> <p>Use this option if you are sure the spectrum or compound exists in the library. The search algorithm uses a fast prescreen search filter.</p>
MS/MS	Searches for an MS/MS spectrum in a library of MS/MS spectra.
In-source HiRes	<p>Searches for an in-source electron ionization (EI) spectrum in a library containing high-resolution in-source EI spectra—that is, a library where the EI-MS spectra are annotated with m/z values to several decimal places. Also searches for an MS/MS spectrum in a library of MS/MS spectra.</p> <p>Unlike an MS/MS search, this search does not compare precursor m/z values.</p>
Similarity	Applies a similarity-search algorithm for the library matching of spectra.
Simple	<p>Applies a simple similarity-search algorithm for library matching of spectra.</p> <p>This option finds a large set of spectra to compare with the submitted spectrum, and is generally slower than an identity search.</p> <p>Use a simple similarity search in these situations:</p> <ul style="list-style-type: none"> You know that the unknown spectrum is not in the library. The spectrum is of poor quality so that a reliable match is unlikely.
Hybrid	<p>Applies a hybrid similarity-search algorithm for library matching of spectra.</p> <p>This option uses a combination of the simple and neutral-loss search strategies. The neutral-loss search requires an estimate of the unknown's molecular weight. If the unknown compound contains chemical structures that generate both characteristic ions and neutral loss patterns, the search algorithm can identify these structures from the match list produced by this search.</p>

Table 49. NIST Search page parameters (Sheet 3 of 5)

Parameter	Description
Neutral Loss	<p>Applies a neutral-loss similarity-search algorithm for library matching of spectra.</p> <p>The neutral losses in a spectrum are the mass differences between the molecular ion and other major ions in the spectrum. Neutral-loss peaks can be very characteristic as spectral features for certain classes of compounds.</p> <p>In a neutral loss search, the algorithm examines the submitted spectrum and identifies the molecular ion. It submits the mass value of the molecular ion to the search algorithm along with the spectrum. The search algorithm calculates the significant neutral losses and compares them with the library data. It returns hits according to matches of the molecular ion and its neutral losses.</p>
MS/MS Hybrid	<p>Applies an MS/MS hybrid similarity-search algorithm for library matching of spectra.</p> <p>This option uses a combination of the simple and neutral-loss search strategies. In this search, the algorithm searches for an MS/MS spectrum in a library of MS/MS spectra.</p>
Options	
Search With MW=	Restricts the library search to entries with a particular molecular weight. Use the associated box to enter the molecular weight.
Reverse Search	Sorts matching library spectra by the reverse search match factor. By default, the algorithm sorts matches by the forward match factor.

Table 49. NIST Search page parameters (Sheet 4 of 5)

Parameter	Description
Penalize Rare Compounds	<p>Limits the impact of rare compounds by reducing the match factor.</p> <p>Use this option only when you select one or more of the NIST databases (such as MAINLIB). It has no effect on spectra in user libraries or other commercial libraries.</p> <p>Each reference spectrum in a NIST library contains a record of other commercial databases containing information about the compound. A compound is considered rare if it is present in a limited number of these databases.</p> <p>If you select the Penalize Rare Compounds option, the application reduces the match factors for matched compounds that are present in few or no databases other than the NIST libraries. The maximum reduction penalty is 50 out of 1000.</p> <p>Selecting this option leads to a relative increase in the match factors of common compounds, placing them higher in the match list than exotic isomers with nearly identical spectra.</p>
Mass Defect	
<p>Use this area to set the parameters for library searches to correct for the differences between the actual masses and the nominal integer masses of the atoms in a molecule. Assign a larger mass defect (in millimass units) for more massive molecules because, in general, they are composed of more atoms than less massive molecules. More massive molecules need a larger correction factor to approximate the linear function that the application uses to calculate masses.</p>	
Enable	Include mass defect values for library searches.
Defect (mmu)	<p>Specify the mass defect (in millimass units).</p> <p>Specify a smaller value for the lower mass ranges in the first box and a larger value for the higher mass ranges in the second box.</p>
At Mass (amu)	<p>Specify the masses at which the application applies the specified mass defect values to calculate mass.</p> <p>Specify a smaller mass value in the first box and a larger mass value in the second box.</p>

Table 49. NIST Search page parameters (Sheet 5 of 5)

Parameter	Description
MS/MS and In-source HiRes Search Options	
Use this area to set the parameters for library searches involving the comparison of peaks in the search spectrum with library spectra whose precursor m/z value or product mass spectral peak might need to fall within a specified tolerance setting.	
Precursor +-	Specify the range of the m/z tolerance for precursor ions, in either ppm or m/z units. Range: 0.015 to 100 000 ppm or m/z 6×10^{-5} to 500; default: 10 ppm
Product Ions +-	Specify the range of mass spectral peak tolerance for product ions, in either ppm or m/z units. Range: 0.015 to 100 000 ppm or m/z 6×10^{-5} to 500; default: 10 ppm
Button	
Make Default	Saves the current settings as the default settings.
Apply	Starts the library search with the entered settings.
Help	Opens the FreeStyle Help to the NIST library search topic.

Modifying an mzVault Search from the mzVault Search Page

Use the mzVault Search page in the Info Bar to search the mzVault library for library entries that match your selected fragmentation spectrum. You can filter the search by compound name, chemical formula, or precursor m/z value.

Follow these procedures:

- [To open the mzVault Search page](#)
- [To modify the display in the mzVault Search Results view](#)
- [To reset the mzVault search parameters and query spectrum](#)

❖ To open the mzVault Search page

Run an mzVault search (see [Performing a Local NIST or mzVault Library Search](#)).

When the search is complete, you can reset the search parameters.

[Figure 143](#) shows the mzVault Search page.

8 Searching Mass Spectrum Libraries

Modifying the Settings for a Local NIST or mzVault Library Search

Figure 143. mzVault Search page

Filter Type	Filter Criterion	Match/Custom

❖ **To modify the display in the mzVault Search Results view**

In the Display list on the mzVault Search page, make a different selection (see [Display](#) in [Table 50](#)).

❖ **To reset the mzVault search parameters and query spectrum**

1. If applicable, click the Browse icon next to the Library list and select another mzVault library.

Note The FreeStyle application lets you select the access-restricted mzVault library.

2. In the Hits box, type the maximum number of compounds that you want the search to return.

3. From the Display list, select how you want to display the matching spectra in the Spectrum view (see [Setting Up the Default Library Search Parameters](#)).
4. (Optional) In the Score Cutoff box in Search Parameters area, specify the minimum score for a match to be returned.

Default: 0

5. Select the type of search to conduct, either **Forward** or **Reverse**.
6. (Optional) For filtered searches, do the following:
 - a. Select the check box on the left to make the filter selections available. You can select up to six filters.
 - b. Select a filter type. You can select the filters from the 12 filter types from the dropdown menu.
 - c. Set up the filter criterion for the selected filter.

For details about selecting the type of filters, see [Table 50](#).

- d. From the Match/Custom dropdown list, select **Match** to update the filter criteria based on the selected scan, or **Custom** to retain the edited filter criteria as you navigate to different scans. The Custom option is applicable only for the following filter types:
 - Collision energy
 - Precursor m/z
 - Retention time
 - Retention time range
7. In the Fragment Tolerance area, select the **Automatic** or **Manual** option to indicate how to specify the fragment tolerance.
8. (Optional) Select whether you want the application to filter out low intensity peaks from the query spectrum before performing the search.
9. (Optional) To have the mzVault application ignore any mass that is within 2.2 Da of the precursor ion for the purpose of calculating scores, select the **Remove Precursor** check box.
10. Click **Apply**.

The application applies the search parameters and displays the mzVault search results for the selected criteria.

8 Searching Mass Spectrum Libraries

Modifying the Settings for a Local NIST or mzVault Library Search

Table 50 describes the parameters for the mzVault Search page in the Info Bar.

Table 50. mzVault Search page parameters (Sheet 1 of 4)

Parameter	Description
Search List	
Use the settings in this area to choose the library for library searches. Specify how many search results to display and how they are displayed.	
Library	Select the library (database file) for the search.
Hits	Specify how many search results to display in the mzVault Search Results View . Range: 1–10 000; default: 10
Display	Select how to display the library spectrum and query spectrum: <ul style="list-style-type: none">Stacked—Displays the query spectrum above the currently selected library spectrum in the match list.Mirror—Plots the query spectrum on the positive y axis from 0 to 100% intensity and the library spectrum on the negative y axis from 0 to –100% intensity.Exclusion—Subtracts the library spectrum from the query spectrum.
Search Parameters	
Score Cut-Off	Filters out results below a percentage of the highest score. For example, if the highest score for a search result is 0.900 and the setting in this box is 60, the search does not return results with scores below 0.540 ($0.9 \times 60\% = 0.540$). Range: 0.00 to 100.00; default: 0.00
Search Type	
Forward	Conducts a forward search, which bases the search on a comparison of the peaks in the unknown spectrum against the peaks in a library spectrum. If the unknown spectrum includes a peak that is not in a given library spectrum, the score for the match is negatively affected. Use a forward search when the unknown spectrum is of high quality—that is, when it has good fragmentation and few low-intensity background peaks.

Table 50. mzVault Search page parameters (Sheet 2 of 4)

Parameter	Description
Reverse	<p>Conducts a reverse search, which bases the search on a comparison of the peaks in a library spectrum against the peaks in the unknown spectrum.</p> <p>If the unknown spectrum does not contain a peak that is in the library spectrum, the score for the match is negatively affected, but the presence of additional peaks in the unknown spectrum has no effect on the score.</p> <p>Use a reverse search if the unknown spectrum includes peaks from several components or has a lot of background noise.</p>
Filters	
Collision Energy	<p>Filters the library entries with scan data acquired at the specified collision energy within the specified tolerance.</p> <p>Valid entries: Numeric value > 0</p>
Compound	<p>Filters the library entries by the name or part of the name of the compound.</p> <p>Valid entries: Alphanumeric and special characters</p>
Compound Class	<p>Filters the library entries for the selected compound class.</p> <p>The Compound Class list includes only compound classes found in the selected library.</p>
Curation Type	<p>Filters the library spectra for the selected curation type.</p> <p>The Curation Type list includes only the curation types in the selected library.</p>
Formula	<p>Filters the library entries by the chemical formula of the library entry.</p> <p>Valid entries: Alphanumeric and special characters</p>
Fragmentation Mode	<p>Filters the library spectra with the same fragmentation mode as the query spectrum.</p> <p>Default: Set to 'Match'</p>
Peptide Sequence	<p>Filters the library entries that include the specified peptide sequence.</p>

8 Searching Mass Spectrum Libraries

Modifying the Settings for a Local NIST or mzVault Library Search

Table 50. mzVault Search page parameters (Sheet 3 of 4)

Parameter	Description
Precursor mass range	Filters the library spectra for the specified range of precursor m/z values. Valid entries: The maximum value must be greater than the minimum value by 0.1amu
Precursor <i>m/z</i>	Filters the library entries by the mass-to-charge ratio of the precursor ion for the library fragmentation scans. When you select this option, the Precursor m/z and the Precursor Tolerance boxes become available. Valid entries: Numeric values up to four decimal places.
Retention Time	Filters the library spectra from the specified retention time within a tolerance. Default: Retention time of the query scan; valid entries: Positive numeric values
Retention time range	Filters the library spectra from the specified retention time range. Valid entries: The maximum value must be greater than the minimum value by 0.1 min.
Scan Filter	Filters the library spectra for the specified scan filter. Valid entries: All or part of a valid scan filter; alphanumeric characters
Filter Criterion	
Defines the filter criteria based on the Filter type.	
Match/Custom	
Match	Updates the filter criteria based on the selected scan.
Custom	Retains the edited filter criteria as you navigate to different scans.
Query Spectrum	
Use the settings in this area to define the mass tolerance and Intensity threshold settings.	
Fragment Tolerance	
Automatic	(Default) Uses an internal algorithm to calculate the mass tolerance.

Table 50. mzVault Search page parameters (Sheet 4 of 4)

Parameter	Description
Manual	<p>Selecting this option activates the mass tolerance box and the corresponding units list. Enter the mass tolerance for the fragment ions.</p> <p>Units: amu, mmu, or ppm</p> <p>Range: 0.001 to 1000.00; default: 10.000</p>
Intensity Threshold	
On or Off	<p>Determines whether the application filters out peaks with an intensity below the Intensity Threshold (%) setting.</p> <ul style="list-style-type: none"> • (Default) On—Filters out peaks that are less than the specified value. Enables the Intensity Threshold% parameters. • Off—Does not filter out peaks on the basis of their intensity. <p>The Intensity Threshold parameter calculates the score that indicates how well a spectrum matches a library spectrum and displays the score in the High Res Score column.</p>
Intensity Threshold%	
Automatic	(Default) Uses an internal algorithm to calculate the threshold value.
Manual	Lets you to specify the threshold value to use for filtering spectrum peaks. The application takes the default value from the scan that you selected from the raw data file.
Additional Query Spectrum parameter	
Remove Precursor Ion	<p>Determines whether the application removes peaks within 2.2 Da of the precursor ion. In an MS/MS experiment, the fragmentation spectrum does not normally contain a peak for the precursor ion, but sometimes it can appear and interfere with the scoring algorithm. This option corrects for the presence of a peak for the precursor ion.</p> <p>Default: Selected</p>
Button	
Make Default	Saves the current settings as the default settings.
Apply	Starts the library search with the entered settings.
Help	Opens the FreeStyle Help to the mzVault search topic.

Searching the Online mzCloud Mass Spectral Database

The mzCloud mass spectral database includes thousands of fragmentation spectra.

❖ **To perform an mzCloud library search for a matching fragmentation spectrum**

1. In a Spectrum view, select a fragmentation spectrum.
2. In the Workspace Processing toolbar, click **mzCloud Search**.

Figure 144 shows the mzCloud Spectrum Search dialog box that opens in your default browser when you run an mzCloud search.

Figure 144. mzCloud search parameters in a browser

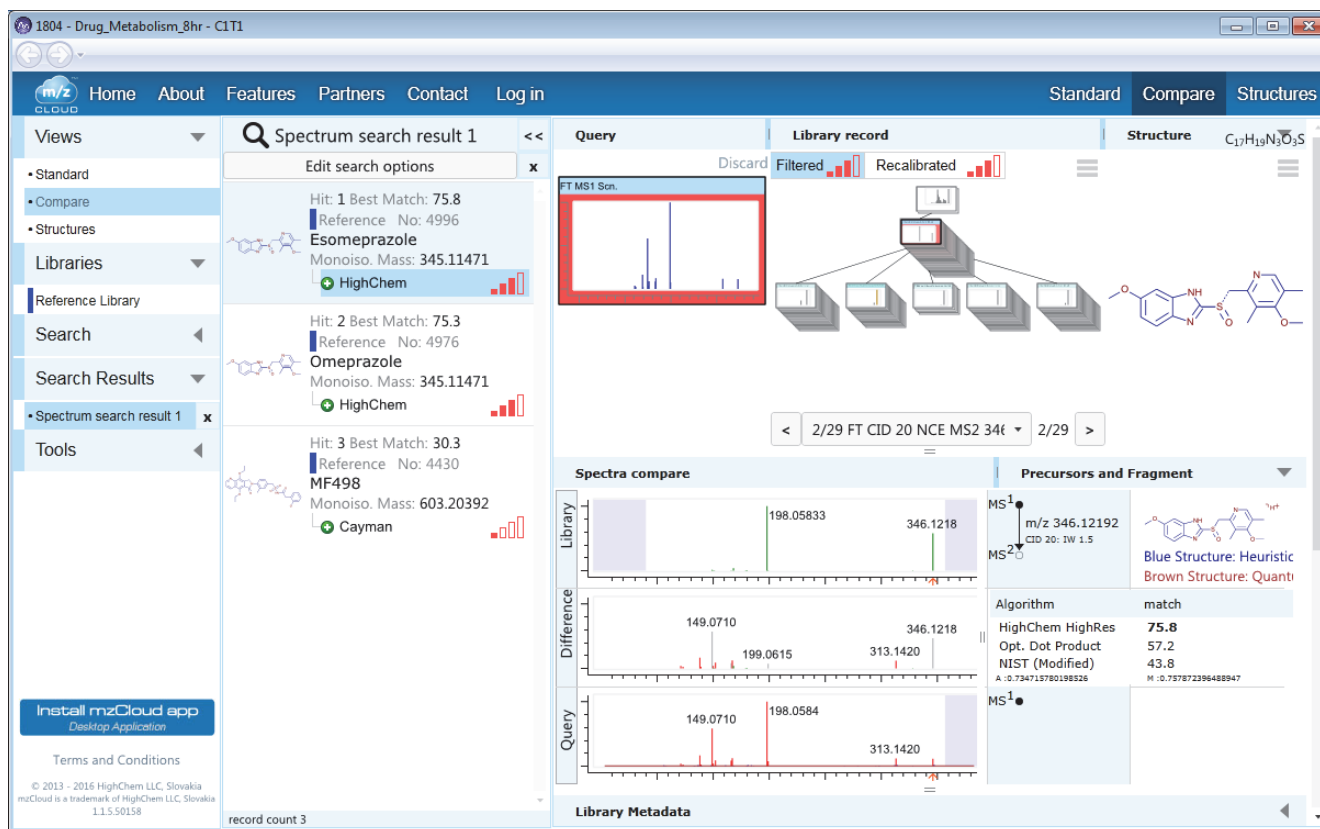
The screenshot displays the 'Spectrum Search' dialog box with the following settings:

- Ionization Mode:** Positive (selected), Negative
- Search Type:** Substructure Identification (Search in MSⁿ) (selected), Compound Identification (Search in MS²)
- Search in:** Filtered Spectra (checked), Recalibrated Spectra (checked), Average Spectra (unchecked), Composite Spectra (unchecked)
- Libraries:** Reference (checked)
- Algorithm:** HighChem HighRes (selected), Opt. Dot Product, NIST (Modified)
- Filters:** Spectrum, Compound
- Query Spectrum:** FTMS + c ESI d Full ms2 346.1219@hcd30.00 [50.0000-357.0000]. The plot shows peaks at m/z 138.0914, 149.0710, 198.0584, and 313.1420.
- Precursor m/z:** m/z 346.12190 ± 0.0040 (selected)
- Accuracy (half width):** From source (selected), Manual 0.001 Th
- Query Spectrum Noise Filter:** Intensity Threshold (%) 2.000 (checked), Num. of highest peaks 10 (unchecked)

Buttons: OK, Cancel

Figure 145 shows the search results for an mzCloud search.

Figure 145. mzCloud search results



Exporting a Mass Spectrum to the NIST MS Search Application

You can export a mass spectrum to the Library Search page of the NIST application.

Note Thermo Fisher Scientific provides the NIST application with the Xcalibur data system.

❖ To export a spectrum to the NIST application

1. Click the Spectrum view or spectrum plot of interest.
2. Click the **Workspace Processing** toolbar tab.
3. In the Library Search area of the toolbar, click **Export to NIST**.

The Library Search page of the NIST application opens with the search spectrum displayed.

Managing Libraries

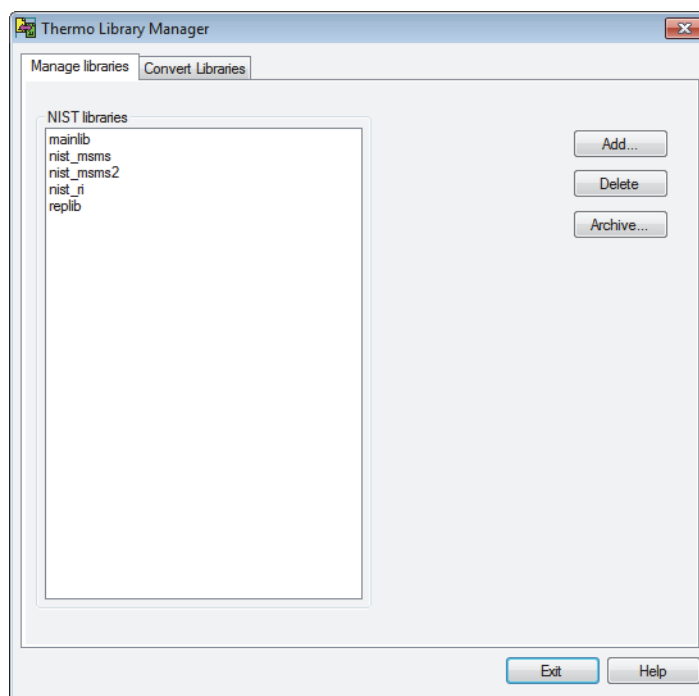
Use the Thermo Library Manager dialog box to add a library, copy a library, or delete a library.

When you add a library to the NIST libraries list, either copy the library file to the local computer or link to the library at a remote location without copying files.

❖ To add a library

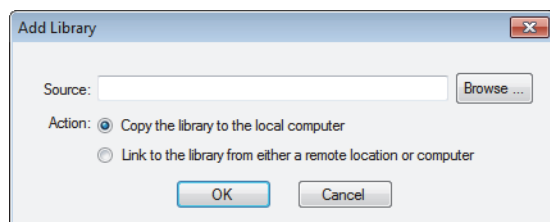
1. Click the **Workspace Processing** toolbar tab.
2. Click **Library Manager** to open the Thermo Library Manager dialog box (see [Figure 146](#)).

Figure 146. Thermo Library Manager dialog box



3. Click **Add** to open the Add Library dialog box (see [Figure 147](#)).

Figure 147. Add Library dialog box



4. Type the path for the new library file in the Source box, or click **Browse** to locate the file.

5. Select one of two options:
 - Copy the library to the local computer.
 - Link to the library from either a remote location or computer.

Tip Because libraries can be large, select the link option rather than the copy option to save time.

6. Click **OK**.

The data system adds the library to the NIST Libraries list in the Thermo Library Manager dialog box.

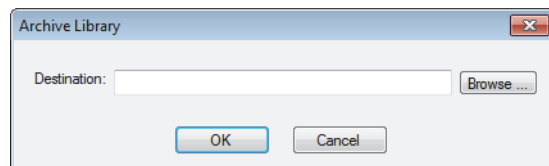
7. If you have no more tasks to complete with the Library Manager, click **Exit**.

❖ To copy a library for archiving

1. In the Thermo Library Manager dialog box (see [Figure 146](#)), select a library in the NIST Libraries list.
2. Click **Archive**.

The Archive Library dialog box opens (see [Figure 148](#)).

Figure 148. Archive Library dialog box



3. In the Archive Library dialog box, type the path for the copied library file in the Destination box, or click **Browse** to find the location.
4. To copy the selected library to the remote location, click **OK**.
5. If you have no more tasks to complete with the Library Manager, click **Exit**.

You can copy a selected library to another directory on the computer or network by using the Archive feature.

❖ To delete a selected library in the NIST Libraries list

1. In the Thermo Library Manager dialog box (see [Figure 146](#)), select a library to delete from the NIST Libraries list.
2. Click **Delete**.

The system prompts you to confirm the deletion.

3. Click **Yes**.

8 Searching Mass Spectrum Libraries

Spectrum Workspace Processing Toolbar – Library Search Buttons

4. If you have no more tasks to complete with the Library Manager, click **Exit**.

IMPORTANT After you delete a library, you cannot bring it back. Make sure that you want to delete a library permanently before you proceed, especially a shared library on the network.

Spectrum Workspace Processing Toolbar – Library Search Buttons

Use the buttons in the Library Search area of the Workspace Processing toolbar to perform library searches.

❖ To display the Spectrum Workspace Processing toolbar

1. Click a [Spectrum View](#) or a [MultiSpectrum View](#) to select it.
2. Click the **Workspace Processing** toolbar tab.

[Figure 149](#) shows the Library Search area in the Workspace Processing toolbar, and [Table 51](#) describes the toolbar buttons.

Figure 149. Library Search areas of the Workspace Processing toolbar

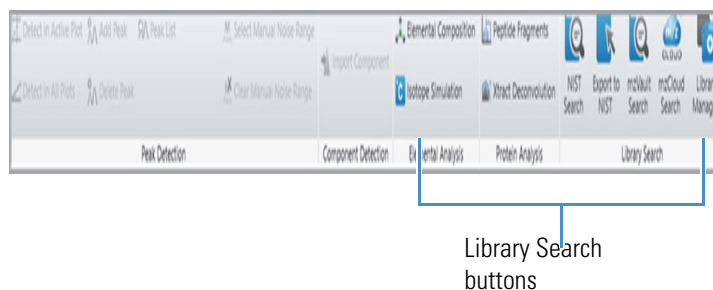


Table 51. Spectra-specific Workspace Processing buttons (Sheet 1 of 2)

Button	Description
Library Search	
NIST Search	Opens the Library Search page in the Info Bar, where you run a NIST library search on the active spectrum and displays the results in the NIST Search Results view. You can modify the library search parameters on the Modifying a NIST Search from the NIST Search Page of the Info Bar. See Performing a Local NIST or mzVault Library Search .
Export to NIST	Exports a spectrum to the NIST application for a search. The NIST application opens and displays the search results.
mzVault Search	Opens the mzVault Search page in the Info Bar, runs an mzVault library search on the active spectrum, and displays the results in the mzVault Search Results view.

Table 51. Spectra-specific Workspace Processing buttons (Sheet 2 of 2)

Button	Description
mzCloud	Uploads a spectrum to mzCloud.org for a search. The website opens and displays the search parameters (see Figure 144). Select the appropriate settings and click OK . The website displays the results (see Figure 145).
Library Manager	Opens the Thermo Library Manager dialog box, where you select or create libraries of spectra. See Managing Libraries .

Simulating Isotope Distributions

The FreeStyle application includes a utility that predicts and displays the isotopic distribution for a specified chemical formula or peptide sequence. The simulated spectrum is based on the elemental composition of the compound or peptide and the charge state distribution of its adduct ions.

To display simulated isotope distribution plots for chemical formulas and peptide sequences, see these topics.

Contents

- [Simulating the Isotope Distribution for a Chemical Formula or Peptide](#)
- [Displaying Centroid and Profile Spectra in the Same Window](#)
- [Isotope Simulation Page](#)

Simulating the Isotope Distribution for a Chemical Formula or Peptide

A simulated isotopic distribution is based on a chemical formula or peptide sequence. For details about the parameter settings, see [Isotope Simulation Page](#).

When you open the FreeStyle application, the Isotope Simulation page opens in the Info Bar. Without opening a raw data file, you can use the Isotope Simulation page to display, in a separate window, the isotope pattern at a specified resolution for compounds or peptides. After opening a raw data file (see [To open a raw data file or a sequence file](#)), you can also use the Isotope Simulation page to insert a simulated isotope pattern or profile spectrum in a Spectrum view and an EIC trace in the Chromatogram view of a Workspace.

❖ To display isotope simulations

1. If the Isotope Simulation page is not visible, click the **Isotope Simulation** tab.
2. For the first simulation, do one of the following:
 - To display the simulation in the Isotope Simulation window, keep the **New** option.

–or–

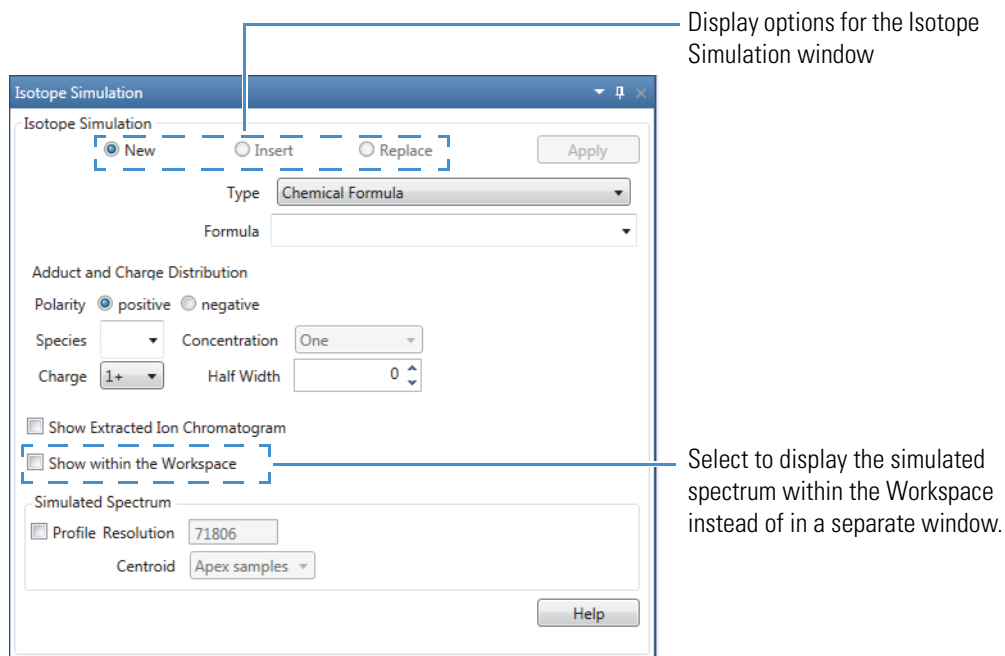
9 Simulating Isotope Distributions

Simulating the Isotope Distribution for a Chemical Formula or Peptide

- To display the simulation in the Spectrum view of a Workspace, open a raw data file, and then on the Isotope Simulation page, select the **Show within the Workspace** check box.

Figure 150 shows the default settings on the Isotope Simulation page.

Figure 150. Isotope Simulation page without an open workspace



3. Enter the formula for a compound or the sequence of a peptide:

- For a compound, select **Chemical Formula** and enter the chemical formula using the chemical symbols in the periodic table.

Type:
Formula: — Elemental composition of caffeine

- To specify a specific isotope, use square brackets before the element name to enclose the isotope mass number, for example, **[13]C**.

Type:
Formula: — Elemental composition of caffeine where one carbon-12 atom has been replaced with carbon-13

Type:
Formula: — Elemental composition of caffeine where all the carbon-12 atoms have been replaced with carbon-13

- To specify repeating moieties, such as those found in polymers, use parentheses.

Type

Formula Polyethylene glycol

- For a peptide, select **Peptide** and enter the peptide sequence using uppercase one-letter symbols for the amino acid residues (see [One- and Three-Letter Abbreviations for Amino Acid Residues](#)).

Type

Formula Kentsin, a tetrapeptide with the following amino acid sequence: Threonine-Proline-Arginine-Lysine

Note If you enter an incorrect formula, the application outlines the Formula box in red. Examples of incorrect entries include the following:

- Peptide—Lowercase letters or three-letter abbreviations
- Chemical formula—Chemical symbols that are not in the periodic table

- Under Adduct and Charge Distribution, do the following:
 - Select the polarity of the ion.
 - Select the adduct species, the charge of the most abundant ion, the concentration of the adduct species, and the half width of the charge distribution.

For example, to simulate a spectrum for both the protonated and sodiated adduct ions of a compound in a high-concentration sodium solution, select the following:

- Polarity: positive
- Species: Na
- Concentration: High
- Half-Width: 0

- To display an EIC trace for the specified ion in the Chromatogram view of a Workspace, open a raw data file if you have not already done so, and then select the **Show Extracted Ion Chromatogram** check box.
- In the Simulated Spectrum area, do one of the following:
 - To display a pattern, leave the **Profile** check box clear.

—or—

 - To display a simulated profile spectrum, select the **Profile** check box.
 - (Optional) Specify the resolution.

9 Simulating Isotope Distributions

Simulating the Isotope Distribution for a Chemical Formula or Peptide

Result:

- If you are working with raw data, the application automatically populates the Resolution box with the experimental resolution in the current raw data file.
- If you have not opened a workspace, the application populates the Resolution box with the default value of 1 000 000. You can type a value from 1 to 2 000 000.

Tip You can set up a simulated profile spectrum to determine whether your instrument is capable of resolving the theoretical isotope pattern for a specific ion, or you can set up several simulations to determine the required resolution.

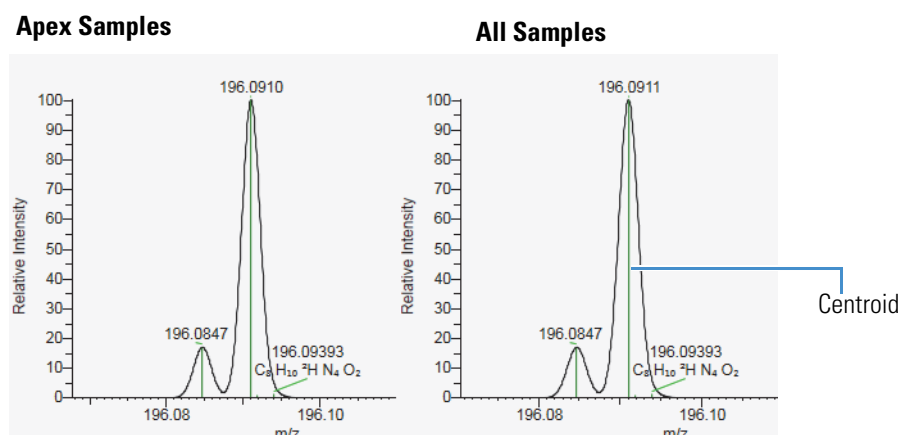
c. Select the centroiding algorithm.

Algorithm	Calculation
Apex Samples	Uses the sampling points nearest the peak apex.
All Samples	Uses all the sampling points in the profile peak, down to the baseline or valley between the resolved peaks.

Figure 151 shows a portion of a simulated profile spectrum. The two graphs show the difference between the two centroiding methods, where the left graph displays the results of the Apex Samples algorithm and the right graph displays the results of the All Samples algorithm. When you zoom in on the peak apices, you can see the slight difference between the centroiding methods.

Note When you display a simulated profile spectrum in the Spectrum view of a Workspace, the application displays the centroids for the Gaussian-shaped peaks as vertical green lines.

Figure 151. A1 peaks in a simulated profile spectrum (in the Spectrum view)

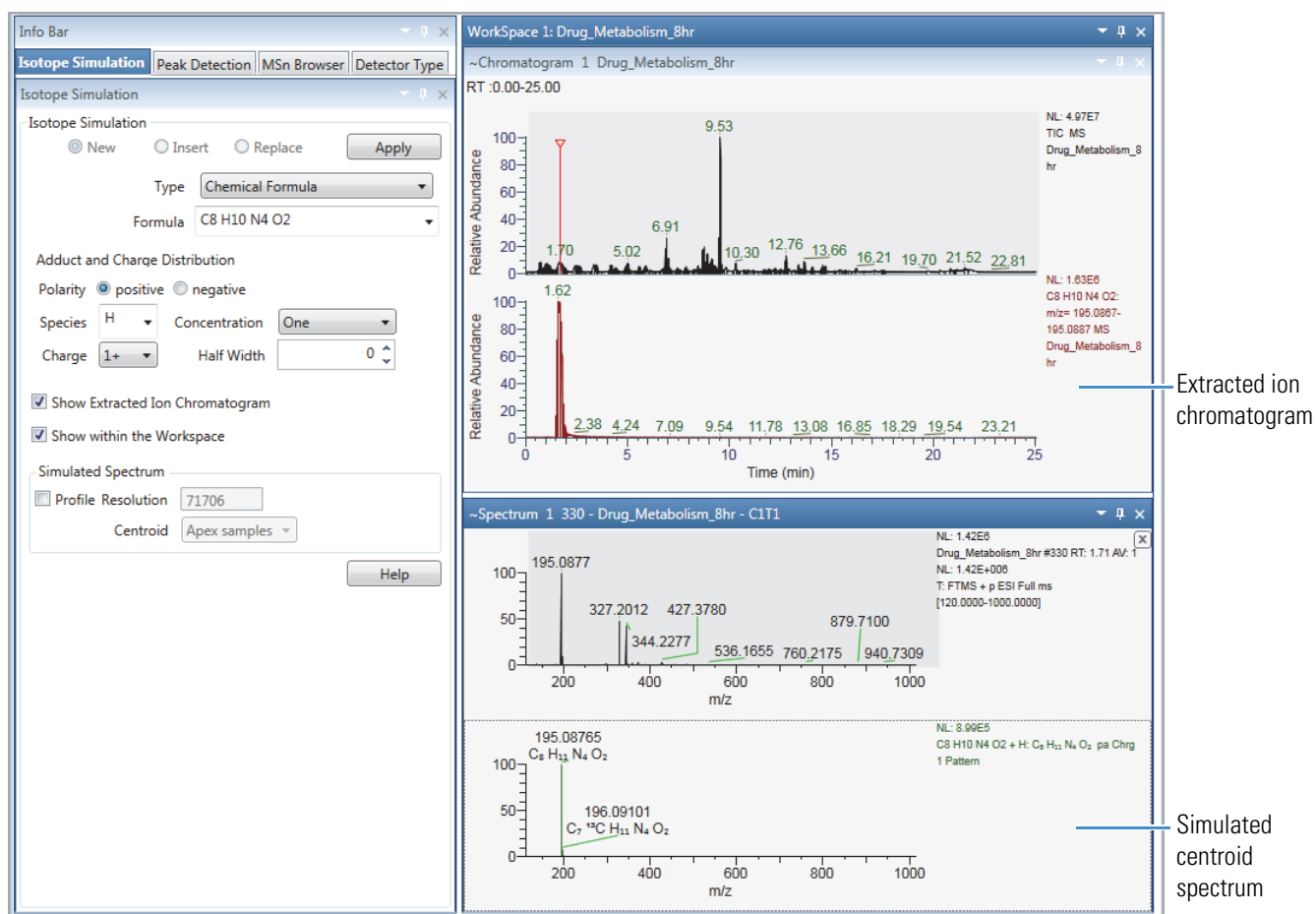


If you leave the Show within the Workspace check box clear, the Insert and Replace options become available after you run the first simulation and the Isotope Simulation window opens.

7. Click **Apply**.
8. For additional simulations, do the following:
 - a. Do one of the following:
 - To display additional simulations in separate Isotope Simulation windows, select **New**.
 - To insert additional simulations into one Isotope Simulation window, select **Insert**.
 - or–
 - To replace the current simulation with another simulation, select **Replace**.
 - b. Modify the compound or peptide information, and then click **Apply**.

Figure 152 shows the result of an isotope simulation for the protonated ion of caffeine.

Figure 152. Simulated centroid spectrum and EIC trace shown within the Workspace



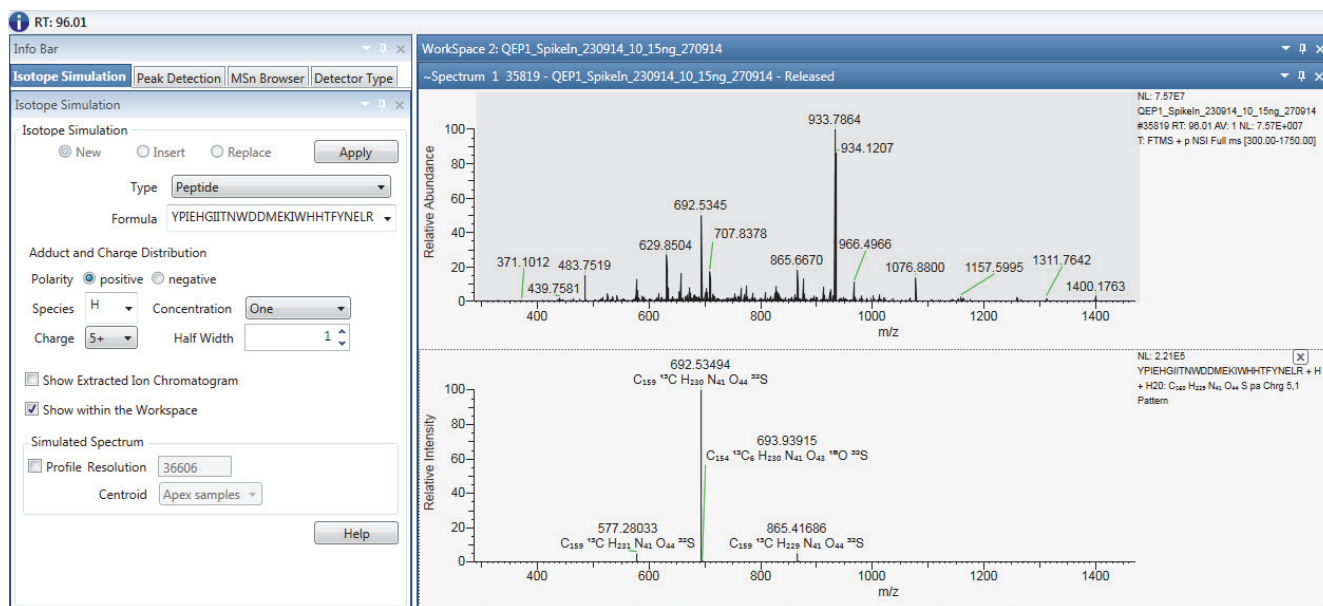
9 Simulating Isotope Distributions

Simulating the Isotope Distribution for a Chemical Formula or Peptide

Note Most peptides have multiple charge sites, which means that their spectra will show multiple charge states.

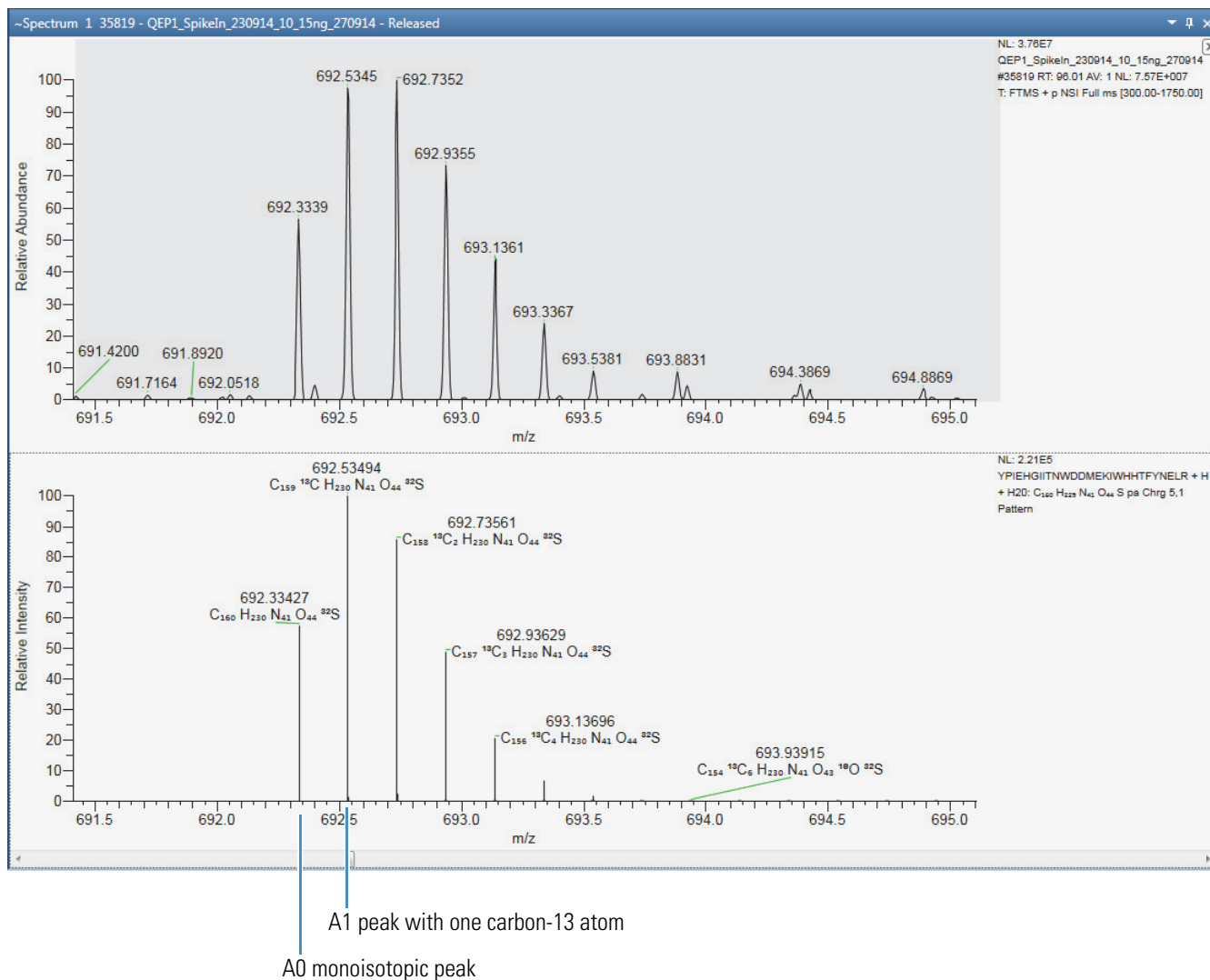
Figure 153 shows an isotope simulation for a peptide with a charge distribution pattern of 5,1—that is, a charge of +5 and a half width of 1. Notice that the monoisotopic peak (A0) is not the most intense peak. In compounds with large numbers of carbon atoms, the probability that one of the atoms will be a carbon-13 atom increases.

Figure 153. Example isotope simulation for a peptide



As shown in Figure 153 and Figure 154, when the number of carbon atom exceeds 91, the probability that the ion includes one carbon-13 atom exceeds 100% and the M+1 peak for the (A1) ion with one carbon-13 atom becomes the most intense peak.

Figure 154. m/z 692 isotope envelope



Displaying Centroid and Profile Spectra in the Same Window

The application does not display formula labels for simulated profile spectra. To view the formulas for the simulated peaks, run the simulation in both the profile and centroid modes within the same Isotope Simulation window.

❖ To view both simulation types in the Isotope Simulation window

1. Set up the initial simulation, make sure that the Show Within the Workspace and Profile check boxes are clear, and click **Apply**.
2. Select the **Insert** option.
3. Set up a second simulation for the same formula and charge distribution.

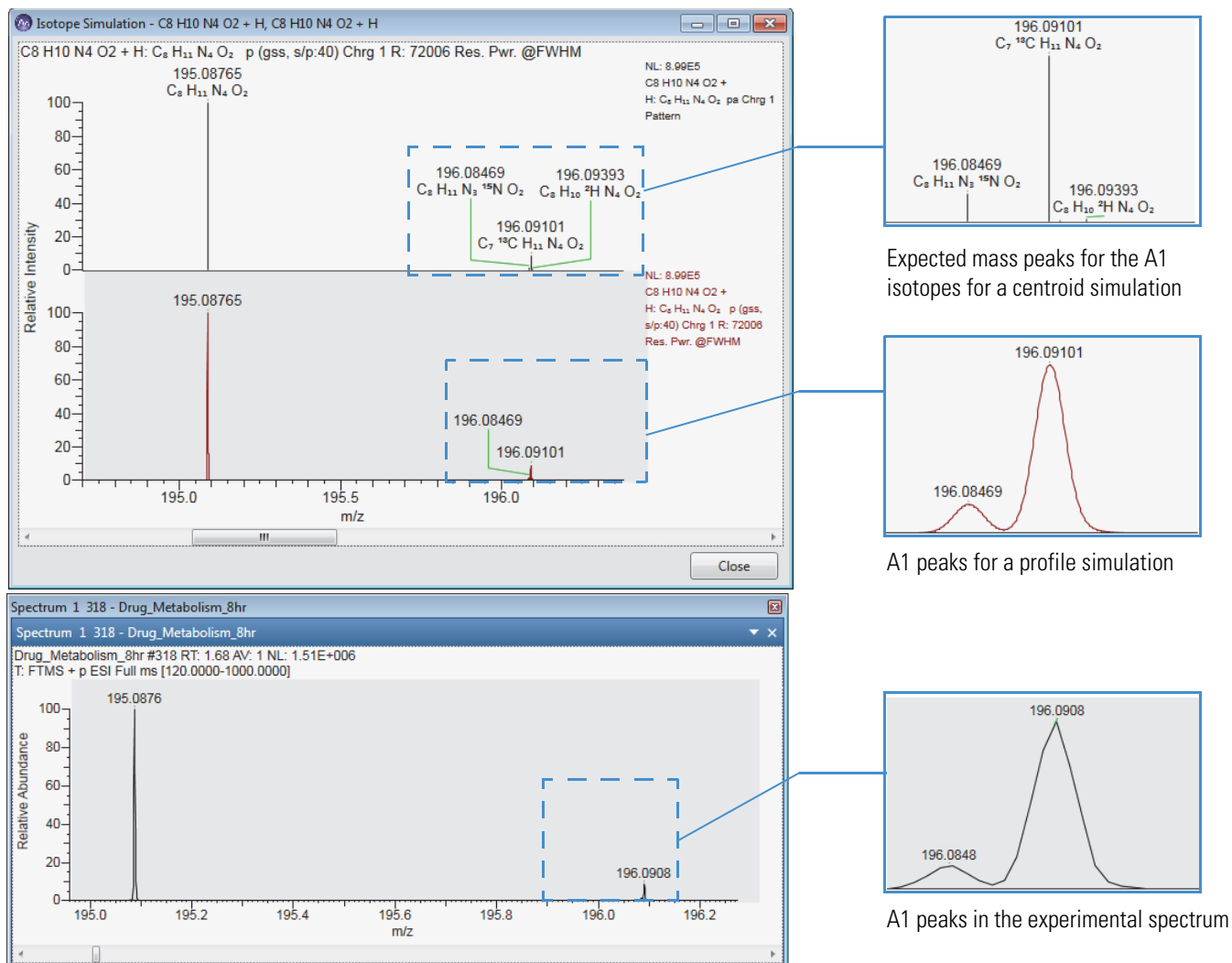
9 Simulating Isotope Distributions

Displaying Centroid and Profile Spectra in the Same Window

4. Select the **Profile** check box, and then select the centroiding algorithm.
5. Click **Apply**.

Figure 155 shows the isotope simulation and the mass spectrum for the protonated caffeine ion. At the experimental resolution, the isotope simulation shows that the A1 peaks at m/z 196.1 will not be completely resolved. The experimental spectrum shows a similar profile at m/z 196.1.

Figure 155. Simulation for caffeine versus the experimental spectrum for a caffeine sample



Isotope Simulation Page

Use the Isotope Simulation page of the Info Bar to create a simulated isotopic distribution plot. The Isotope Simulation page is always available. Without opening a workspace, you can only display a simulated spectrum in the stand-alone Isotope Simulation window. After you open a workspace, you can display an EIC trace in the Chromatogram view and add an isotope simulation plot to the Spectrum view.

For information about running the isotope simulation utility and reviewing the resulting spectrum, see [Simulating the Isotope Distribution for a Chemical Formula or Peptide](#).

Table 52 describes the parameters for the Isotope Simulation page.

Table 52. Isotope Simulation page parameters (Sheet 1 of 4)

Parameter	Description
New	Opens a separate Isotope Simulation window each time you click Apply.
Insert	Adds repeated simulations to the same Isotope Simulation window. Available when an Isotope Simulation window is open.
Replace	Replaces the current simulated spectrum with the latest simulation. Available when an Isotope Simulation window is open.
Type	Specify that the formula is either a Chemical Formula or a Peptide.
Formula	Enter the formula or peptide sequence for the compound or peptide that you want to create a simulated spectrum for. After you enter a formula or peptide sequence, it appears in a selection list for this box.
• Chemical Formula	Uses the custom periodic table elements for isotope simulation, if the custom periodic table is selected for elemental composition. For entering a chemical formula, do the following: <ul style="list-style-type: none">• Use the IUPAC symbol for each element—that is, match the notation, including the capitalization, in the periodic table of elements.• To enter a specific isotope, use square brackets before the element name to enclose the isotope mass number, for example, [13]C.• Use parentheses to specify repeating moieties such as those found in polymers, for example, HO(C2H4O)5H.

Table 52. Isotope Simulation page parameters (Sheet 2 of 4)

Parameter	Description
<ul style="list-style-type: none"> Peptide 	<p>For a peptide sequence, do the following:</p> <ul style="list-style-type: none"> Use the single capital letter abbreviations for amino acids. You can specify mixtures of substances by using additional symbols “+” (addition) and “x” (multiplication). Both symbols are required to specify a mixture. A valid mixture has the format substance x quantity + substance x quantity, for example, A x 2 + C x 5. <p>See One- and Three-Letter Abbreviations for Amino Acid Residues. The maximum molecular weight for the formula is less than 600 000 amu.</p>
Adduct and Charge Distribution	
Polarity	<p>Select the Positive or Negative option.</p> <p>Default: Positive</p>
Species	<p>Select, if necessary, the adduct for the simulated spectrum.</p> <ul style="list-style-type: none"> For positive polarity, you can select H, K, Na, NH₄, or no adduct. Or, you can type the chemical symbol or formula for a positive adduct such as Ca for calcium. For negative polarity, you can select H, Cl, OH, HCOO, or no adduct. Or, you can type the chemical symbol or formula for a negative adduct.
Charge	<p>Specify the charge of the adduct ion or peptide.</p> <ul style="list-style-type: none"> (Positive polarity) Range: 1–25; default: 1 (Negative polarity) Range: –1 to –25; default: –1
Half Width	<p>Simulate the number of additional charges on each side of the most abundant ion. Select a value from 0 to 99.</p> <p>For example: If you simulate charge 5 and a half width of 2, then the data system draws charges 5 ± 2, giving 7, 6, 5, 4, 3 (with the largest mass peak (normalized to 100% relative intensity) at charge 5).</p> <p>For most small molecules, simulate charge 1 with a half width of 0.</p> <p>To simulate an intensity distribution, the peaks at the edge of the distribution are shown at 5% of the height of the most abundant peak.</p>

Table 52. Isotope Simulation page parameters (Sheet 3 of 4)

Parameter	Description
Concentration	There are four possibilities for how the adduct is added to the ion: One, Low, High, and 100%.
<ul style="list-style-type: none"> One adduct 	<p>For positive charge states</p> <p>The application creates the specified charge state by adding the specified species for the first charge and H⁺ ions for subsequent charges. For example, a charge distribution of Charge 2+ (most abundant) and Half Width = 1 (2,1) and a Na⁺ adduct shows these ions and the simulated spectrum includes three isotope clusters:</p> <ul style="list-style-type: none"> – Charge +1) <i>Chemical_Formula</i>Na⁺ – Charge +2) <i>Chemical_Formula</i>HNa⁺² – Charge +3) <i>Chemical_Formula</i>H₂Na⁺³ <p>For negative charge states</p> <p>When you select H as the species, each charge state shows a loss of H⁺ to attain the specified charge—that is, a loss of one hydrogen ion for a charge of negative one, a loss of two hydrogen ions for a charge of negative two, and so on. If the molecule does not contain hydrogen, nothing is removed.</p> <p>For the neutral compound C₈H₈:</p> <ul style="list-style-type: none"> – Charge -1) C₈H₇ – Charge -2) C₈H₆ – Charge -3) C₈H₅ <p>For negative species, such as Cl⁻, the -1 charge state shows the addition of the negative species to attain the specified charge. Higher charge states show the addition of one Cl⁻ and the loss of one or more H⁺ to attain the specified charge.</p> <p>For the neutral compound C₈H₈:</p> <ul style="list-style-type: none"> – Charge -1) C₈H₈Cl – Charge -2) C₈H₇Cl – Charge -3) C₈H₆Cl
<ul style="list-style-type: none"> Low 	The ion with no adduct is included at 100% intensity, one adduct at 25% intensity, two adducts at 11% intensity, and so on (to the limit of the charge distribution).
<ul style="list-style-type: none"> High 	The ion with N adducts are included at 100% intensity, N-1 adduct at 25% intensity, N-2 adducts at 11% intensity, and so on, where N is the absolute value of the maximum charge simulated (to the limit of the charge distribution).

Table 52. Isotope Simulation page parameters (Sheet 4 of 4)

Parameter	Description
<ul style="list-style-type: none"> • 100% 	The ion of charge N contains M adducts (where M is the absolute value of N).
Show Extracted Ion Chromatogram	Select to display an EIC trace for the monoisotopic ion in the Chromatogram view.
Show within the Workspace	Select to display the isotope simulation in the Spectrum view of a Workspace.
Simulated Spectrum	
Profile and Resolution	<p>When the Profile check box is clear, the application generates a theoretical isotope pattern with infinite resolution—that is, it displays the peaks as one-dimensional sticks rather than as Gaussian peaks similar to the intensity versus frequency data that the instrument generates.</p> <p>When the Profile check box is selected, for each peak in the simulated pattern, the application uses a Gaussian peak-shape model that bases the peak width on the specified resolution. The application then sums the simulated Gaussian peaks to create a simulated profile spectrum for the specified resolution. The application displays the centroid for each profile peak as a green vertical line.</p> <p>When you open a raw data file in the workspace, the Resolution box displays the resolution value for the most intense peak (full width at half maximum [FWHM]) in the selected scan. You can overwrite this value.</p> <p>Tip To view the resolution values for each mass spectrum peak, click Peak Resolution in the Spectrum – Display Options toolbar.</p>
Centroid	<ul style="list-style-type: none"> • (Default) Apex Samples—Selects the sampling points nearest the peak apex to calculate the centroid of each simulated profile peak. • All Samples—Uses all the sampling points across the profile peak, down to the baseline or valley between resolved peaks.
Button	
Apply	Calculates the isotope distribution and displays the simulated spectrum by using the parameter settings.
Help	Opens the FreeStyle Help to the isotope simulation topic.

Working with CV Plots for FAIMS Data

The Thermo Scientific high-field asymmetric waveform ion mobility spectrometry (FAIMS) interface increases sensitivity by reducing chemical noise and matrix interferences. Use the CV Plot view to determine the optimal compensation voltages for your FAIMS-MS experiments.

Follow these topics.

Contents

- [CV Traces Dialog Box](#)
- [CV Plot View](#)

CV Traces Dialog Box

The CV Traces dialog box includes check boxes for the selected traces in the Chromatogram Ranges view.

❖ To select the traces for display in the CV Plot view

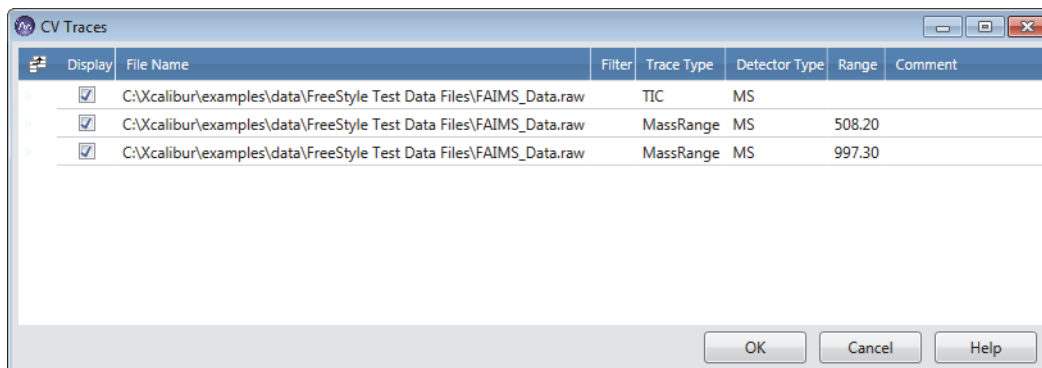
1. Open an Xcalibur RAW file with data acquired from a mass spectrometer with the FAIMS interface.
2. In the Chromatogram Ranges view, set up the traces of interest.
3. In the Workspace Options toolbar, click **CV Plot**.

The CV Traces dialog box opens ([Figure 156](#)).

4. Select the check boxes for the traces of interest.
5. Click **OK**.

The CV Plot view opens. For information about working with the CV Plot, see the next topic [CV Plot View](#).

Figure 156. CV Traces dialog box



CV Plot View

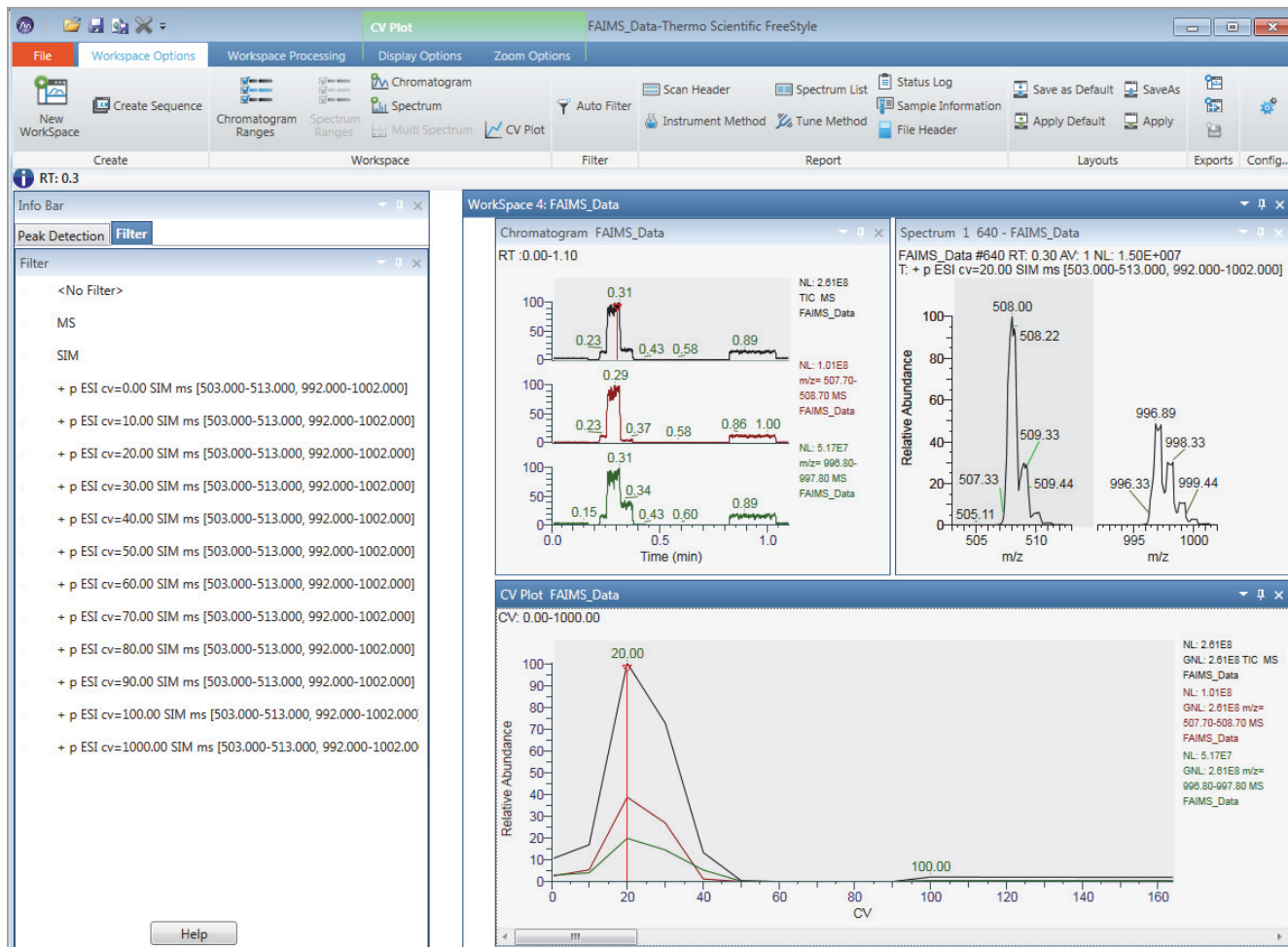
Use the CV Plot view to determine the optimum compensation voltages for your LC/MS/FAIMS experiments.

To open the CV Plot view, see [CV Traces Dialog Box](#). You can reopen the CV Traces dialog box to add or remove traces from this view.

[Figure 157](#) shows a CV Plot view with three overlaid traces: a TIC MS trace in black, a mass range m/z 507.2 ± 0.5 trace in red, and a mass range m/z 997.3 ± 0.5 trace in green.

The Filter page on the left shows scans acquired at 12 stepped compensation voltages from 0.00 to 1000 and two mass ranges. The CV Plot view, which plots the relative abundance versus the compensation voltage, shows a maximum response at a compensation voltage of 20 for both mass ranges.

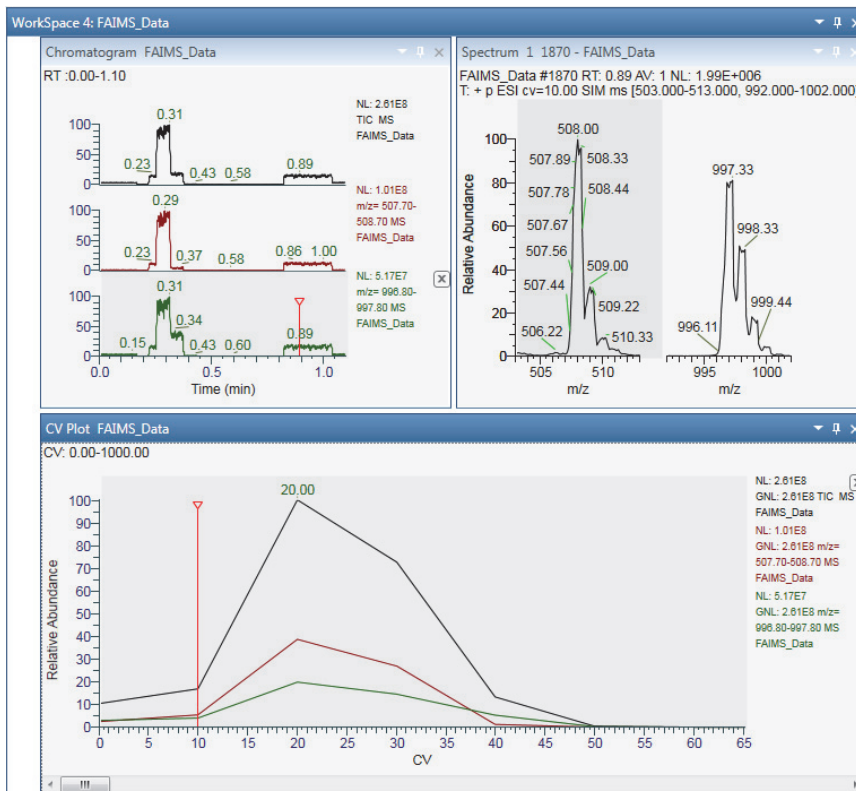
Figure 157. Workspace showing a selected CV Plot view



The CV Plot view, Chromatogram view, and Spectrum view are interactive—that is, as you move the red vertical bar across the CV voltage range on the x axis, the traces update in the Chromatogram and Spectrum views. Use this interactive behavior to determine the optimum compensation voltage for your analytes—that is, monitor the relative abundance of the ions in the Spectrum view as you move across the CV range in the CV Plot view.

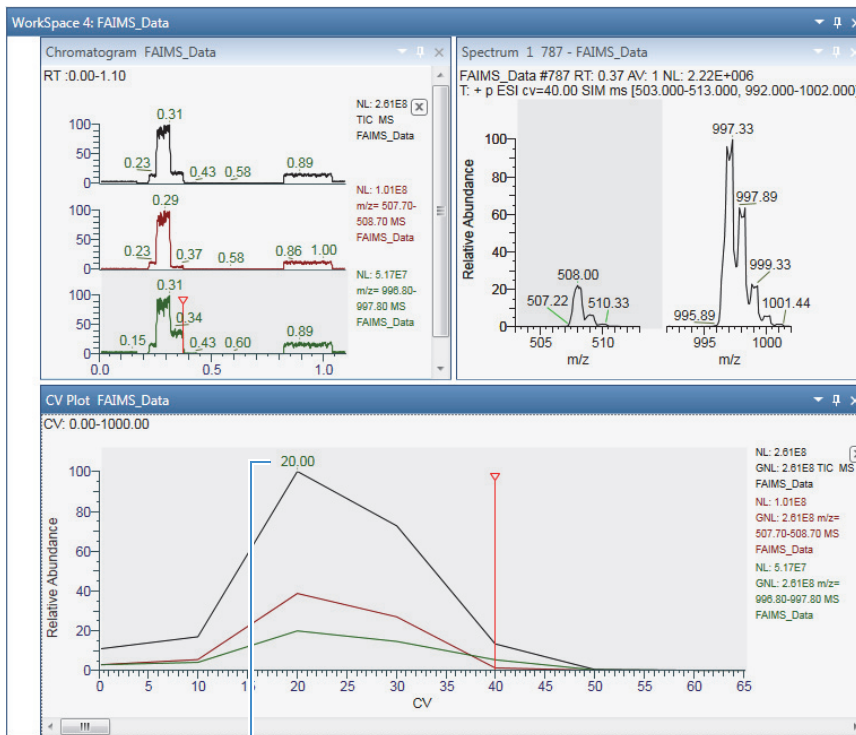
At a CV of 10, the ion at m/z 508.00 is relatively more abundant than the ion at m/z 997.33 (Figure 158).

Figure 158. Compensation voltage 10



At a CV of 40, the ion at m/z 997.33 is relatively more abundant than the ion at m/z 508.00 (Figure 159).

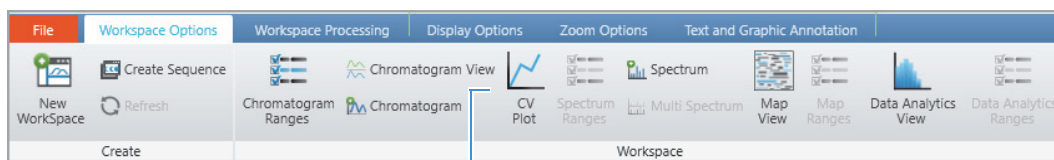
Figure 159. Compensation voltage 40



CV label

The display options for the CV Plot view are similar to those for the Chromatogram view (see “Chromatogram – Display Options Toolbar” on page 95). Use the CV Plot button in the Display Options toolbar to add the compensation voltage label to the y -axis maxima in the CV Plot view (Figure 160).

Figure 160. Display Options toolbar for the CV Plot view



CV Plot button

Peptide Fragment Annotation

The Thermo Scientific mass spectrometers support different strategies for acquiring peptide fragment spectra. The ETD or CID activation types performed in mass spectrometers provide different fragment scans of the same peptide.

This chapter explains how Peptide Fragment annotation in Freestyle application annotates the mass scans of the selected peptide sequence.

Contents

- [Overview of Peptide Fragments](#)
- [Fragmentation Methods](#)
- [Annotating Peptide Fragments](#)
- [Peptide Fragments Info Bar](#)
- [Peptide Results View](#)

Overview of Peptide Fragments

The fragment ions of peptides are produced by a CID or ETD activation process in which a peptide ion is fragmented in a collision cell. The fragment ion spectra contain peaks of the fragment ions formed by cleavage of the peptide bond and are used to determine the amino acid sequence. A fragment must have at least one charge for it to be detected. If this charge is retained on the N-terminal fragment, the ion is classed as a, b, or c. If the charge is retained on the C-terminal fragment, the ion type is x, y, or z. A subscript indicates the number of residues in the fragment.

The Peptide Fragment annotation displays all the fragmented ion series in the spectrum. The fragment ion series includes a&cx, c&z, and b&cy.

Fragmentation Methods

The Peptide Fragmentation in Freestyle application helps you to annotate the fragment ions for user defined peptide sequence in the spectrum. The spectrum header displays the following details: name of the raw file, activation type (CID or ETD) used in fragmentation, collision energy, and if the peptide is intact or fragmented. The peptide sequence is fragmented in one of the following fragmentation methods:

CID—Uses the collision-induced dissociation method of fragmentation, where molecular ions are accelerated to high kinetic energy in the vacuum of a mass spectrometer and then allowed to collide with neutral gas molecules such as helium, nitrogen, or argon. The collision breaks the bonds and fragments the molecular ions into smaller pieces. CID produces b and y ions.

ETD—Uses the electron transfer dissociation method of fragmentation, where singly charged reagent anions transfer an electron to multiply protonated peptides within an ion trap mass analyzer to induce fragmentation. ETD cleaves along the peptide backbone while side chains and modifications, such as phosphorylation, are left intact. This method is used to fragment peptides and proteins. ETD produces primarily c and z ions.

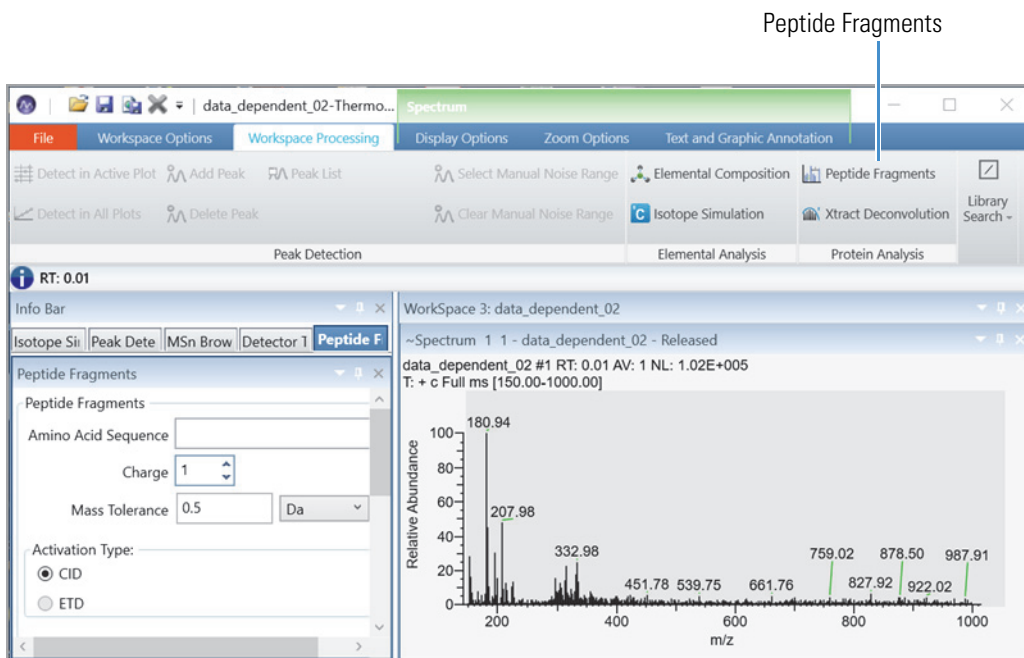
The fragment ions produced are identified according to where they are fragmented in the peptide. The a, b, and c fragment ions have a charge on the N-terminal side, and x, y, and z fragment ions have a charge on the C-terminal side. Fragment ions a*, b*, and y* are ions that have lost ammonia (–17 Da), and fragment ions a^o, b^o, and c^o are ions that have lost water (–18 Da). The subscript next to the letter indicates the number of residues in the fragment ion.

Annotating Peptide Fragments

The spectrum view shows the annotated spectra of the selected peptide sequence. When you use a CID or ETD fragmented peptide spectra, the view shows peaks of fragmented ions formed by the cleavage of peptide bonds. These fragments are used to confirm the amino-acid sequence.

Each alphabet represents an amino acid in the peptide sequence; for example, R for Arginine, H for histidine, G for Glycine, and so on. The carboxyl group of one amino acid links to the amine group of another amino acid to form an amino-acid sequence.

Figure 161. Workspace Processing – Peptide Fragments



❖ **To apply an amino acid sequence to a fragmented spectra**

1. Open a raw data file.
2. Click the Spectrum view to make it active.
3. From the ribbon menu, click **Workspace Processing**.
4. In the Protein Analysis area, click **Peptide Fragments**.

The Peptide Fragments page opens in the Info Bar.

5. Specify the parameter settings that you want to apply.

[Table 53](#) describes the parameters for the Peptide Fragments page.

6. Click **Apply**.

The application runs and annotates the selected fragment spectra.

Peptide Fragments Info Bar

Use the Peptide Fragments page in the Info Bar to specify the parameter settings.

❖ **To display the Peptide Fragments page**

1. Select the spectrum view to make it active.
2. Click the **Workspace Processing** tab.

3. Click **Peptide Fragments**.

Figure 162. Peptide Fragments Info Bar

Table 53 describes the parameters for the Peptide Fragments page.

Table 53. Peptide Fragments Page

Parameter	Description
Peptide Fragments	
Amino Acid Sequence	Specifies the amino acid sequence that you want to use to annotate the spectrum.
Charge	Specifies the charge state of the precursor ion. You can either select or manually enter a value.
Mass Tolerance	Specifies the mass tolerance. The application annotates the spectrum with ions within this mass tolerance. You can use daltons (Da) or millimass units (mmu).
Activation Type	
CID	Collision Induced Dissociation. The application automatically selects the option based on the spectrum selected.

Table 53. Peptide Fragments Page

Parameter	Description
ETD	Electron Transfer Dissociation. The application automatically selects the option based on the spectrum selected. IMPORTANT By default, Freestyle selects CID as the activation type for all spectrum types other ETD (for example, HCD, PQD, and so on).
Display Options	
Ion Series	Specifies the types of ion series to annotate the spectra. The application automatically selects the b and y ion series for CID activation type and c and z ion series for ETD activation type.
Charge State	Displays the charge state of the precursor ion as stored in the scan header of the raw data file. You can change the charge state to filter the fragment ions.
Show peptide results table	Select or clear to show or hide the Peptide Fragments table.

Peptide Results View

The Peptide Results table shows the analyzed spectra of the selected peptide sequence. It displays the major fragment ion series in the spectrum. The fragment ion series includes a, b, c, x, y, and z ions, as well as water and ammonia losses from b and y fragment ions. You can use the Peptide Results table to identify the spectral peaks that matches the amino-acid sequence provided.

❖ To display or hide the Peptide Fragments Table

1. In the Peptide Fragments page of the Info bar, select the **Show Peptide Results Table** check box.

The Peptide Fragments table appears.

Figure 163. Peptide Fragments Table

Peptide Fragments - Spectrum 1								
	(NH3+) - A>	S>	<T	T>	<V	V>	<R	R - (COOH)
b (+1)	72.04	159.08		260.12		359.19		515.29
b*(+1)	55.02	142.05		243.1		342.17		498.27
b°(+1)	54.03	141.07		242.11		341.18		497.28
y (+1)	462.27	375.24		274.19		175.12		533.3
y*(+1)	445.24	358.21		257.16		158.09		516.28
y°(+1)	444.26	357.22		256.18		157.11		515.29

2. To hide the table, clear the check box.

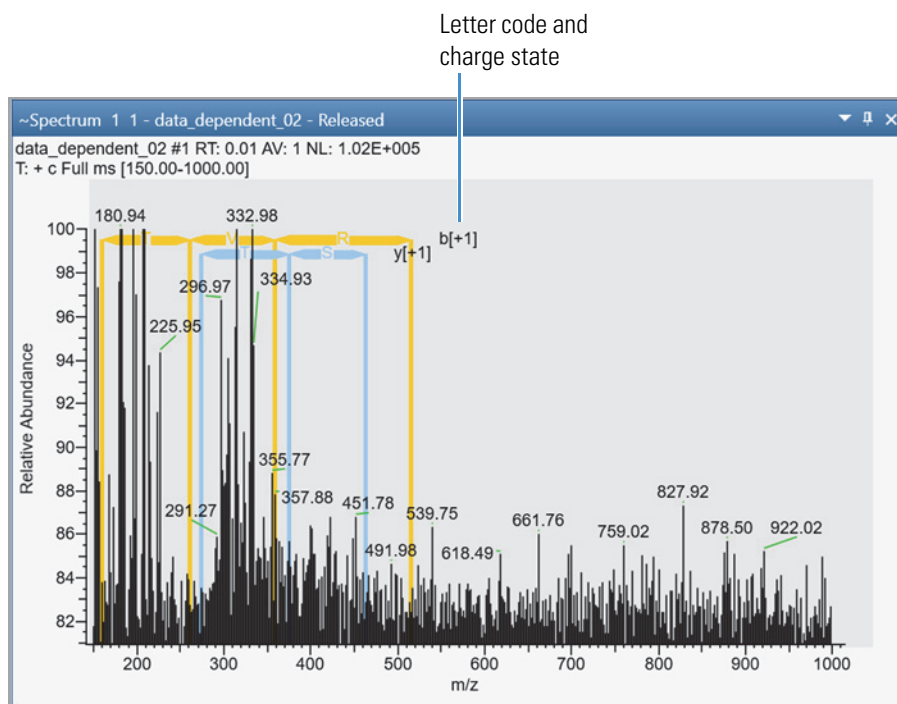
11 Peptide Fragment Annotation

Peptide Results View

The header of the Peptide Fragments table displays the terminal amino acid(s) or N-terminus for each ion; for example, (NH₃⁺) - A>. Each entry in the grid displays the m/z value of the indicated ion.

The Spectrum view displays vertical and horizontal color-coded lines to indicate the position of amino acid ions. The vertical color-coded bars indicate the position of matched ions and horizontal lines indicate the amino acids matched between adjacent ions. Each horizontal line also displays the ion series letter code and selected charge state.

Figure 164. Spectrum view



The following table describes the fragment ion series used in the activation type to annotate the spectra.

Table 54. Fragment ion series (Sheet 1 of 2)

Activation type	Ion series	Description
CID	<i>b</i>	b ion with charge on the N-terminal side
CID	<i>b</i> [*]	b ion that has lost ammonia (-17 Da)
CID	<i>b</i> ⁰	b ion that has lost water (-18 Da)
CID	<i>y</i>	y ion with charge on the C-terminal side

Table 54. Fragment ion series (Sheet 2 of 2)

Activation type	Ion series	Description
CID	y*	y ion that has lost ammonia (-17 Da)
CID	y ^o	y ion that has lost water (-18 Da)
CID	Immonium	Immonium ion
ETD	a•	a ion
ETD	z•	z ion
ETD	c'	c ion
ETD	y'	y ion

In Peptide Fragments table, colors have the following significance:

- White indicates that the peptide was not found.
- Yellow represents a series fragment ions.
- Gold represents b series fragment ions.
- Blue represents y series fragment ions.
- Orange represents c series fragment ions.
- Green represents z series fragment ions.

By default, the Peptide Fragments table displays all ion types searched in the analysis. You can remove any of these by clearing the appropriate ion check box on the right.

Filtering the Fragmented Ions

After you apply the parameters given in the Peptide Fragments Info Bar page, the Peptide Fragments table and Spectrum view displays all ion types based on the parameters defined. You can use the Display Options to filter the data.

❖ To apply filters in the display options

1. In the Ion Series area, select or clear the check box for the appropriate ions in the Ion Series area. In CID activation type, b and y ions are selected by default. In ETD activation type, c and z ions are displayed by default.
2. In the Charge State box, select **All** or specify a charge state from the dropdown to filter the charge states.

The Spectrum view and Peptide Fragments table display changes based on the filters selected.

Viewing Experiment and Instrument Information

In addition to containing the acquired data, the raw data file contains information about the scan header, instrument status, instrument method, tune method, and sample. Use the FreeStyle application to extract this information from the raw data files. Viewing the experiment information for a particular acquisition can help you interpret the results of that acquisition. Viewing the mass spectrometer readback (MS Trending detector type) information for the acquisition can help you troubleshoot anomalies in the data.

To review the experiment and instrument information in your raw data files, follow these topics.

Contents

- [Workspace Options Toolbar – Report Area](#)
- [Spectrum List – Display Options Toolbar](#)
- [Scan Header View](#)
- [Status Log View](#)
- [Instrument Method View](#)
- [Tune Method View](#)
- [Sample Information View](#)
- [File Header View](#)
- [Error Log View](#)

Workspace Options Toolbar – Report Area

Use the buttons in the Report area of the Workspace Options toolbar to view information that is contained in the raw data file.

❖ To display the Workspace Options toolbar

Click the **Workspace Options** toolbar tab.

Figure 165 shows the Report area, and Table 55 describes the toolbar buttons, which display information-specific views.

Figure 165. Report area of the Workspace Options toolbar

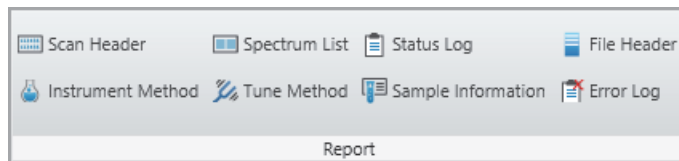


Table 55. Workspace Options toolbar – Report area buttons

Button	Description
Scan Header	<p>Displays the Scan Header View, showing the scan header of the active raw data file.</p> <p>Click the chromatogram trace to display the scan header for the chosen retention time and scan number.</p>
Spectrum List	<p>Displays the Spectrum List View, showing spectral peak information in a table.</p> <p>Click the chromatogram trace to display the spectrum list for the chosen retention time and scan number.</p>
Status Log	<p>Displays the Status Log View, showing mass spectrometer readbacks.</p> <p>Click the chromatogram trace to display the status log for the chosen retention time and scan number.</p>
Instrument Method	<p>Displays the Instrument Method View, showing the instrument method parameters that the instrument used to obtain the raw data file.</p>
Tune Method	<p>Displays the Tune Method View, showing the tune parameters that the instrument used to obtain the raw data file.</p>
Sample Information	<p>Displays the Sample Information View, showing sample-specific information.</p>
File Header	<p>Displays the File Header View.</p>
Error Log	<p>Displays the Error Log View, showing errors recorded during data acquisition.</p>

Spectrum List – Display Options Toolbar

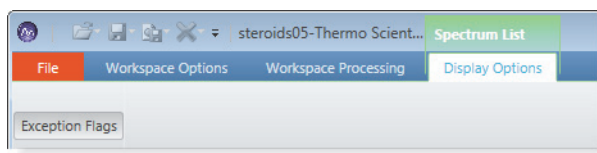
Use the Exception Flags button in the Spectrum List Display Options toolbar to add or remove the Flags column.

❖ To display the Spectrum List Display Options toolbar

1. Click the **Spectrum List** view to select it.
2. Click the **Display Options** toolbar tab.

Figure 166 shows the Spectrum List Display Options toolbar.

Figure 166. Spectrum List Display Options toolbar



The Exception Flags button is the only button in the Spectrum List Display Options toolbar. It displays or hides the Flags column in the Spectrum List view.

Scan Header View

Use the Scan Header view to display scan parameters and instrument data for the retention time and scan number that you select in the Chromatogram view.

❖ To open the Scan Header view

1. Click the **Workspace Options** toolbar tab.
2. Click **Scan Header**.

Figure 167 shows an example of a Scan Header view.

Figure 167. Scan Header view

Name	Value
Total Ion Current:	177517000
Scan Low Mass:	100
Scan High Mass:	1000
Scan Start Time (Min):	0
Scan Number:	1
Base Peak Intensity:	52672800
Base Peak Mass:	195.09
Scan Mode:	FTMS (1,1) + p ESI Full lock ms [100.00-1000.00]
===== Instrument data =====	
AGC	True
Micro Scan Count	1
Scan Segment	1
Scan Event	1
Ion Injection Time (ms)	21.9808921813965
Max. Ion Time (ms)	50
FT Resolution	50000
AGC Target	1000000
Analyzer Temperature	33.7397919725252
=== Mass Calibration: ===	

The Scan Header view can include the following information:

- Scan parameters
- Instrument data
- Mass calibration parameters
- Ion optics readbacks
- Diagnostic data

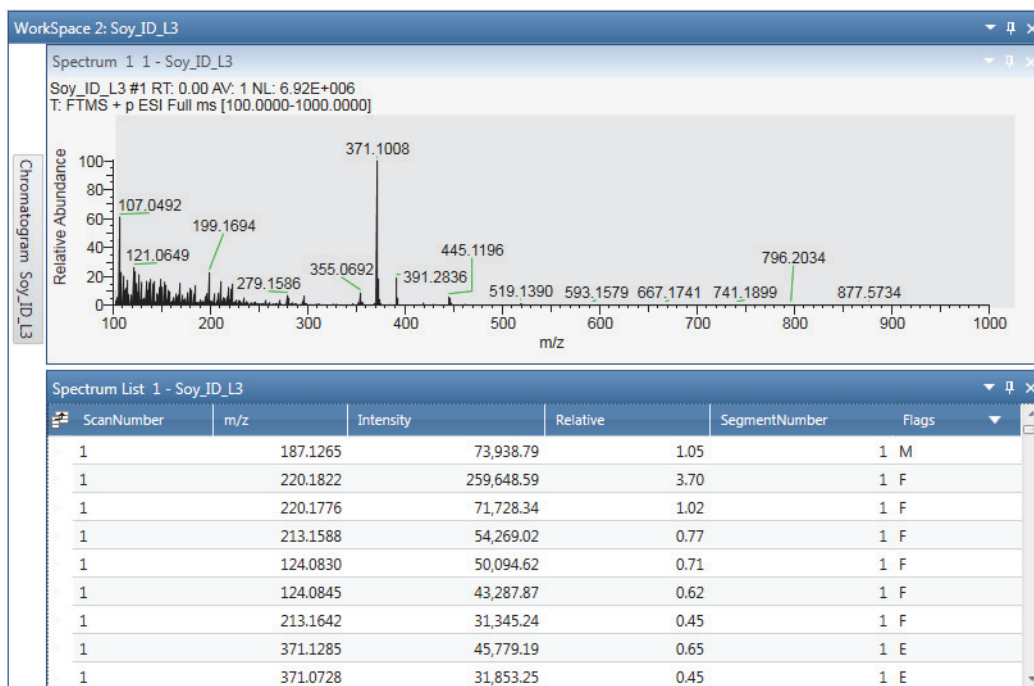
Spectrum List View

The Spectrum List view lists in tabular form the peak positions, intensities, and relative intensities of the peaks in the Spectrum view for the retention time that you select in a Chromatogram view.

❖ To open the Spectrum List view

1. Click the **Workspace Options** toolbar tab.
2. Click **Spectrum List**.

Figure 168 shows an example of a Spectrum List view and the associated mass spectrum.

Figure 168. Spectrum List view (bottom) associated with the mass spectrum (top)

Some raw data files contain flags that provide supplemental information about the peak data. The possible flags are as follows:

- R: Reference peaks are peaks from a reference compound used for an internal recalibration of a scan.
- E: Exception peaks are also peaks from a reference compound, but they are not used for recalibration. These are typically small isotopes or fragments of the main references.
- M: Merged peaks are peaks where the centroider combined two nearby peaks.
- F: Fragmented peaks are peaks separated into multiple peaks by the centroiding activity.

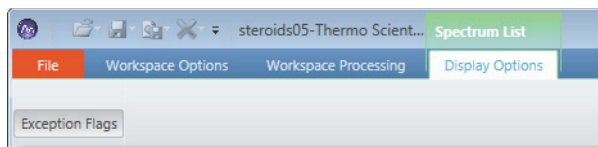
If the Spectrum List view does not show the Flags column (Figure 168), you can click Exception Flags in the Display Options toolbar for the Spectrum List view to see this column and the letters representing the flags.

❖ To display flags in the spectrum list

Do one of the following:


- Click the **Spectrum List** view to select it.
- Click the **Display Options** toolbar tab.

- c. Click **Exception Flags**.



The Flags column appears as the last column. To hide the Flags column, click **Exception Flags** again.

–or–

- a. Click the **Field Chooser** icon, , to the left of the first column heading.
- b. In the Field Chooser dialog box, select the **Flags** check box.

Status Log View

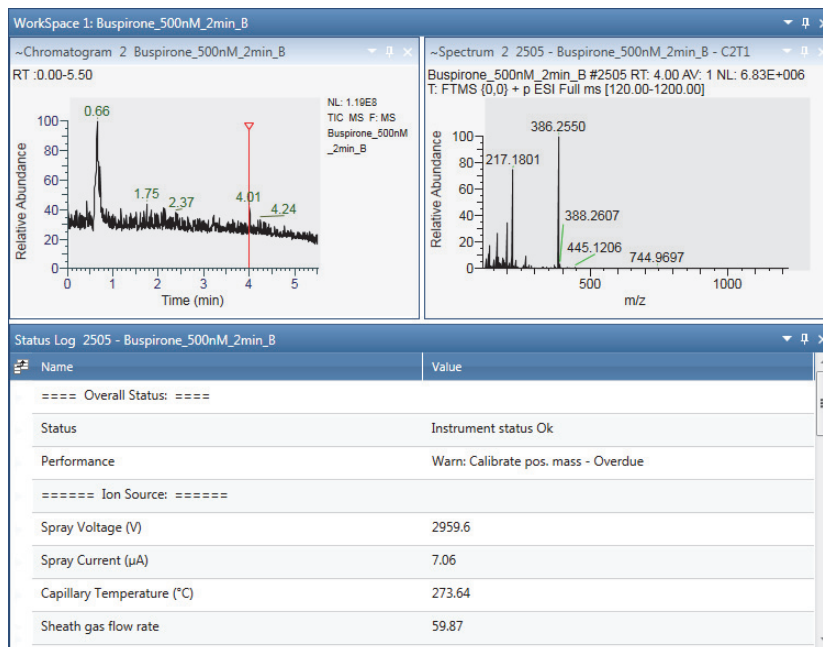
The Status Log view displays the mass spectrometer component readbacks that the system recorded at the retention time selected in the Chromatogram view.

❖ To open the Status Log view

1. Click the **Workspace Options** toolbar tab.
2. Click **Status Log**.

Figure 169 shows an example of the Status Log view. The view's title bar includes the scan number that corresponds to the selected retention time.

Figure 169. Status Log view



You can display individual instrument status readbacks, as a function of the retention time, in the Chromatogram view (see [Setting up Instrument Status Traces](#)). The readback display can aid in troubleshooting anomalies in the data.

Instrument Method View

The Instrument Method view displays the instrument method. An instrument method is a set of experiment parameters that define operating settings for the autosampler, liquid chromatograph (LC), mass spectrometer, divert valve, syringe pump, and so on. Instrument methods are saved as file type .meth.

For an LC/MS system, the Instrument Method view displays the instrument method settings for the autosampler, LC pump, and mass spectrometer on separate pages. This information does not change from scan to scan.

❖ To open the Instrument Method view

1. Click the **Workspace Options** toolbar tab.
2. Click **Instrument Method**.

For example, the Autosampler page lists the injection-cycle information.

[Figure 170](#) shows an example of the Autosampler page of an Instrument Method view.

Figure 170. Autosampler page of the Instrument Method view

Click these links to see more information about these instruments.

WorkSpace 1: Apple_01

Instrument Method - Apple_01

Thermo Pal Accela 1250 Pump Exactive - Orbitrap MS

Description:
Duplication of PAL local LC-Inj cycle.
This is only a sample method for short runs. For analytical work the method parameters must be set according to the specific needs.

Syringe: 100ul

Macro Sequence

LC-Inj (Step 1 of 1)
Air Volume (ul): 0
Pre Clean with Solvent 1: 0
Pre Clean with Solvent 2: 0
Pre Clean with Sample: 0
Filling Speed (ul/s): 1
Filling Strokes: 3
Inject to: LC Vlv1
Injection Speed (ul/s): 10
Pre Inject Delay (ms): 500
Post Inject Delay (ms): 500
Post Clean with Solvent 1: 3
Post Clean with Solvent 2: 3
Valve Clean with Solvent 1: 3
Valve Clean with Solvent 2: 3

Recommended Injection Volume
Volume: 1.25

Injection Look Ahead
Delay Time: 0.00

Macros used within method

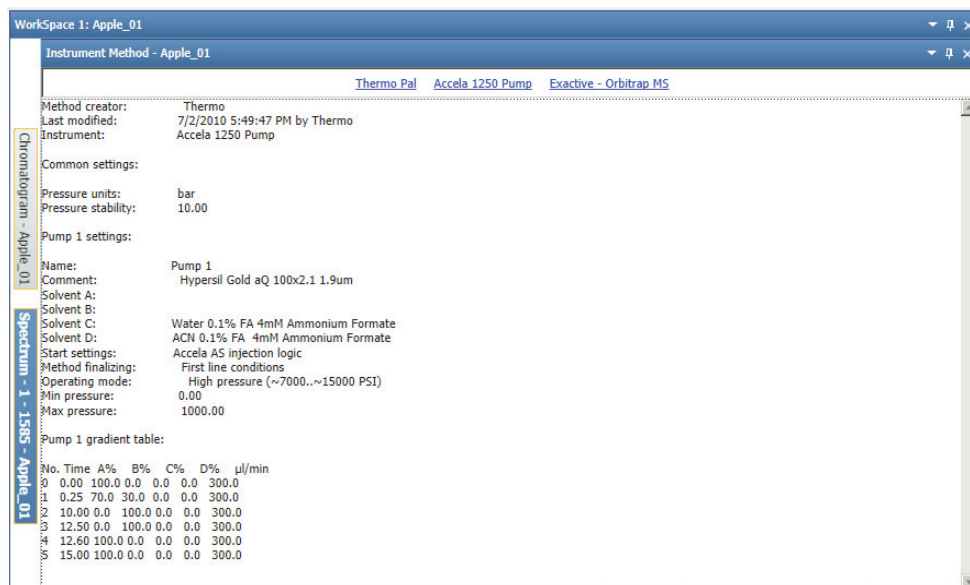
Macro name: LC-Inj
COMPLETE CYCLE

Duplication of PAL's local LC injection cycle.
Tray Name, Sample Number and Sample Volume are taken from the sample list.

The LC page can include the solvent composition and the gradient information for the acquisition.

Figure 171 shows an example of the LC page of an Instrument Method view.

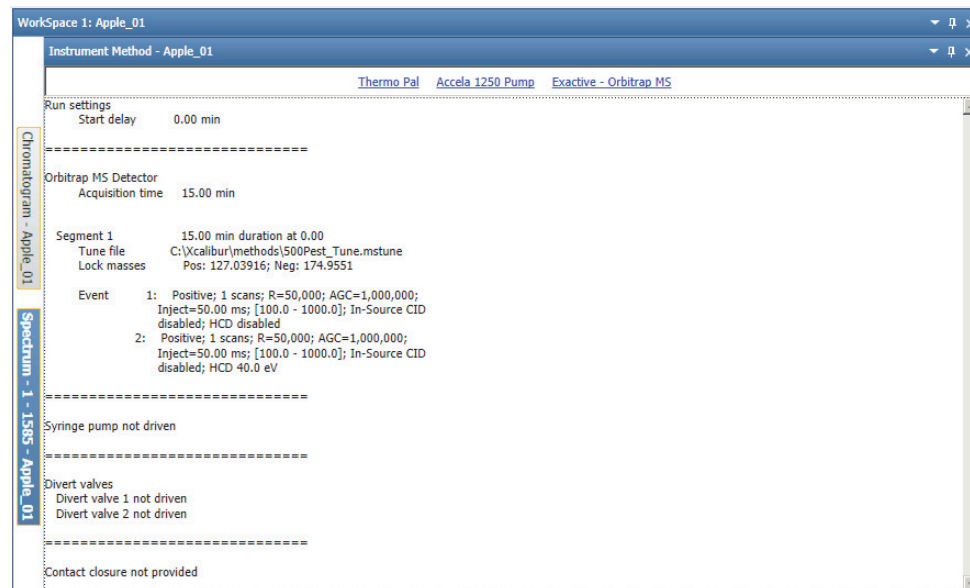
Figure 171. LC page of the Instrument Method view



The Mass Spectrometer page might list the tune file, the lock masses, the scan event descriptions, the activation types, the syringe pump settings, and the divert valve settings.

Figure 172 shows an example of the Mass Spectrometer page of the Instrument Method view.

Figure 172. Mass Spectrometer page of the Instrument Method view



Tune Method View

The Tune Method view displays the settings in the tune method. A tune method is a defined set of mass spectrometer tune parameters for the ion source, ion optics, and mass analyzer.

❖ To open the Tune Method view

1. Click the **Workspace Options** toolbar tab.
2. Click **Tune Method**.

Figure 173 shows an example of a Tune Method view.

Figure 173. Tune Method view

Name	Value
Capillary Temp (C):	175.00
APCI Vaporizer Temp (C):	500.00
Source Voltage (kV):	8.00
Source Current (uA):	6.00
Sheath Gas Flow (l):	90.00
Aux Gas Flow (l):	10.00
Capillary Voltage (V):	5.00
Tube Lens Offset (V):	55.00
Octapole RF Amplifier (Vp-p):	400.00
Octapole 1 Offset (V):	-2.25
Octapole 2 Offset (V):	-5.50
InterOctapole Lens Voltage (V):	-16.00
Trap DC Offset Voltage (V):	-10.00
Multiplier Voltage (V):	0.00
Maximum Ion Time (ms):	400.00
Ion Time (ms):	5.00
Data Type:	Centroid

Sample Information View

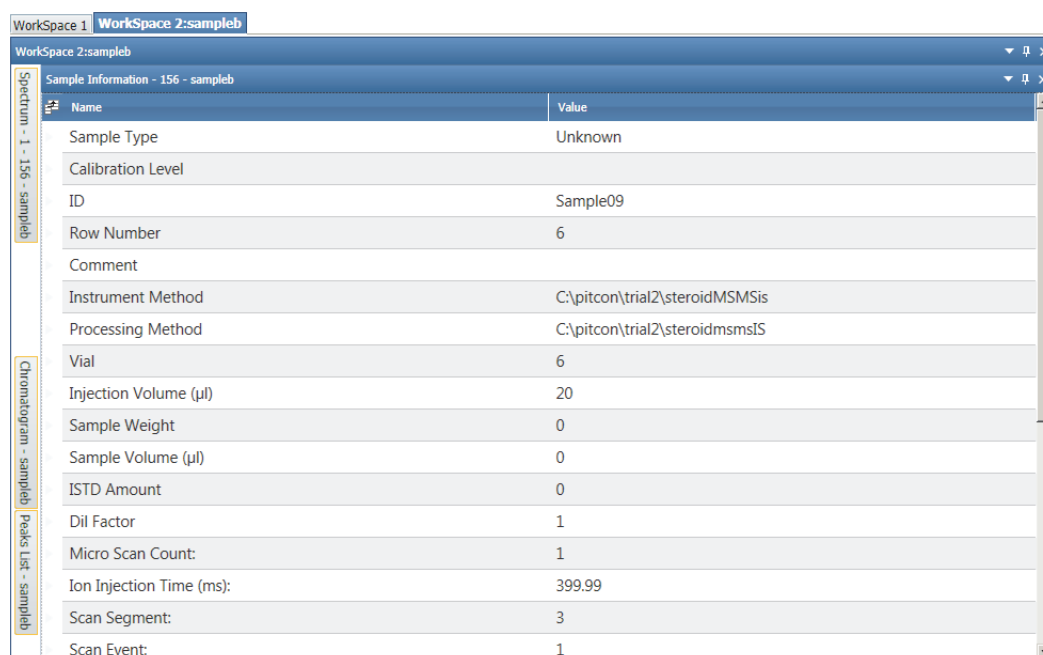
Use the Sample Information view to view sample-specific autosampler, LC pump, and mass spectrometer information in a raw data file. This information does not change from scan to scan.

❖ To open the Sample Information view

1. Click the **Workspace Options** toolbar tab.
2. Click **Sample Information**.

Figure 174 shows an example of a Sample Information view.

Figure 174. Sample Information view



Name	Value
Sample Type	Unknown
Calibration Level	
ID	Sample09
Row Number	6
Comment	
Instrument Method	C:\pitcon\trial2\steroidMSMSis
Processing Method	C:\pitcon\trial2\steroidmsmsIS
Vial	6
Injection Volume (µl)	20
Sample Weight	0
Sample Volume (µl)	0
ISTD Amount	0
Dil Factor	1
Micro Scan Count:	1
Ion Injection Time (ms):	399.99
Scan Segment:	3
Scan Event:	1

File Header View

Use the File Header view to view acquisition information about the current raw data file. The File Header view displays information from the acquisition sequence, the autosampler, and the mass spectrometer.

❖ To open the File Header view

1. Open a raw data file.
2. In the Workspace Options toolbar, click **File Header**.

The File Header view opens above the other views.

Figure 175. File Header view

Label1	Value1	Value2
drugx_06.raw	5/4/1996 11:21:33 PM	9
drugx_06.raw	StdUpdate	STD=100 pg/mL;US=100 pg/mL

The view has three columns, each displays a custom label name and value as configured in the Heading Editor dialog box.

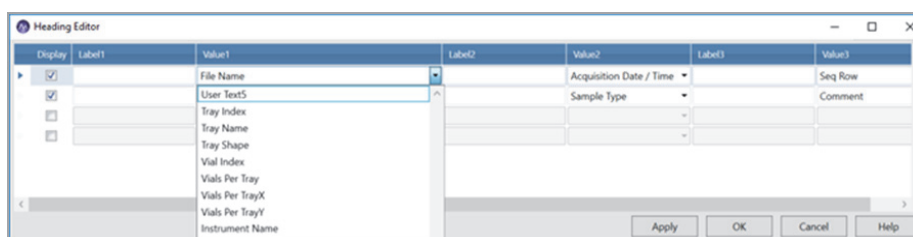
Editing the File Header Values

Use the Heading Editor dialog box to edit the values in the File Header view.

❖ To open the Heading Editor

1. Right-click the File Header, and click **Heading Editor**.

The Heading Editor dialog box opens.

Figure 176. Heading Editor dialog box

2. In the Heading Editor dialog box, choose a label name and value from the dropdown list for each column.

3. Click **Apply**.

The selected values are populated in the File Header.

4. Click **OK**.

Table 56 describes the information in the Heading Editor dialog box.

Table 56. Heading Editor (Sheet 1 of 3)

Read-only parameter	Description
Acquisition Date/Time	Displays the date and time that the displayed data was acquired.
Sample Name	Displays the text in the Sample Name column.
Comment	Displays the text in the Comment column.
Seq Row	Displays the number of the sequence row.

Table 56. Heading Editor (Sheet 2 of 3)

Read-only parameter	Description
Sample Type	Displays the sample type selection in the Sample Type column.
Path	Displays the directory and file name in the Path column, which is the file's original storage location and file name.
Cal Level	Displays the calibration level in the Level column for calibration standards.
Injection Volume	Displays the injection volume, in microliters, in the Inj. Vol. column.
Sample Weight	Displays the numeric value in the Sample Wt column. The application does not use this value for any internal calculations.
Sample Volume	Displays the numeric value in the Sample Vol column. The application does not use this value for any internal calculations.
Sample ID	Displays the sample identification text in the Sample ID column.
ISTD Amount	Displays the numeric value in the ISTD Corr Amt column.
CD Factor	Displays the numeric value in the Dil Factor column. The application does not use this value for any internal calculations.
Bar Code	Displays the bar code information acquired by the autosampler's bar code reader during an injection sequence.
Bar Code Status	Indicates whether the autosampler read the bar code for an injection vial.
Inst Method	Displays the instrument method in the Inst Meth column.
Proc Method	Displays the processing method in the Proc Meth column.
User Text 1 to User Text 5	Displays the text in the respective sequence table columns. An Xcalibur sequence table can include up to five additional text columns for user-specific information. The default headings for these columns are (1) Study, (2) Client, (3) Laboratory, (4) Company, and (5) Phone. You can customize the column headings in the Xcalibur data system's User Labels dialog box.
The following values are from the autosampler. Not all autosamplers supply this information to the data system.	
Tray Index	Displays text that identifies the autosampler tray.
Tray Name	Displays text that identifies the type of autosampler tray.
Tray Shape	Displays the shape of the autosampler tray (for example, rectangular).
Vial Index	Displays a numeric value.
Vials Per Tray	Displays a numeric value.

Table 56. Heading Editor (Sheet 3 of 3)

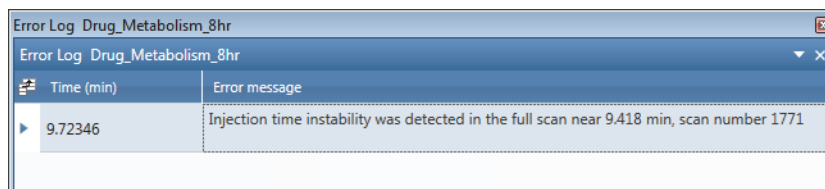
Read-only parameter	Description
Vials Per Tray X	Displays the number of vials across the autosampler tray.
Vials Per Tray Y	Displays the number of vials that the autosampler is deep.
Instrument Name	Displays the name of the mass spectrometer series (for example, Orbitrap Fusion).
Instrument Model	Displays the model name of the mass spectrometer (for example, Orbitrap Fusion).
Instrument Number	Displays the serial number of the mass spectrometer.
Instrument Software Version	Displays the version number of the mass spectrometer's instrument control software that is installed on the data system computer.
Instrument Hardware Version	Displays the version number of the hardware components that are installed on the mass spectrometer.
Flags	Displays additional information about the scan data.
Mass Tolerance	Displays the default mass tolerance setting on the Mass Options page of the Xcalibur Configuration dialog box. The FreeStyle application does not use this setting.
<p>Tip To open the Xcalibur Configuration dialog box, from the Xcalibur RoadMap view, choose Tools > Configuration.</p>	
Created by	Displays the user name.
File Name	Displays the name and path of the file storing the displayed data.

Error Log View

Use the error log to view errors recorded during data acquisition.

❖ To open the Error Log view

1. Open a raw data file.
2. Click the **Workspace Options** toolbar tab.
3. In the Reports area, click **Error Log**.

Figure 177. Error Log view

Deconvolving and Deisotoping Spectra with the Xtract Algorithm

The Xtract algorithm of the FreeStyle application uses a fitting scheme to deconvolve and deisotope isotopically resolved mass spectra of peptides, proteins, and nucleotides.

To deisotope and deconvolve the spectra for peptides, proteins, and nucleotides, follow these topics.

Contents

- [Understanding the Xtract Algorithm](#)
- [Applying the Xtract Algorithm to an Isotopically Resolved Spectrum](#)
- [Xtract Page](#)
- [Deconvolved Spectrum View](#)
- [Xtract Results View](#)

Understanding the Xtract Algorithm

The Xtract algorithm first examines a cluster of isotopically resolved peaks and uses the peak spacing of a cluster to determine an initial estimate of the mass of the relevant component. For peptides and proteins, it fits an average² distribution to the observed peak profile in that cluster to determine the monoisotopic mass that best reproduces that profile. (The monoisotopic mass is the mass of an ion for a given molecular formula, which is calculated by using the exact mass of the most abundant isotope of each element, for example, C = 12.000000, H = 1.007825, and O = 15.994915.)

² Senko, M.W.; Beu, S.C.; McLafferty, F.W. Determination of monoisotopic masses and ion populations for large biomolecules from resolved isotopic distributions. *J. Am. Soc. Mass Spectrometry*. **1995**, Vol. 6, 226–233.

13 Deconvolving and Deisotoping Spectra with the Xtract Algorithm

Applying the Xtract Algorithm to an Isotopically Resolved Spectrum

For nucleotides, the algorithm fits a distribution that is typical for nucleotides. Finally, the Xtract algorithm combines the results for all observed charge states for each mass component to produce a single mass value for that component. The resulting spectrum shows only the monoisotopic masses and their ion populations for the components that the algorithm identified.

When used properly, the Xtract algorithm reduces spectral noise and provides a high-intensity mass spectrum of the monoisotopic peaks. You can use the Xtract Results table, which contains the monoisotopic mass list of the deconvolved mass-spectral peaks or the extracted deconvolved spectrum, as the input to various search engines.

Figure 178 shows an isotopically unresolved mass spectrum. The mass-spectral peaks represent different charge states.

Figure 178. Isotopically unresolved mass spectrum of myoglobin

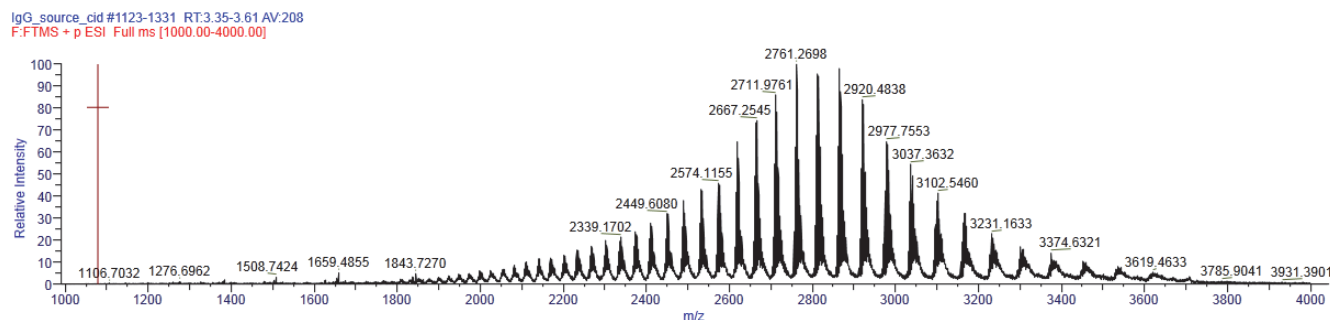
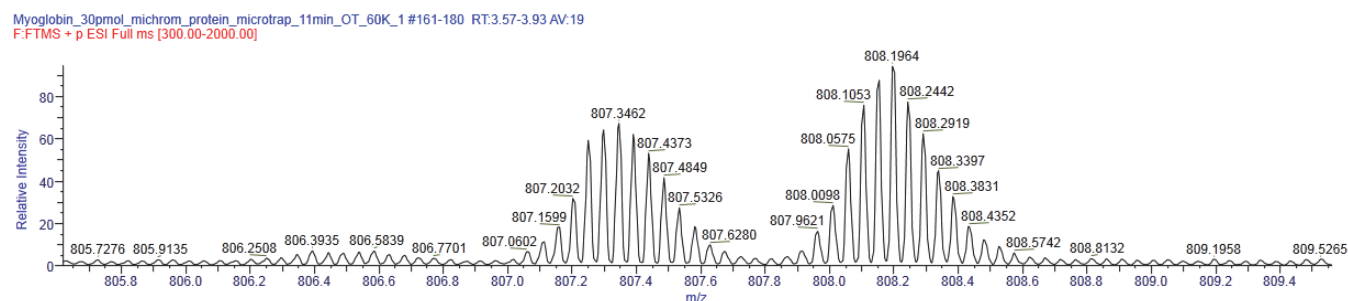


Figure 179 shows an isotopically resolved mass spectrum for one charge state.

Figure 179. Isotopically resolved mass spectrum of myoglobin for one charge state showing two components



Applying the Xtract Algorithm to an Isotopically Resolved Spectrum

❖ To apply the Xtract algorithm to an isotopically resolved spectrum

1. Review the default settings for the Xtract algorithm on the Xtract Options page of the Default Options Configuration dialog box as follows:
 - a. In the Workspace Options toolbar, click **Default Options**.
 - b. In the left pane, select **Xtract Options**.

- c. If needed, modify the settings and click **Save**.
2. Open a raw data file.
3. Select a mass spectrum.
4. Click the **Workspace Processing** tab.
5. Click **Xtract**.

The Xtract page opens in the Info Bar. The m/z Range boxes are populated with the m/z range of the selected spectrum.

6. If necessary, modify the m/z range by entering a range within the selected spectrum.
7. Review and modify the settings in the Deconvolution Parameters area as appropriate.
8. Click **Apply**.

The application runs the Xtract algorithm.

Xtract Page

Use the Xtract page in the Info Bar to specify the parameter settings for the Xtract algorithm.

❖ To display the Xtract page

1. Click the **Workspace Processing** toolbar tab.
2. Click **Xtract**.

[Figure 180](#) shows the default settings on the Xtract page. The m/z range depends on the selected spectrum.

Figure 180. Default settings on the Xtract page

The screenshot shows the Xtract Parameters dialog box with the following settings:

- Data Selection:** M/Z Range: 300.00 to 2000.00
- Deconvolution Parameters:**
 - Output Mass: M MH+
 - Adduct Element: H+ (1.00727663) K+ (38.9631585) Na+ (22.9892213) Custom
 - Charge Range: 5 to 50
 - Analyzer Type: OT
 - Rel. Abund. Threshold (%): 0
 - Isotope Table: Protein
 - Negative Charge:
 - Min Num Detected Charge: 3

Buttons at the bottom: Load Defaults, Apply

Note The default parameters on the Xtract page are for protein and peptide deconvolution in the positive polarity mode.

Table 57 describes the parameters for the Xtract page.

Table 57. Xtract page parameters (Sheet 1 of 5)

Parameter	Description
Data Selection	
<i>m/z</i> Range	<p>Specifies the portion of the input spectrum that the Xtract algorithm processes. The options are as follows:</p> <ul style="list-style-type: none"> Min: Specifies the lowest end of the input spectrum Max: Specifies the highest end of the input spectrum <p>For example, if the total mass range of the spectrum is mass 100–2000, a setting of 300–500 for the <i>m/z</i> Range parameter means that the Xtract algorithm processes only peaks with masses between <i>m/z</i> 300 and 500.</p> <p>Range: minimum <i>m/z</i> of the input spectrum to the maximum <i>m/z</i> of the input spectrum</p> <p>Default: minimum <i>m/z</i> of the input spectrum to the maximum <i>m/z</i> of the output spectrum</p>

Table 57. Xtract page parameters (Sheet 2 of 5)

Parameter	Description
Deconvolution	
Output Mass	<p>Determines whether the Xtract algorithm returns a single peak at either the monoisotopic mass or the monoisotopic MH^+ mass for each of the detected components. The options are as follows:</p> <ul style="list-style-type: none"> • M: Specifies that the results file contains a single peak for the monoisotopic mass for each of the detected components. This option generates masses without adducts. • MH^+: Specifies that the results file contains a monoisotopic MH^+ mass for each of the detected components. This option generates masses with adducts. <p>Default: M</p>
Adduct Element	<p>Specifies the adduct ions used during ESI processing. Adduct ions bring the charge to the molecule, and this charge converts it to an ion. The options are as follows:</p> <ul style="list-style-type: none"> • H^+ (1.00727663): Specifies that the adduct is hydrogen. • K^+ (38.9631585): Specifies that the adduct is potassium. • Na^+ (22.9892213): Specifies that the adduct is sodium. • Custom: Specifies that the adduct is a charge carrier other than hydrogen, potassium, or sodium. When you select this option, a box opens so that you can type the mass of the custom charge carrier. <p>Default: H^+</p>
Charge Range	<p>Specifies the lowest and highest charge state to be deconvolved. The options are as follows:</p> <ul style="list-style-type: none"> • Low: Specifies the lowest charge state. • High: Specifies the highest charge state. <p>For example, if you set this parameter range to 1–5, the Xtract algorithm considers only charge states 1 through 5 for deconvolution. It ignores charge states 6 and higher.</p> <p>Default range: 5–50</p>

Table 57. Xtract page parameters (Sheet 3 of 5)

Parameter	Description
Analyzer Type	<p>Specifies the type of mass analyzer that was used to obtain the mass spectral data. Different types of mass analyzers have different mass accuracies. The options are as follows:</p> <ul style="list-style-type: none"> • FT: Specifies a Fourier transform - ion cyclotron resonance (FT-ICR) mass analyzer • OT: Specifies an Orbitrap™ mass analyzer • Sector: Specifies a magnetic-sector mass analyzer • Quad: Specifies a quadrupole mass analyzer <p>Default: OT</p>
Rel. Abund. Threshold (%)	<p>Specifies a threshold (as a percentage) below which the Xtract algorithm filters out data for data reporting.</p> <p>This option sets a relative threshold as a percentage of the most abundant component in the spectrum. The most abundant peak in the deconvolved spectrum has a relative abundance of 100 percent, and all other peaks are calculated relative to that one.</p> <p>In the Xtract Results table of the Xtract Results view, the application shows only those components that are greater than or equal to this relative abundance threshold in the deconvolved spectrum. For example, if the highest peak has an absolute abundance of 1000, the relative abundance is 1 percent, and no peaks below an absolute abundance of 10 appear in the deconvolved spectrum.</p> <p>Zero (0) displays all results, and 100 displays only the most abundant component.</p> <p>Range: 0–100; default: 0</p>

Table 57. Xtract page parameters (Sheet 4 of 5)

Parameter	Description
Isotope Table	<p>Specifies the type of isotope table to use.</p> <p>Isotope tables simulate the distribution of isotopic peaks, in m/z, for different choices of the monoisotopic mass. The Xtract algorithm chooses the monoisotopic mass with the best fit between the theoretical and the observed isotope distribution.</p> <p>To generate an isotope table, the Xtract algorithm uses a chemical formula to describe the type of molecule. You can choose one of the following formulas:</p> <ul style="list-style-type: none"> • Protein: Uses an average formula typical for peptides and proteins to generate the isotope table • Nucleotide: Uses an elemental formula typical for nucleotides to generate the isotope table <p>Default: Protein</p>
Negative Charge	<p>Indicates whether the data was acquired in positive charge mode or negative charge mode during the ESI process.</p> <p>You might want to use this option when you process compounds that contain nucleotides like those found in DNA and RNA. When you acquire these compounds in negative mode, the resulting mass spectra are often clearer. Deprotonation of nucleotides, which are acidic, occurs when the compound is dissolved in a basic solution and negative voltage is applied to produce negatively charged ions. The options are as follows:</p> <ul style="list-style-type: none"> • Selected: The data was acquired in negative charge mode. • Cleared: The data was acquired in positive charge mode. <p>Default: Cleared</p> <p>IMPORTANT Do not select Negative Charge if your data was acquired in positive mode. The results will not be usable.</p>
Min Num Detected Charge	<p>Specifies the minimum number of charge states required to produce a component. No components with less than this minimum number appear in the deconvolved spectrum.</p> <p>Valid values: an integer greater than or equal to 1 Default: 3</p>

13 Deconvolving and Deisotoping Spectra with the Xtract Algorithm

Xtract Page

Table 57. Xtract page parameters (Sheet 5 of 5)

Parameter	Description
Button	
Load Defaults	Resets all the parameters on the Xtract page to the default settings.
Apply	Runs the Xtract algorithm on the selected spectrum.

Deconvolved Spectrum View

The Xtract algorithm transforms the source mass spectrum into a deconvolved mass spectrum and displays it in the Deconvolved Spectrum view, labeled with mass units rather than with the mass-to-charge ratio on the x axis. The spectrum that the application displays in the Deconvolved Spectrum view is a mass spectrum that consists of the monoisotopic peaks of the components that the Xtract algorithm detected. The range of the mass spectrum is from the mass of the component with the lowest monoisotopic mass to the mass of the component with the highest monoisotopic mass. Information about each component can be found in the Xtract Results view.

❖ To run the Xtract algorithm and display the Deconvolved Spectrum view

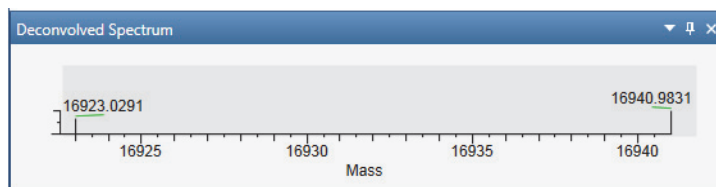
1. Select a mass spectrum for the Xtract algorithm to deconvolve.
2. Click the **Workspace Processing** toolbar tab.
3. Click **Xtract**.

The Xtract page opens in the Info Bar.

4. Set the deconvolution parameters on the Xtract page.
5. Click **Apply**.

Figure 181 shows the deconvolved spectrum for the myoglobin example.

Figure 181. Deconvolved Spectrum view showing the two monoisotopic peaks of myoglobin (located at the endpoints)



The spectrum in the [Spectrum View](#) or [MultiSpectrum View](#) shows the multiple masses (see [Figure 184](#)) used to create the deconvolved peak displayed in the Deconvolved Spectrum view.

Xtract Results View

The Xtract Results view displays the experimentally determined results for each monoisotopic mass and charge state.

Xtract deconvolves isotopically resolved mass spectra—that is, spectra in which it is possible to distinguish separate peaks for different isotopic compositions of the same component. When the Xtract algorithm finishes processing, it displays the deconvolved spectrum in the [Deconvolved Spectrum View](#) as a set of peaks in mass and relative intensity.

13 Deconvolving and Deisotoping Spectra with the Xtract Algorithm

Xtract Results View

It also displays the component list in the Xtract Results view as a list of masses, intensities, charge state information, and quality scores (Figure 182). The values in the columns represent the outputs of the deconvolution. Each peak in the list is composed of isotopic clusters. Each isotopic cluster in the original spectrum provides evidence for the peak in the deconvolved spectrum.

❖ To run the Xtract algorithm and display the Xtract Results view

1. Display the mass spectrum for the Xtract algorithm to deconvolve in the Spectrum or Multi Spectrum view.
2. Click the **Workspace Processing** toolbar tab.
3. Click **Xtract**.
4. Set the parameters on the [Xtract Page](#).
5. Click **Apply**.

Figure 182 shows the Xtract Results view for the two-component myoglobin example.

Figure 182. Xtract Results view in two parts (left and right) showing the 16 923 Da and 16 941 Da components

Xtract Results					
No.	Monoisotopic Mass	Sum Intensity	Number of Charge States	Average Charge	
+	1	16923.0291	572,776.58	9	20.96
+	2	16940.9831	850,259.11	12	20.56

Delta Mass	Relative Abundance	Fractional Abundance	RT Range
-17.95	67.3649	40.2503	3.654
0.00	100.0000	59.7497	3.654

Table 58 describes the parameters for the Xtract Results view.

Table 58. Xtract Results view parameters (Sheet 1 of 2)

Parameter	Description
No.	An integer, starting at one, that labels the components in the order of increasing monoisotopic mass.
Monoisotopic Mass	The weighted average of the monoisotopic masses of each charge state:

$$\text{Monoisotopic Mass} = \frac{\sum_i (\text{Monoisotopic Mass of This Charge} \times \text{Charge Normalized Intensity})}{\text{Sum Intensity}}$$

where i is the sequential order of the charges in the Charge column of the expanded Xtract table.

Table 58. Xtract Results view parameters (Sheet 2 of 2)

Parameter	Description
Sum Intensity	The sum of the intensities of all the charge states and the isotopes of a component.
Number of Charge States	The number of charge states that the Xtract algorithm used to calculate the monoisotopic mass.
Average Charge	The average of the charges weighted by intensity.
Delta Mass	The difference in mass (in daltons) of a component from the mass of the most intense component.
Relative Abundance	<p>Displays the components that are above the relative abundance threshold set by the Relative Abundance Threshold parameter on the Xtract page. The application assigns the largest peak in a deconvolved spectrum a relative abundance of 100 percent. An abundance number in the Relative Abundance column represents the intensity in the same row of the Sum Intensity column divided by the greatest intensity in the Sum Intensity column multiplied by 100.</p> <p>For example, if the largest peak in a deconvolved spectrum has an intensity of 1000, the application assigns it a relative abundance of 100 percent. If the next most abundant peak has an intensity of 500, the application assigns it an abundance of 50 percent:</p> $\frac{500}{1000} \times 100\% = 50\%$
Fractional Abundance	The fractional abundance of a component, which is the abundance for that peak relative to the total abundance of all peaks in the spectrum, expressed as a percentage. The sum of all fractional abundances of all peaks in a deconvolved spectrum is 100 percent.
RT Range	The retention time (in minutes), or the retention time range (if the retention time was averaged), that produced the mass spectrum that the Xtract algorithm deconvolved.

Click **+** to expand the Xtract Results view for a component. The expanded list shows the experimentally-determined parameters for the individual charge states of a component.

In the expanded list (see [Figure 183](#)), six new columns appear: Charge State, Calculated Monoisotopic m/z, Monoisotopic Mass for This Charge, Mostabund m/z, Charge Normalized Intensity, and Fit %. These values represent the isotopic clusters with different charge states from the source spectrum that were used to produce the peak in the deconvolved spectrum.

Figure 183 shows the Xtract Results view of Figure 182 with the 16 923 Da component expanded to show the individual charge states.

Figure 183. Xtract Results view with the 16 923 Da component expanded to show results for the nine individual charge states

Charge State	Calculated Monoisotopic m/z	Monoisotopic Mass for This Charge	Mostabund m/z	Charge Normalized Intensity	Fit %
18	941.1756	16923.0428	941.7327	83,056.26	89.10
19	891.6931	16923.0477	892.2206	105,669.90	89.20
20	847.1588	16923.0393	847.6607	90,557.22	87.60
21	806.8658	16923.0641	807.3449	73,625.01	91.80
22	770.2359	16923.0486	770.6924	64,146.69	84.10
23	736.7912	16923.0498	737.2283	64,615.90	92.20
24	706.1335	16923.0670	706.5534	47,677.56	86.80
25	677.9285	16923.0196	678.3286	27,615.76	86.40
26	651.8930	16923.0233	652.2789	15,812.27	86.10

Table 59 describes the parameters for the expanded Xtract Results view.

Table 59. Expanded Xtract Results view parameters

Parameter	Description
Charge State	The charge states of the individual isotopic clusters that constitute the total number shown in the Number of Charge States column.
Calculated Monoisotopic m/z	The mass-to-charge ratio of the monoisotopic mass that the Xtract algorithm calculated from the isotopic peak envelope for a specific charge state.
Monoisotopic Mass For This Charge	The mass (in daltons) that is equal to the calculated monoisotopic m/z multiplied by the charge state, z .
Mostabund m/z	The mass-to-charge ratio of the most abundant isotope, or the height of the tallest peak in the isotopic distribution.
Charge Normalized Intensity	The sum of the intensities of the isotopic peaks for a particular charge state.
Fit %	The quality of the match between a measured isotope pattern and an average distribution of the same mass. This column displays a value between 0 and 100 percent. <ul style="list-style-type: none"> 0 percent indicates a poor fit between the measured pattern and the average pattern. 100 percent indicates that the observed peaks in the measured isotope profile are absolutely identical to those in a theoretical average distribution and that any missing peaks fall below a restrictive threshold.

When you select any charge state row in the expanded Xtract Results view, the application automatically highlights in blue the following items:

- The blue spectral line in the [Spectrum View](#) or [MultiSpectrum View](#) corresponding to the most abundant m/z for the selected row, as shown in [Figure 184](#) and [Figure 185](#).

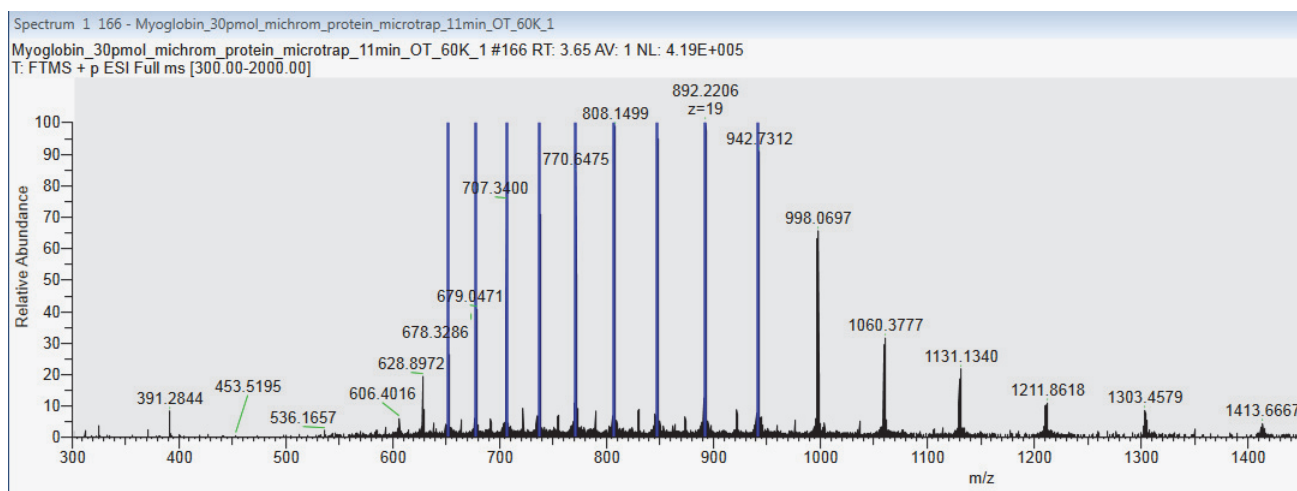
For example, in [Figure 185](#), a blue spectral line marks the value of 770.6924 corresponding to the [Mostabund \$m/z\$](#) value of the selected row.

- The deconvolved peak in the [Deconvolved Spectrum View](#).

You might have to right-click and choose Reset to Scale to see the blue spectral lines.

[Figure 184](#) shows the mass spectrum from [Figure 178](#) with nine blue bars that indicate the positions of the theoretical most abundant m/z for each of the nine charge states that the Xtract algorithm used to deconvolve the spectrum.

Figure 184. Mass spectrum of myoglobin showing the nine charge states (blue bars) that the Xtract algorithm used to calculate the 16923 Da monoisotopic mass

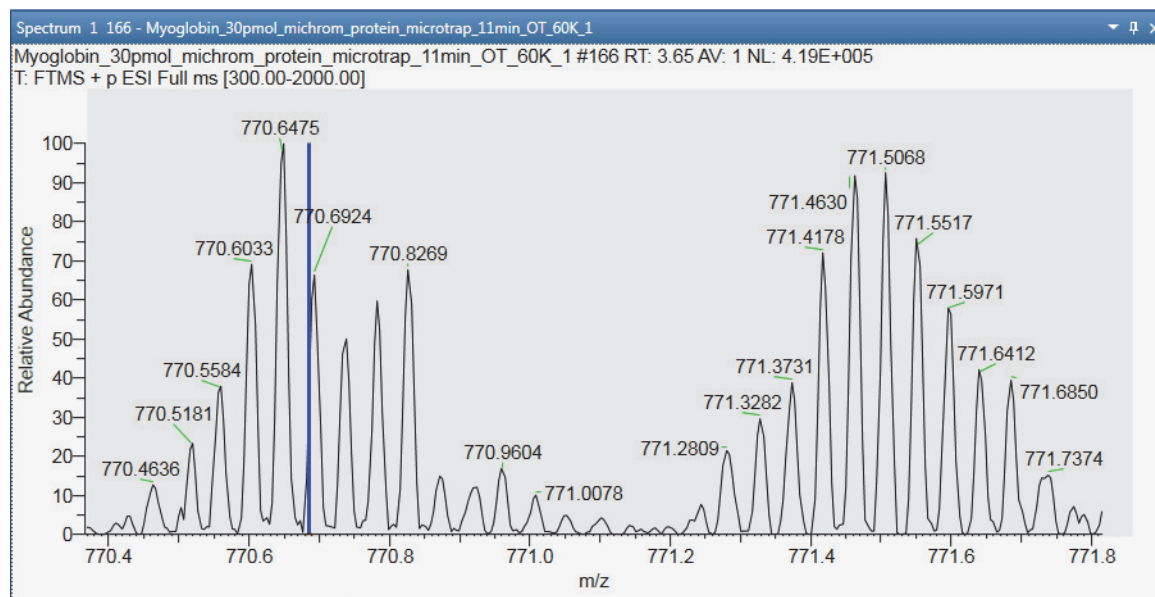


[Figure 185](#) shows the theoretical most abundant m/z for the $z=22$ charge state of the 16923 Da component.

13 Deconvolving and Deisotoping Spectra with the Xtract Algorithm

Xtract Results View

Figure 185. Mass spectrum at the $z = 22$ charge state showing the location (blue bar) of the theoretical most abundant m/z for the 16 923 Da monoisotopic mass



FreeStyle Default Settings

Use the pages of the Default Options Configuration dialog box to specify the initial default settings that the FreeStyle application uses when you start the application.

Follow these topics.

Contents

- [Default Mass Precision Page](#)
- [Default Peak Detection Page](#)
- [Default Recently Used Files Page](#)
- [Default Raw Data Files Directory Page](#)
- [Default Elemental Composition Page](#)
- [Factory Default Template Page](#)
- [Default Library Search Page](#)
- [Default Workspace Options Page](#)
- [Default Xtract Options Page](#)

Modifying, Saving, and Resetting the Default Configuration Options

Use the pages of the Default Options Configuration dialog box to modify some of the parameter settings for the FreeStyle application.

Follow these procedures:

- [To open the Default Options Configuration dialog box](#)
- [To display a specific page of the dialog box](#)
- [To open the Help topic for a specific page](#)
- [To save your new parameter settings](#)

A FreeStyle Default Settings

Modifying, Saving, and Resetting the Default Configuration Options

- To cancel all your changes to the parameter settings
- To return all the configuration parameters to their factory default settings
- To apply the factory default layout to the active Workspace view

❖ To open the Default Options Configuration dialog box

- Open the [File Menu](#) and click **Options** at the bottom right.

–or–

- In the [Workspace Options Toolbar](#), click **Default Options**.

The dialog box opens to the Default Mass Precision page.

❖ To display a specific page of the dialog box

In the left pane of the dialog box, click the page link.

❖ To open the Help topic for a specific page

Click **Help** at the bottom right of the page.

❖ To save your new parameter settings

Click **Save**.

The Default Options dialog box closes, and the application saves all the modified parameter settings.

❖ To cancel all your changes to the parameter settings

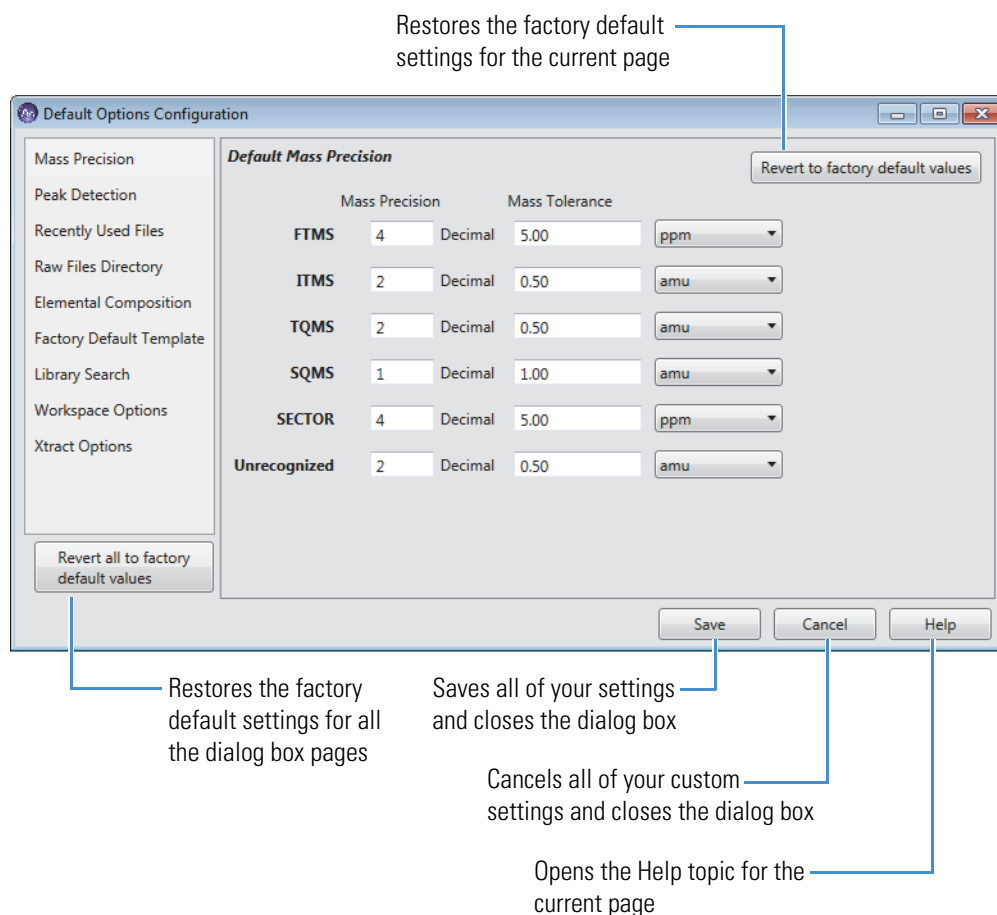
Click **Cancel**.

The Default Options Configuration dialog box closes without changing the parameter settings.

❖ To return all the configuration parameters to their factory default settings

At the bottom left, click **Revert All to Factory Default Values**.

Figure 186. Location of the Revert All to Factory Default Values button



❖ **To apply the factory default layout to the active Workspace view**

1. In the left pane of the dialog box, click **Factory Default Layout**.
2. At the top right of the [Factory Default Template Page](#), click **Revert to Factory Default Values**.

Default Mass Precision Page

Use the Default Mass Precision page to specify how many decimal places the application displays, by default, and also the default mass tolerance.

❖ **To display the Default Mass Precision page**

1. Open the Default Options Configuration dialog box (see [Modifying, Saving, and Resetting the Default Configuration Options](#)).
2. In the left pane, click **Mass Precision**.

Figure 186 shows the Default Mass Precision page with its original default values, and Table 60 lists the default Mass Precision page parameters.

Table 60. Default Mass Precision page parameters (Sheet 1 of 2)

Column	Description
Mass Precision	<p>Specify the default number of decimal places in the mass-to-charge ratios that the application displays in the Spectrum view and the MultiSpectrum view.</p> <p>Default:</p> <ul style="list-style-type: none"> • FTMS: 4 • ITMS: 2 • TQMS: 2 • SQMS: 1 • Sector: 4 • Unrecognized: 2 <p>Range: 0–5</p>
Mass Tolerance	<p>Specify the default mass tolerance.</p> <p>The mass tolerance defines how close a measured mass must be to the mass in the mass list to be considered the same mass.</p> <p>Default:</p> <ul style="list-style-type: none"> • FTMS: 5.00 ppm • ITMS: 0.50 amu • TQMS: 0.50 amu • SQMS: 1.00 amu • Sector: 5.00 ppm • Unrecognized: 0.50 amu <p>Range: 0.00–1000.00 ppm or 0.00–10.00 amu</p>
Default Mass Tolerance Units	<p>Specify the default mass tolerance units as follows:</p> <ul style="list-style-type: none"> • ppm: In parts per million. Select this option for FTMS and sector mass analyzers because the peak width is proportional to the mass-to-charge ratio. When you select ppm, the size of the window around the reference mass is proportional to the mass-to-charge ratio of the reference mass. • amu: In units of the mass-to-charge ratio. When you select amu, the size of the window around the reference mass remains the same regardless of the mass-to-charge ratio of the reference mass. • mmu: In units of one one-thousandth of an amu.

Table 60. Default Mass Precision page parameters (Sheet 2 of 2)

Column	Description
Button	
Revert to Factory Default Values	Returns the parameter settings on this page to their factory default values.
Help	Opens the Help topic for this page.

Default Peak Detection Page

Use the Default Peak Detection page to select the default peak detection algorithm and its default parameters. For a description of these parameters, see these topics: [Avalon Peak Detection Page](#), [Genesis Peak Detection Page](#), [ICIS Peak Detection Page](#), or [PPD Peak Detection Page](#).

❖ To display the Default Peak Detection page

1. In the Workspace Options toolbar, click **Default Options**.

The Default Options Configuration dialog box opens.

2. In the left pane, click **Peak Detection**.

To save your custom settings, see [Modifying, Saving, and Resetting the Default Configuration Options](#).

Default Recently Used Files Page

Use the Default Recently Used Files page to set the number of recently used files that the application displays in the Recent Items area of the File menu. By default, this number is set to 10. You can change the setting to a value from 0 to 10.

❖ To display the Default Recently Used Files page

1. In the Workspace Options toolbar, click **Default Options**.

The Default Options Configuration dialog box opens.

2. In the left pane, click **Recently Used Files**.

The Default Number of Recently Used Files page appears ([Figure 187](#)).

Figure 187. Default Number of Recently Used Files page

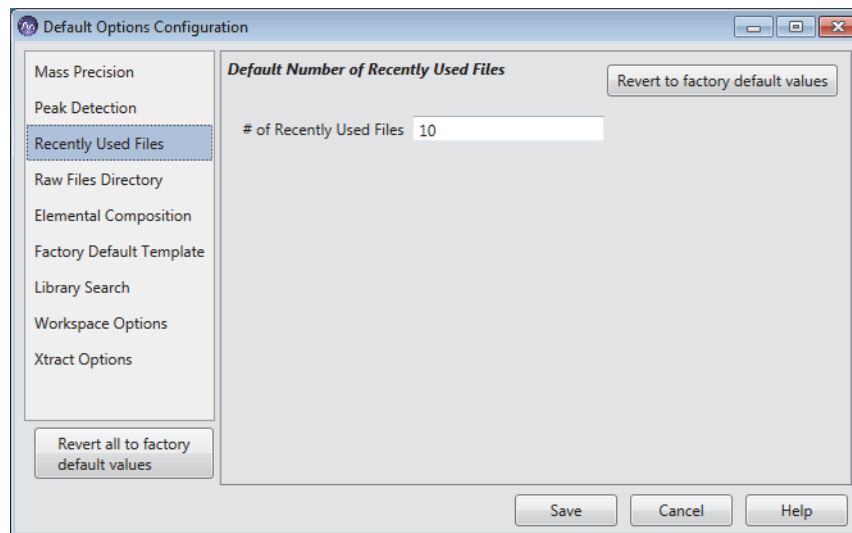


Table 61 describes the parameter and buttons on the Default Number of Recently Used Files page. To save a custom setting, see [Modifying, Saving, and Resetting the Default Configuration Options](#).

Table 61. Default Number of Recently Used Files page parameter and buttons

Parameter	Description
# of Recently Used Files	Specifies the maximum number of files that the application displays in the Recent Items area of the File menu. Range: 0 to 10; default: 10
Button	
Revert to Factory Default Values	Returns the parameter settings on this page to their factory default values.
Help	Opens the Help to the topic for this page.

Default Raw Data Files Directory Page

Use the Default Raw Data Files Directory page to specify the default location where the FreeStyle application looks for raw data files.

❖ To display the Default Raw Data Files Directory page

1. In the Workspace Options toolbar, click **Default Options**.
The Default Options Configuration dialog box opens.
2. In the left pane, click **Raw Files Directory** ([Figure 188](#)).

Figure 188. Default Raw Files Directory page

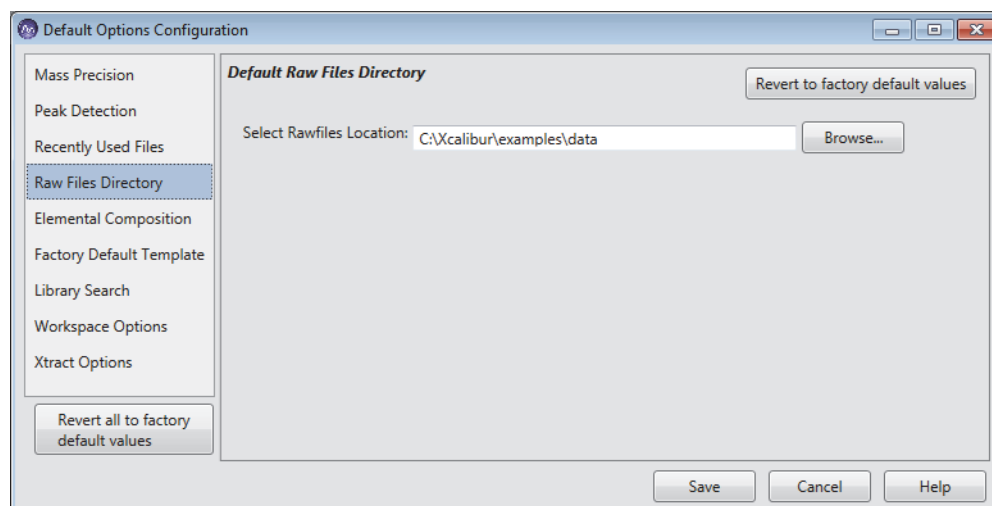


Table 62 describes the parameter and buttons on the Default Raw Files Directory page. To save a custom setting, see [Modifying, Saving, and Resetting the Default Configuration Options](#).

Table 62. Default Raw Files Directory page parameter and buttons

Parameter	Description
Select RawFiles Location	Default: C:\Xcalibur\examples\data
Button	
Revert to Factory Default Values	Returns the parameter settings on this page to their factory default values.
Help	Opens the Help to the topic for this page.

Default Elemental Composition Page

Use the Default Elemental Composition page to specify the default elemental composition search parameters. For information about the basic parameter settings and performing an elemental composition analysis, see [Elemental Composition Page](#). To save your custom settings, see [Modifying, Saving, and Resetting the Default Configuration Options](#). To use a custom periodic table for elemental composition analysis, see [Using Custom Periodic Table for Analysis](#).

❖ To display the Default Elemental Composition page

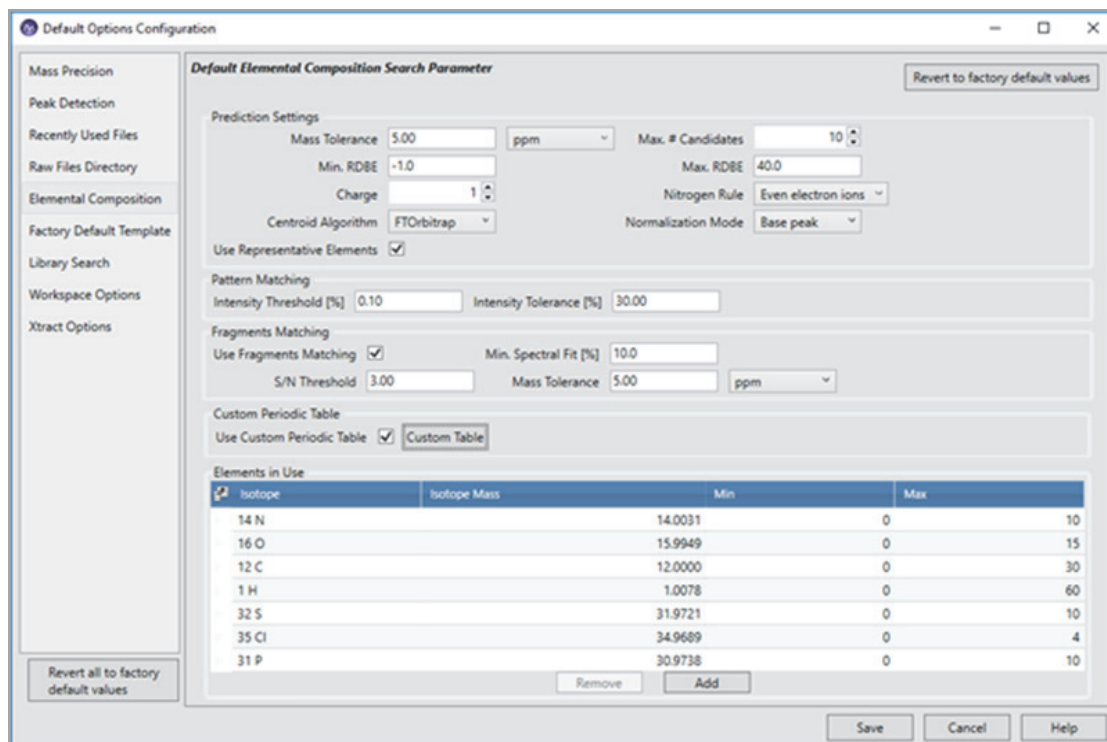
1. In the Workspace Options toolbar, click **Default Options**.

The Default Options Configuration dialog box opens.

2. In the left pane, click **Elemental Composition**.

Figure 189 shows the Default Elemental Composition Search Parameter page. For information about the parameter settings, see [Elemental Composition Page](#).

Figure 189. Default Elemental Composition Search Parameter page



Factory Default Template Page

Use the Factory Default Template page to apply the factory default layout to the active Workspace view (see [Modifying, Saving, and Resetting the Default Configuration Options](#)). The settings on this page are read only.

❖ To display the Factory Default Template page

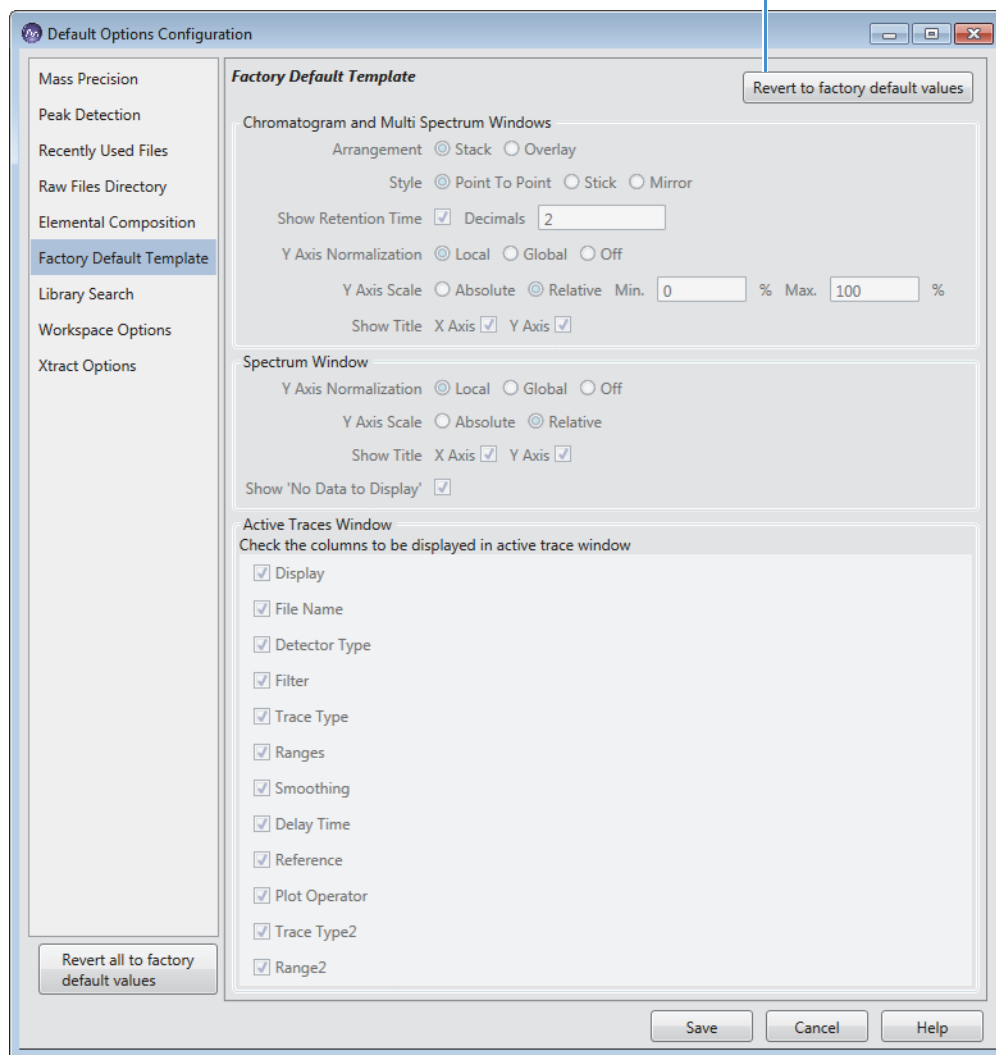
1. In the Workspace Options toolbar, click **Default Options**.

The Default Options Configuration dialog box opens.

2. In the left pane, click **Factory Default Template** (Figure 190).

Figure 190. Factory Default Template page

Click to apply the factory default layout to the active Workspace view.



Default Library Search Page

Use the Default Library Search page to specify the default library search parameters for the NIST libraries and the *mzVault* libraries. For a description of these parameters, see [Modifying a NIST Search from the NIST Search Page](#) and [Modifying an *mzVault* Search from the *mzVault* Search Page](#). To save your custom settings, see [Modifying, Saving, and Resetting the Default Configuration Options](#).

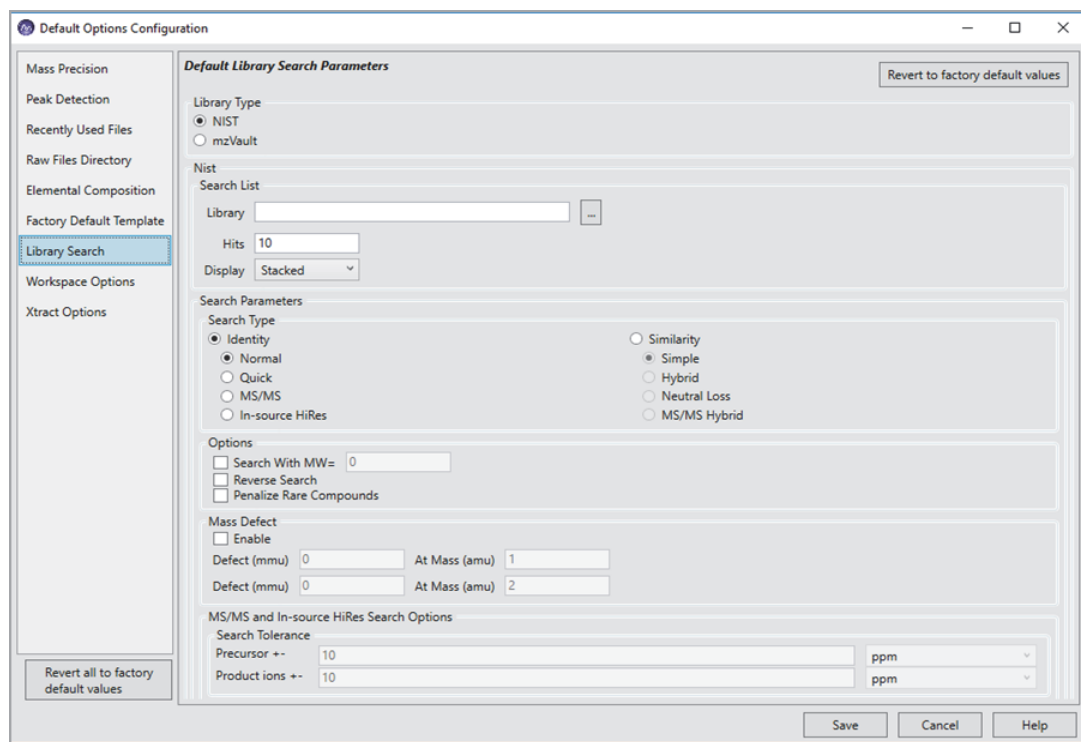
❖ To display the Default Library Search page

1. In the Workspace Options toolbar, click **Default Options**.

The Default Options Configuration dialog box opens.

[Figure 191](#) shows the factory default settings for the Library Search page.

Figure 191. Default Library Search page



2. In the left pane, click **Library Search**.

❖ To display the NIST library parameters

On the Default Library Search page, select the **NIST** option in the Library Type area.

❖ To display the *mzVault* library parameters

1. On the Default Library Search page, select the ***mzVault*** option in the Library Type area.

2. Make sure that the Library list in the Search List area includes the mzVault databases that you want to search (see [To specify the location of your local mzVault database files](#)).

Default Workspace Options Page

Use the Default Workspace Options page to specify the following:

- The minimum trace height (in centimeters) for the [Chromatogram View](#), [Spectrum View](#), or [MultiSpectrum View](#)
- The maximum number of scan filters to display in the Chromatogram view when applying the auto filter command (see [Adding Chromatogram Traces with the Auto Filter Feature](#))

Note When you adjust the height of these views, if the height becomes smaller than the set minimum value, a scrollbar automatically appears to the right of the views.

- How to open the Chromatogram view and Spectrum view toolboxes

❖ To display the Default Workspace Options page

1. In the Workspace Options toolbar, click **Default Options**.

The Default Options Configuration dialog box opens.

2. In the left pane, click **Workspace Options**.

[Figure 192](#) shows the factory default settings for the Workspace Options page.

Figure 192. Default Workspace Options page

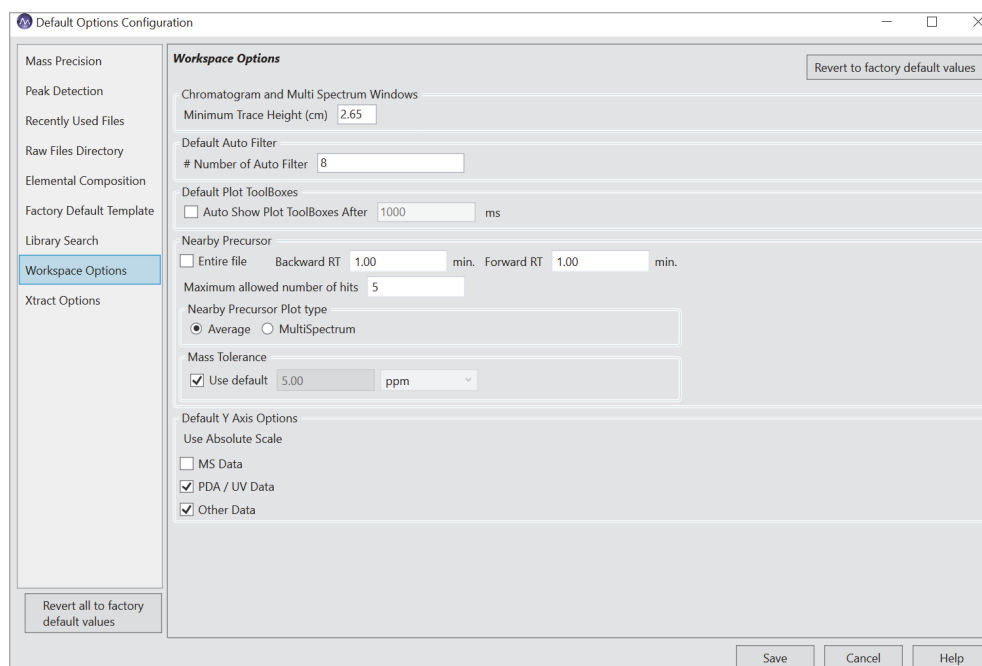


Table 63 describes the parameters on the Workspace Options page of the Default Options Configuration dialog box.

Table 63. Workspace Options page parameters (Sheet 1 of 3)

Parameter	Description
Chromatogram and Multi Spectrum Windows	
Minimum Trace Height (cm)	Specifies the minimum height, in centimeters, for the traces in the Chromatogram and Multi Spectrum views. Default: 2.65; range: 0.5 to 15.0
Default Auto Filter	
# Number of Auto Filter	Specifies the maximum number of filtered chromatograms that the application displays in the Chromatogram Ranges and Chromatogram views. Default: 8; range: 1 to 500
Default Plot Toolboxes	
Auto Show Plot Toolboxes After	<ul style="list-style-type: none"> • Select to turn on the automatic display of the view's toolbox with these mouse pointer actions: <ul style="list-style-type: none"> – Point to the Chromatogram view for the specified time, in milliseconds. – Point to an <i>m/z</i> value in the Spectrum view for the specified time, in milliseconds. • Clear to turn off the automatic display of the view's toolbox. <p>For more information, see Selecting How You Open the Floating Toolboxes.</p> <p>Tip By default, there are two ways to open the toolbox:</p> <ul style="list-style-type: none"> • Right-click the view and choose Show Toolbox. <p>–or–</p> <ul style="list-style-type: none"> • Use the mouse pointer. <p>Clearing this check box turns off the mouse pointer action.</p>
Time box (ms)	Available when the Auto Show Plot Toolboxes After check box is selected. Default: 1000 ms; range: 100 to 2000 ms

Table 63. Workspace Options page parameters (Sheet 2 of 3)

Parameter	Description
Nearby Precursor	
Use this feature to label the spectrum peaks with related data-dependent scans within the specified time range of the current scan. A related data-dependent scan is an MS _n +1 stage data-dependent scan for a precursor ion that matches the <i>m/z</i> value of the current peak within the specified mass tolerance. See Using the Nearby Precursors Marker .	
Entire File	When selected, sets the search to every scan in the file (no time range limit). Default: Clear
Backward RT	Specifies the first time point for the search: $RT_{\text{current scan}} - \text{Specified time in minutes}$ Default: 1.00 min; range: 0.00 to 10.00 min Not available when the Entire File check box is selected.
Forward RT	Specifies the last time point for the search: $RT_{\text{current scan}} - \text{Specified time in minutes}$ Default: 1.00 min; range: 0.00 to 10.00 min Not available when the Entire File check box is selected.
Maximum Allowed Number of Hits	Specifies the maximum number of matching data-dependent scans to display or average. If the application finds more than this maximum number, it displays the data-dependent scans with the nearest RT to the precursor “survey” scan. Default: 5; range: 1 to 50
Nearby Precursor Plot Type	Specifies whether double-clicking a marker in the precursor “survey” scan opens an average spectrum of the matching data-dependent scans in another Spectrum view or opens a MultiSpectrum view with a separate plot for each matching data-dependent scan. Options: Average (default) or MultiSpectrum

Table 63. Workspace Options page parameters (Sheet 3 of 3)

Parameter	Description
Mass Tolerance	<p>Specifies how close the m/z value of the precursor ion for the data-dependent scan must be to the m/z value of the current spectrum peak.</p> <p>Default: Use Default is selected—the application uses the mass tolerance of the mass spectrometer.</p> <p>Use Default check box is clear—the settable range is from 0.00 to 1000.00 (amu, mmu, or ppm).</p>
Default Y Axis Options	
<p>Use this feature to set the Y-Scale units as Relative or Absolute for each detector type. Change the Y Axis units of individual traces of the Chromatogram view by enabling the Separate Axis button in the Y Axis area in Display Options toolbar.</p>	
MS Data	<p>The default Y-Scale unit for MS Data is Relative.</p> <p>Select the check box for MS Data to change the default setting to Absolute scale.</p>
PDA/UV Data	<p>By default, the Y-Scale unit for PDA and UV Data is set to Absolute.</p> <p>Clear the default check box to view the PDA and UV data in Relative scale.</p>
Other Data	<p>By default, the Y-Scale unit for Other Data is set to Absolute.</p> <p>Clear the default check box to view the Other data in Relative scale.</p>

Default Xtract Options Page

Use the Default Xtract Options page to specify the default Xtract parameters. For more information, see [Chapter 13, “Deconvolving and Deisotoping Spectra with the Xtract Algorithm.”](#)

❖ To display the Default Xtract Options page

- In the Workspace Options toolbar, click **Default Options**.
The Default Options Configuration dialog box opens.
- In the left pane, click **Xtract Options**.

[Figure 193](#) shows the Default Xtract Options page with its original default values.

Figure 193. Default Xtract Options page

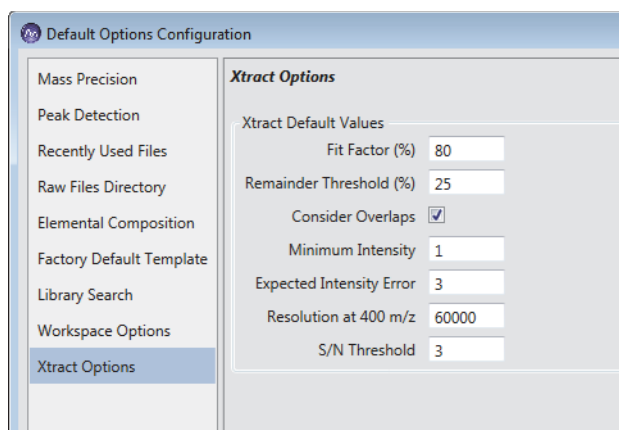


Table 64 lists the default Xtract parameters.

Table 64. Default parameters for the Xtract algorithm (Sheet 1 of 3)

Parameter	Description
Fit Factor (%)	<p>Measures the quality of the match between a measured isotope pattern and an average distribution of the same mass, as a percentage.</p> <ul style="list-style-type: none"> 0% requires a low fit only. 100% means that the measured isotope profile is identical to the theoretical average isotope distribution. <p>Range: 0–100; default: 80</p>
Remainder Threshold (%)	<p>Specifies the height of the smaller overlapping isotopic cluster (as a percentage) with respect to the height of the most abundant isotopic cluster when the Xtract algorithm attempts to resolve overlapping isotopic clusters.</p> <p>For example, if one isotopic cluster in a spectrum has an abundance of 100 and you set the Remainder Threshold parameter to 30 percent, the Xtract algorithm ignores any overlapping clusters with an abundance less than 30.</p> <p>Range: 0–100; default: 25</p>

Table 64. Default parameters for the Xtract algorithm (Sheet 2 of 3)

Parameter	Description
Consider Overlaps	<p>Determines whether the Xtract algorithm is more tolerant of errors when the spectrum intensity is significantly higher than expected for the theoretical isotopic cluster.</p> <ul style="list-style-type: none"> Selected: The Xtract algorithm is more tolerant of errors when the spectrum intensity is significantly higher than expected for the theoretical isotopic cluster. Because this option can lead to increased false positives, select it only in cases where you expect overlapping isotopic clusters in a data set. Cleared: The Xtract algorithm is less tolerant of errors when the spectrum intensity is significantly higher than expected for the theoretical isotopic cluster. <p>Default: Selected</p>
Minimum Intensity	<p>Specifies a minimum intensity threshold to filter out possible background noise, even when you set the S/N Threshold parameter to zero.</p> <p>Range: 0–9999; default: 1</p>
Expected Intensity Error	<p>Specifies the permissible percentage of error allowed in calculating the ratio of the most abundant isotope to the next isotope that is higher in mass in the isotope series.</p> <p>Range: 0–9999; default: 3</p>
Resolution at 400 m/z	<p>Defines the resolution of the source spectrum at m/z 400.</p> <p>This parameter is not needed if the Xtract algorithm deconvolves FTMS, Orbitrap, or Exactive data because the data contains the instrument information in the spectrum. You must set this parameter for all other spectrum types and for exported spectrum files (in -qb.raw data file format), which lack instrument information.</p> <p>Range: 6000–240 000; default: 60 000</p>

Table 64. Default parameters for the Xtract algorithm (Sheet 3 of 3)

Parameter	Description
S/N Threshold	<p data-bbox="781 338 1450 474">Specifies a signal-to-noise (S/N) threshold, x, above which the Xtract algorithm considers a measured peak to be a real (accepted) peak. The Xtract algorithm ignores peaks below this threshold.</p> <p data-bbox="781 506 1386 611">Any spectral peak must be x times the intensity of the calculated noise for that spectrum before the Xtract algorithm considers it.</p> <p data-bbox="781 642 1078 674">Range: 0–9999; default: 3</p>

Scan Filters and Scan Headers

The scan method used for each scan is recorded as a scan event. The data system creates scan filters from the scan-event settings. You can select or create scan filters to apply processing to a subset of the scans in the raw data file. Scan headers provide important information about the scan.

Contents

- [Scan Filter Parameters](#)
- [Scan Headers and Scan Header Abbreviations](#)

Scan Filter Parameters

Table 65 lists the scan filter parameters. Use only the fields that apply to your mass spectrometer. You can define additional scan filters by adhering to the following scan filter format.

Note Not all features are applicable for every mass spectrometer.

Table 65. Scan filter format (Sheet 1 of 4)

Feature	Option	Interpretation
Polarity	+, –	Positive or negative.
Data type	p, c	Profile or centroid.
Dependent scans	d, !d	Include dependent scans or exclude dependent scans.
TurboScans	t, !t	Include TurboScan scans or exclude TurboScan scans.
Source CID	cid, !cid	Include Source CID scans or exclude Source CID scans. (Source CID scans are scans for ions that are produced by collision-induced dissociation in the ion source.)
Scan type	FULL, Z, SIM, SRM, CRM, Q1MS, Q3MS	Full scan, ZoomScan, selected ion monitoring (SIM), selected reaction monitoring (SRM), consecutive reaction monitoring (CRM), Q1 quadrupole analysis, or Q3 quadrupole analysis.

Table 65. Scan filter format (Sheet 2 of 4)

Feature	Option	Interpretation
Scan mode	ms, ms2, ms3, ... MS10	MS ⁿ where $n = 1$ to 10 Each order can be followed by the appropriate number of parents. The parents can also be omitted. Example: “ms3 345.3, 253.2” indicates an MS ³ scan with parents with m/z 345.3 and 253.2.
	pr	Parent (followed by the product mass)
	cng	Constant neutral gain (followed by the mass of the neutral)
	cnl	Constant neutral loss (followed by the mass of the neutral)
Mass Analyzer	ITMS, TQMS, SQMS, TOFMS, FTMS, Sector	Ion trap, triple-quadrupole, single-quadrupole, time-of-flight, Fourier transform, or magnetic sector mass spectrometry.
Photo Ionization	pi, !pi	Include photo ionization scans or exclude photo ionization scans.
Compensation Voltage	cv, !cv	Include compensation voltage scans or exclude compensation voltage scans.
Detector Valid	det, !det	Include detector valid scans or exclude detector valid scans.
Enhanced	E, !E	Include enhanced scans or exclude enhanced scans.
Wideband	w, !w	Include wideband scans or exclude wideband scans.
Supplemental Activation	sa, !sa	Include supplemental activation scans or exclude supplemental activation scans.
Multistate Activation	msa, !msa	Include multistate activation scans or exclude multistate activation scans.
Product masses or mass range of scan	[m1a–m1b, m2a–m2b, m3a–m3b, ...]	Scans with a specific mass range or mass ranges, such as SIM, SRM, and CRM. Example: [50.00–1500.00] for a scan from m/z 50.00 to 1500.00
		If a scan is exactly 1 Da wide, it appears as a single value (the center mass). This is typical for SIM, SRM, and CRM. Filters for precursor ions in dependent scans are matched with a tolerance of m/z 1.0 so that minor differences in precursor mass measurements from scan to scan do not give different filters.
Segment/scan event number pairs	{segment, scan number}	Example: “{3, 4} + c ms” indicates segment 3, scan event 4 for a positive centroid MS scan The curly brackets { } are required.

Table 65. Scan filter format (Sheet 3 of 4)

Feature	Option	Interpretation
Ionization mode	APCI, ESI, EI, CI, NSI, FAB, TSP, FD, MALDI, GD, PSI	Atmospheric pressure chemical ionization (APCI), electrospray (ESI), electron ionization (EI), chemical ionization (CI), nanoelectrospray ionization (NSI), fast atom bombardment (FAB), thermospray ionization (TSP), field desorption (FD), matrix-assisted laser desorption ionization (MALDI), or glow discharge (GD), Paper Spray Ionization (PSI). Example: “+ c ESI ms” indicates a positive centroid electrospray MS scan.
Corona on/off	corona, !corona	Corona on or corona off. Example: “+ APCI !corona ms” indicates a positive centroid APCI scan with the corona off.
Detector value	“det=## .##”	Detector value is ## .## with no spaces. Example: “+ ESI det= -800.0” indicates a positive electrospray scan at -800.0 detector units (usually volts).
MS/MS and MS ⁿ CID energies	mass@energy	Mass is the precursor mass and energy is the CID relative energy (no units). Example: “- c ms2 196.1@25.0” indicates a negative centroid MS/MS scan of <i>m/z</i> 196.1 at 25.0 units of CID energy.
Quadrupole identification	Q1MS, Q3MS	Q1 quadrupole or Q3 quadrupole. Example: “+ c ESI Q3MS” indicates a positive centroid electrospray MS scan using Q3 quadrupole.
Accurate mass	AM, !AM, AMI, AME	Include accurate mass scans, exclude accurate mass scans, include accurate mass internal, or include accurate mass external.
Ultra	u, !u	Include ultra scans or exclude ultra scans.
Sector	BSCAN, !BSCAN	Include magnetic sector scans or exclude magnetic sector scans.
	ESCAN, !ESCAN	Include electric sector scans or exclude electric sector scans.
LOCK	lock, !lock	Include lock scans or exclude lock scans.
Multiplex	msx, !msx	Include multiplexing scans or exclude multiplexing scans.
Electron capture dissociation	ecd, !ecd	Include electron capture dissociation or exclude electron capture dissociation.
Multi-photon dissociation	mpd, !mpd	Include photo dissociation or exclude photo dissociation.

B Scan Filters and Scan Headers

Scan Headers and Scan Header Abbreviations

Table 65. Scan filter format (Sheet 4 of 4)

Feature	Option	Interpretation
Pulsed dissociation	pqd, !pqd	Include pulsed dissociation scans or exclude pulsed dissociation scans.
Electron transfer dissociation	etd, !etd	Include electron transfer dissociation scans or exclude electron transfer dissociation scans.
NPTR	nptr, !nptr	Include NPTR, Exclude NPTR
Higher-energy CID	hcd, !hcd	Include higher energy CID scans or exclude higher energy CID scans.
Source SID	sid, !sid	Include source SID scans or exclude Source SID scans. (Source SID scans are surface-induced scans.)
Isolation width	iw ##, ##	Isolation width value at ##, ##

Scan Headers and Scan Header Abbreviations

Use the Scan Header dialog box to select what parameters appear in the scan header of the [Spectrum View](#) or [MultiSpectrum View](#).

❖ To display the Scan Header dialog box

1. Click the [Spectrum View](#) or [MultiSpectrum View](#) to select it.
2. Click the **Display Options** toolbar tab.
3. Click **Scan Header**.

[Figure 194](#) shows the default scan header selections.

Figure 194. Default scan header selections

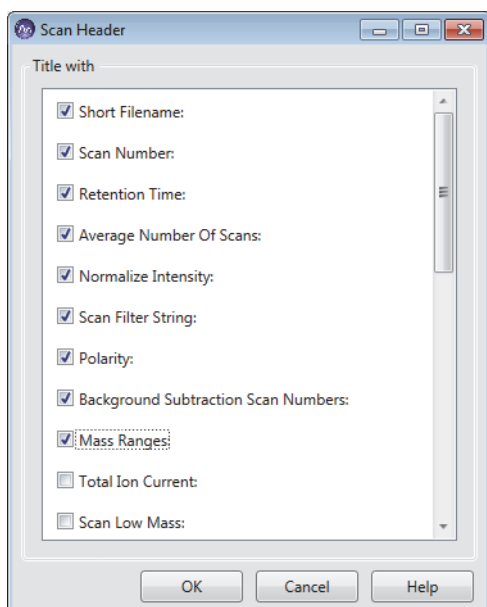


Table 66 lists the parameters that you can display in the scan header. The parameters displayed in the Scan Header dialog box depend on the data acquisition settings for the mass spectrometer used to acquire the raw data file.

Table 66. Scan header parameters and their abbreviations (Sheet 1 of 2)

Parameter	Abbreviation before the value
Short filename	NA
Scan Number	#
Retention time	RT
Average Number of Scans	AV
Normalize Intensity	NL
Scan Filter String	T:
Polarity	NA (P: when the Scan Filter String is not selected)
Background Subtraction Scan Numbers	SB
Mass Ranges	[...] (MR: [...] when the Scan Filter String is not selected)
Total Ion Current	TIC:
Scan Low Mass	SLM:

Table 66. Scan header parameters and their abbreviations (Sheet 2 of 2)

Parameter	Abbreviation before the value
Scan High Mass	SHM:
Scan Start Time (Min)	SST (Min):
Base Peak Intensity	BPI:
Base Peak Mass	BPM:
Multiple Injection	Multiple Injection::
Multi Inject Info	Multi Inject Info::
AGC (automatic gain control—On, Off, or predicted)	AGC::
Micro Scan Count	MSC:
Ion Injection Time (ms)	IIT
Scan segment	SS:
Scan Event	SE:
Master Index	Master Index::
Charge State	CS:
Reagent Ion AGC	Full parameter name
Reagent Ion Injection Time (msec)	Full parameter name
Elapsed Scan Time (sec)	EST
API Source CID Energy	Full parameter name
Average Scan By Instrument	ABSI
Charge State	CS
Monoisotopic Ion M/Z	Full parameter name
MS ⁿ Isolation Width ($n = 2$ to 10) (m/z)	Full parameter name
FT analyzer settings	Full parameter names
FT analyzer message	Text message
FT resolution	Full parameter names
Conversion parameters	Full parameter names

One- and Three-Letter Abbreviations for Amino Acid Residues

Table 67 lists common one- and three-letter abbreviations for amino acid residues that you enter in the peptide/protein Formula box on the [Isotope Simulation Page](#).

Table 67. Common one- and three-letter abbreviations for amino acid residues (Sheet 1 of 2)

One letter	Name	Formula	Three letter
A	Alanine	C ₃ H ₅ NO	Ala
C	Cysteine	C ₃ H ₅ NOS	Cys
D	Aspartate	C ₄ H ₅ NO ₃	Asp
E	Glutamate	C ₅ H ₇ NO ₃	Glu
F	Phenylalanine	C ₉ H ₉ NO	Phe
G	Glycine	C ₂ H ₃ NO	Gly
H	Histidine	C ₆ H ₇ N ₃ O	His
I	Isoleucine	C ₆ H ₁₁ NO	Ile
K	Lysine	C ₆ H ₁₂ N ₂ O	Lys
L	Leucine	C ₆ H ₁₁ NO	Leu
M	Methionine	C ₅ H ₉ NOS	Met
N	Asparagine	C ₄ H ₆ N ₂ O ₂	Asn
O	Ornithine	C ₅ H ₁₁ N ₂ O	Orn
P	Proline	C ₅ H ₇ NO	Pro
Q	Glutamine	C ₅ H ₈ N ₂ O ₂	Gln
R	Arginine	C ₆ H ₁₂ N ₄ O	Arg
S	Serine	C ₃ H ₅ NO ₂	Ser

Table 67. Common one- and three-letter abbreviations for amino acid residues (Sheet 2 of 2)

One letter	Name	Formula	Three letter
T	Threonine	C ₄ H ₇ NO ₂	Thr
V	Valine	C ₅ H ₉ NO	Val
W	Tryptophan	C ₁₁ H ₁₀ N ₂ O	Trp
Y	Tyrosine	C ₉ H ₉ NO ₂	Tyr

Table 68 lists less common three-letter abbreviations for amino acid residues to enter in the peptide/protein formula for the simulated spectrum.

Table 68. Less common three-letter abbreviations for amino acid residues (Sheet 1 of 3)

Three letter	Name	Formula
Abu	2-Aminobutyric acid (2-aminobutanoic acid)	C ₄ H ₇ NO
Aec	Aminoethylcysteine	C ₅ H ₁₀ N ₂ OS
Aib	Aminoisobutyric acid	C ₄ H ₇ NO
Aln	--	C ₁₃ H ₁₁ NO
Aly	Alveolysin	C ₁₂ H ₂₂ N ₂ O ₆
Amc	--	C ₆ H ₁₀ N ₂ O ₂ S
Bcy	--	C ₁₀ H ₁₁ NOS
Bgl	--	C ₁₂ H ₁₃ NO ₃
Bly	--	C ₁₆ H ₂₆ N ₄ O ₃ S
Bse	--	C ₁₀ H ₁₁ NO ₂
Bth	--	C ₁₁ H ₁₃ NO ₂
Cmc	Carboxymethylcysteine	C ₅ H ₇ NO ₃ S
Cml	--	C ₈ H ₁₄ N ₂ OS
Cph	Chlorophenylalanine	C ₉ H ₈ NOCl
Cya	Cysteic acid	C ₃ H ₅ NO ₄ S
Dha	Dehydroalanine	C ₃ H ₃ NO
Dhb	Dehydro-2-aminobutyric acid	C ₄ H ₅ NO
Dpr	D-proline	C ₅ H ₅ NO
Dty	Diiodotyrosine	C ₉ H ₇ NO ₂ I ₂
Fcy	--	C ₁₈ N ₂₉ NOS
Fph	--	C ₉ H ₈ NOF
Ftr	--	C ₁₂ H ₁₀ N ₂ O ₂

Table 68. Less common three-letter abbreviations for amino acid residues (Sheet 2 of 3)

Three letter	Name	Formula
Gaa	--	C ₄ H ₇ NO
Gcg	--	C ₅ H ₅ NO ₄
Gla	Carboxyglutamic acid	C ₆ H ₇ NO ₅
Glp	--	C ₅ H ₅ NO ₂
Hse	Homoserine	C ₄ H ₇ NO ₂
Hsl	Homoserine lactone	C ₄ H ₅ NO
Hya	Beta-hydroxyaspartate	C ₄ H ₅ NO ₄
Hyg	Hydroxyglycine	C ₅ H ₇ NO ₄
Hyl	Hydroxylysine	C ₆ H ₁₂ N ₂ O ₂
Hyp	Hydroxyproline	C ₅ H ₇ NO ₂
Ils	Isolysine	C ₉ H ₁₈ N ₂ O
Ity	Iodotyrosine	C ₉ H ₈ NO ₂ I
Iva	Isovaline	C ₅ H ₉ NO
Mar	--	C ₇ H ₁₄ N ₄ O
Mas	--	C ₅ H ₇ NO ₃
Mbt	--	C ₁₇ H ₁₇ NO ₂
Mes	--	C ₅ H ₉ NO ₃ S
Mga	--	C ₆ H ₁₀ N ₂ O ₂
Mgl	--	C ₆ H ₉ NO ₃
Mhi	--	C ₇ H ₉ N ₃ O
Mls	--	C ₇ H ₁₄ NO
Mme	--	C ₆ H ₁₁ NOS
Mph	--	C ₁₀ H ₁₁ NO
Mso	Methioninesulfoxide	C ₅ H ₉ NO ₂ S
Mty	C ₁₀ H ₁₁ NO ₂	
Nle	Norleucine	C ₆ H ₁₁ NO
Nls	Norlysine	C ₁₂ H ₁₅ N ₃ O ₂
Pal	--	C ₈ H ₈ N ₂ O
Pcy	--	C ₁₉ H ₃₅ NO ₂ S
Pec	--	C ₁₀ H ₁₂ N ₂ OS
Pip	2-Piperidinecarboxylic acid	C ₆ H ₉ NO

Table 68. Less common three-letter abbreviations for amino acid residues (Sheet 3 of 3)

Three letter	Name	Formula
Psr	Phosphoserine	C ₃ H ₆ NO ₅ P
Pth	Phosphothreonine	C ₄ H ₈ NO ₅ P
Pty	Phosphotyrosine	C ₉ H ₁₀ NO ₅ P
Pyr	Pyroglutamic acid	C ₅ H ₅ NO ₂
Sar	Sarcosine	C ₃ H ₅ NO
Sas	--	C ₈ H ₈ NO ₅
Tml	E-amino trimethyl-lysine	C ₉ H ₁₉ N
Tys	Tyrosinesulfonic acid Tyr (SO ₃ H)	C ₉ H ₉ NO ₅ S

Common Isotopes

Table 69 lists the exact mass and natural abundance for some common isotopes.

Table 69. Common elements with multiple stable isotopes

Element	Isotope symbol	Mass (Da) ^a	%Abundance ^b
Bromine	⁷⁹ Br	78.918336	50.69
	⁸¹ Br	80.916290	49.31
Carbon	¹² C	12.000000	98.93
	¹³ C	13.003354	1.07
Calcium	⁴⁰ Ca	39.962591	96.95
	⁴² Ca	41.958622	0.65
	⁴⁴ Ca	43.955485	2.086
Chlorine	³⁵ Cl	34.968853	75.77
	³⁷ Cl	36.965903	24.23
Potassium	³⁹ K	38.963708	93.20
	⁴¹ K	40.961825	6.73
Nitrogen	¹⁴ N	14.003074	99.63
	¹⁵ N	15.00011	0.37
Oxygen	¹⁶ O	15.994915	99.76
	¹⁸ O	17.999159	0.20
Sulfur	³² S	31.972071	95.02
	³³ S	32.971459	0.75
	³⁴ S	33.967867	4.21

^a Mass values are for reference only.

^b Isotopes with a natural abundance below 0.20% are not listed. Percent abundance values are for reference only.



Index

Numerics

10x Absolute button 98, 170

A

active views 19

Add Peak button 100

adding

chromatogram traces 53

chromatographic peaks manually 88

custom text and graphic annotations 28

EIC plots 67

filtered chromatograms with the Auto Filter
button 334

labels, chromatogram trace 73

MultiSpectrum views 142

mzVault database files to the Library list 242

scan header information 164

spectrum peak labels 166

spectrum plots 142–143

Spectrum views 140

text labels 29

adduct ions

isotope simulation 278

Xtract algorithm 313

adduct species 271

adductions 5

All Samples, centroiding algorithm 280

amino acid residues

common one-letter abbreviations (table) 347

common three-letter abbreviations (table) 347

less common three-letter abbreviations (table) 348

analog-to-digital converter 118

annotations, adding 28

Apex Samples, centroiding algorithm 280

Auto Filter button, using 53

auto filter feature, default setting 334

Auto Show Plot Toolboxes After check box 37

Automatic setting 257

Avalon Peak Detection page 130

average spectrum

creating 147

displaying 161

MSn Browser page 158, 188

procedure 146

toolbox icon 115

averagine isotope distribution 320, 337

B

background subtraction, spectra

procedure 149

toolbox icon 115

Behind Graph icon, using 29

C

Calculated Monoisotopic m/z column 319

CAS ID column 249

centroiding algorithm 280

centroids for isotope simulation in Spectrum view 272

Charge Normalized Intensity column 319

Charge State column 319–320

Chemical Formula column, Chromatogram Ranges view 106

chemical formula labels 167

chemical formula type (isotope simulation) 277

chemical formulas used to filter library searches 259

chemical ionization 5

Chromatogram button, Workspace Options toolbar 52

Chromatogram Display Options toolbar 95

Chromatogram Ranges view

adding chromatogram traces 52

comparing chromatogram traces 93

Display column 55

entering a chemical formula 106

figure 104

fill down feature 108

parameters 105

- chromatogram traces
 - adding 53
 - base peak 51
 - Chromatogram Display Options toolbar 95
 - comparing 93
 - delay time 107
 - deleting 54
 - detector type, selecting 105
 - extracted ion 51
 - filtering 105, 114
 - formats 95
 - labeling 96
 - mass range 51
 - neutral fragment 51
 - Normalization area, Display Options toolbar 97
 - normalizing 76
 - plot operators 107
 - smoothing 106
 - total ion current (TIC) 51
 - trace type, selecting 105
 - y-axis scaling 98
 - Chromatogram view
 - comparing chromatogram traces 93
 - description 102
 - displaying 52
 - figure 102
 - pointer actions 27
 - toolbox 114–115
 - Chromatogram View button, Workspace Options toolbar 40
 - Clear Manual Noise Range button 100
 - Clipboard, copying images to 34
 - commands
 - Create Sequence 13
 - Report an Issue 14
 - Save As (template) 13
 - Save as Default (template) 13
 - communicator bar 14
 - composite spectra, viewing 162
 - Compound column 248
 - Compound ID column 248
 - compound name, displayed 86
 - computer requirements x
 - concentration effect on adduct formation 271
 - Consider Overlaps parameter 338
 - contacting us xii
 - context menus
 - Chromatogram view 103, 182
 - Elements in Use Table 240
 - MSn Browser page 188
 - Copy to Clipboard command 34
 - Copy to Clipboard/Export dialog box 35
 - copying
 - graphical views 35
 - report views 34
 - Create Sequence button, Workspace Options toolbar 53
 - Create Sequence command, File menu 13
 - creating a library entry with the spectral data 265
 - CSV files, copying data to and from 59
 - Current Trace button 101
 - CV button 285
 - CV Plot view 282
 - CV Traces dialog box 281
- ## D
- data-dependent scans
 - Nearby Precursor marker 155
 - precursor markers for 152
 - viewing 152
 - decimal places, numeric chromatogram labels 74
 - Deconvoluted Spectrum view 185
 - deconvolution results 318
 - Deconvoluted Spectrum view 317
 - default data visualization application, choosing 36
 - defaults, setting
 - elemental composition 329
 - library search parameters 332
 - mass precision 325
 - minimum trace height 334
 - number of recently used files 327
 - peak detection algorithm 327
 - raw data file directory 328
 - workspace options 333
 - Xtract parameters 336
 - deisotoping 309
 - Delete Peak button 100
 - deleting
 - chromatogram traces 54, 102
 - elements from the Elements in Use table 228
 - events, Avalon Peak Detection 133
 - libraries 264
 - peaks from the Peaks List 90
 - rows, Chromatogram Ranges view 110
 - text annotations 45
 - Delta Mass button, using 168
 - Delta Mass column 249
 - Detect in Active Plot button 100
 - Detect in All Plots button 100
 - Detector Type page 117
 - documentation
 - accessing x
 - additional x
 - double bonds 236
 - drawing lines and boxes 28

E

- Edit_text 30
- elemental composition
 - default settings, specifying 329
 - toolbox icon 185
- Elemental Composition button, using 167
- Elemental Composition check box, Peaks List view 112
- Elemental Composition page
 - description 234
 - parameters 235
- Elemental Composition Results view
 - description 231
 - parameters 231
- Elements in Use table 227
- EMF files, saving images as 34
- Error Log view 307
- Exactive mass spectrometers, resolution information 338
- Expected Intensity Error parameter 338
- experiment information, acquisition-specific 295
- exported spectrum files (in -qb raw data file format) 338
- exporting
 - graphical views as EMF images 35
 - tabular data to a CSV file 35
- extracted ion current (EIC) traces, displaying 67

F

- F1 key, using to access context-sensitive Help x
- factory default layout 17
- factory default layout, applying 325
- factory default template settings 330
- File Header view 304–305
- files, default number of recently used files 327
- Fill Down dialog box, Chromatogram Ranges view 108
- Filter column 249
- Fit % column 319
- Fit Factor (%) parameter 337
- floating toolboxes 37
- floating views 21
- Formula column 248
- formula labels, adding 167
- formula types
 - chemical (isotope simulation) 277
 - peptide (isotope simulation) 278
- Fragment Tolerance parameter 257
- fragmentation patterns 4
- FreeStyle application
 - default program for RAW files 36
 - default settings 323
 - FreeStyle window (figure) 18
 - Info Bar pages 47

- spectra-specific views 180
- toolbars 38
- Workspace Options toolbar 38
- FTMS mass spectrometers 338

G

- Genesis Peak Detection page 122–123
- global normalization, chromatograms 77
- global normalization, spectrum plots 177
- green, vertical lines 272

H

- half-width of charge distribution 271
- height, trace 334
- Help, accessing x
- hiding views 21
- High Res Score column 249
- hover box
 - Chromatogram view 114
 - Spectrum view 185
- hydrogen 313

I

- ICIS Peak Detection page 127–128
- images of views, copied to Clipboard 34
- Info Bar
 - location (figure) 18
 - pages, hiding 48
 - pages, list 47
 - pages, rearranging 48
- Info Bar pages
 - Avalon Peak Detection page 130
 - Detector Type page 117
 - Elemental Composition page 234
 - Genesis Peak Detection page 122
 - ICIS Peak Detection page 127
 - Isotope Simulation page 277
 - mzVault Search page 255
 - NIST Search page 250
 - PPD Peak Detection page 133
 - Scan Filters page 119
 - Sequence page 136
 - Trace Type page 118
- information, acquisition-specific 295
- information, general 14
- Insert Chromatogram command 53
- instrument
 - information 295
 - readbacks, displaying 301
- Instrument Method view

Index: L

- Autosampler page (figure) 301
 - description 301
 - displaying 301
 - LC page (figure) 302
 - Mass Spectrometer page (figure) 302
- integration markers 89
- integration method, manual 89
- ionization modes, effect 4
- isotope intensity threshold 238
- Isotope Simulation page
 - figure 270
 - parameter descriptions 277
- isotope simulations
 - chemical formula or peptide 277
 - Isotope Simulation window 273
 - labeled compounds 270
- isotopic
 - clusters 318, 337
 - patterns for proteins 274
- isotopically resolved mass spectra 309

L

- Label Threshold box
 - chromatograms 75
 - spectra 166
- labeled compounds, isotope simulations 270
- labeling
 - intensity threshold, chromatograms 75
 - x axis, Chromatogram view 99
 - y axis, Chromatogram view 98
- labels
 - compound name 86
 - formula 167
 - integrated chromatograms 74
 - spectrum peaks 166
- layouts
 - applying 41
 - factory default 17
 - Factory Default Template page 330
 - saving 41
- libraries
 - adding 264–265
 - copying, from NIST libraries list 265
 - creating entries in 264–265
 - searching NIST or mzVault 241
 - searching, default settings 332
- linking a spectrum to a chromatogram 146
- local normalization, chromatograms 76, 97
- local normalization, spectrum plots 177

M

- manual integration method 89
- Manual Noise Region check box 87
- Manual option, fragment tolerance 257
- mass chromatograms 7
- mass peaks, labeling 166
- mass precision, defaults 325
- mass tolerance
 - Chromatogram Ranges view 58
 - Default Mass Precision page 325
 - Elemental Composition page 235
 - File Header view 307
 - mzVault Search page 260
- Mass Tolerance column, Chromatogram Ranges view 107
- mass-to-charge labels, Spectrum view 68
- maximum number, autofiltered chromatograms 53
- Microsoft Office applications, copying images for 34
- Minimum Intensity parameter 338
- Minimum Trace Height parameter 334
- monoisotopic mass 310
- Monoisotopic Mass for This Charge column 319–320
- Monoisotopic Mass parameter 318
- Mostabund m/z column 319–320
- mouse pointer actions 27
- MS Trending detector type 67
- MSn Browser page 158
- MSn experimental data, analyzing and displaying 186
- Multi Spectrum view 183
- mzVault option 332
- mzVault Search page 255, 258

N

- Name column, Peaks List 113
- Nearby Precursors marker 155
- neutral loss 3
- Neutral Mass column 249
- New Trace button 53, 101
- New Workspace icon 16
- NIST Mass Spectral Search application, purchasing 5
- NIST Search page 250–251
- NIST Search Results view 245–246
- Nitrogen Rule 237
- noise region, selecting 87
- normalization, chromatograms
 - global 77
 - local 76
 - off 78
 - options 97
- normalization, spectrum plot 177

nucleotides, deconvolving spectra with the Xtract algorithm 309

O

One Page option, printing 36

Orbitrap resolution of source spectrum 338

P

PDF file, creating 36

peak area label 74

peak detection

algorithm default settings 327

Avalon 130

Genesis 122

ICIS 127

manual 88

PPD 133

peak height label 74

Peak List button 100

Peaks List view 110, 112

Peptide Fragments page 289

peptides

deconvolving spectra with the Xtract algorithm 309

formula type (isotope simulation) 278

pointers

Add Peak pointer 88

Average Spectrum pointer 147

Background Subtraction pointers 149

Delete Peak pointer 90

zooming in on a chromatogram 83

zooming in on a spectrum 172

potassium 313

PowerPoint presentations, copying images for 34

PPD Peak Detection page 133–134

Precursor m/z column 249

precursor markers 152

presentations, copying images for 34

printing

contents of a view or workspace 35

MSn tree 189

procedures

background spectra, subtracting 149

chromatogram peaks

adding 88

deleting 90

integrating 84

chromatogram traces

adding 53

comparing 93

normalization 76

composite spectra, viewing 162

detecting and integrating chromatographic peaks 84

instrument readbacks, displaying 301

sequences, creating 92

simulating an isotope pattern 269

spectra

averaging 146

displaying 140

templates

applying a saved template 27

applying default 26

restoring factory default 14, 27

saving as default 13, 26

saving as XML file 13, 26

views, arranging 23

Profile check box 276

proteins

deconvolving spectra with the Xtract algorithm 309

isotope patterns 274

publications, copying images for 34

purple circle spectrum peak marker 156

R

raw data files

Chromatogram Ranges view 105

default directory, selecting 328

displaying name and URL 249

opening 15

Peaks List view 113

Raw File URL column 249

readback parameters 67

reagent gas 5

recently used files 327

Released description 146

releasing a spectrum 55

Remainder Threshold (%) parameter 337

Remove Precursor check box 257

Remove Precursor Ion parameter, mzVault library search 261

repeating moieties, isotope simulations 277

Replace All Traces button 101

Report an Issue command 14

requirements, system x

Reset icon, Peak List view 112

Reset Scaling command 83, 172

Resolution at 400 m/z parameter 338

resolution, effect of 272

Results table, Xtract

columns in 319

results of deconvolution in 318

Retention Time column 249

retention time label, formatting 74

retention time, mzVault spectrum information 249

right-click menus
 Chromatogram view 103, 182
 Elements in Use table 240
 MSn Browser page 188
 ring equivalents 236

S

S/N Threshold parameter 339
 Sample Information view
 description 304
 displaying 304
 figure 304
 Save As command 13
 Save as Default command 13
 saving template 13, 26
 scan filter 249
 Scan Filter page 119
 Scan Filters button 40, 71
 scan filters, formats (table) 341
 scan header information 182
 Scan Header view 297–298
 scan headers
 abbreviations 344
 parameters 298
 Scan Header dialog box 178
 scan number 249
 Scan Number column 249
 Score Cutoff parameter 257
 Search Results view 245
 seconds, RT label 74
 Select Manual Noise Range button 100
 Select Trace to Link 145
 selected ion monitoring 7
 selected views 19
 Selection As button 35
 Separate Page option 36
 Sequence File tab 53
 Sequence page 136–137
 Sequence toolbar 100–101
 sequences
 comparing chromatogram traces 93
 creating 92
 shortcut menus
 Chromatogram view 103, 182
 Elements in Use table 240
 MSn Browser page 188
 Show Extracted Ion Chromatogram check box 280
 Show Toolbox command 37, 103, 114, 185, 334
 Show within the Workspace check box 280
 signal-to-noise ratio label, chromatograms 74

signal-to-noise threshold 339
 simulated isotopic distribution spectrum,
 creating 173, 277
 sodium 313
 software requirements x
 spectra
 averaging
 command 115
 procedure 146
 background subtraction toolbox icon 115
 color 175
 data-dependent scans, viewing 152
 deconvolving with Xtract 309
 deisotoping with Xtract 309
 display options 175, 179
 displaying 140
 elemental composition icon 185
 isotopically resolved (figure) 310
 isotopically unresolved (figure) 310
 labels 175
 normalization 177
 Spectrum Display Options toolbar 174
 subtracting background 149
 working with 139
 Spectrum Display Options toolbar 174–175, 179
 Spectrum ID column 249
 Spectrum List Display Options toolbar 297
 Spectrum List view 298–299
 Spectrum Ranges dialog box 141
 Spectrum view
 about 180
 highlighting most abundant isotope 321
 pointer actions in Chromatogram view 27
 scan header information 182
 toolbox 185
 spreadsheet files 59
 Status Log view 300
 status readbacks 67
 system requirements x

T

templates
 applying a saved template 27
 applying the default template 26
 description 26
 restoring factory default template 14, 26–27
 saving as default template 13, 26
 saving as XML file 13, 26
 text labels, adding 29
 Theo. button, using 168
 tilde description 19

time range marker, spectrum averaging 148
 timebase, updating 27
 To Clipboard in EMF option 35
 To CSV File option 35
 tool, view arranger 23
 toolbars
 list 38
 location (figure) 18
 Sequence toolbar 100
 spectra-specific toolbars 172
 Spectrum Display Options toolbar 174
 Spectrum List Display Options toolbar 297
 Workspace Options toolbar 38, 295
 Workspace Processing toolbar 99, 173, 266
 Zoom Options toolbar 43
 toolboxes
 Chromatogram view 114
 Spectrum view 185
 trace analysis 8
 Trace Title icon 99
 Trace Type page 118–119
 triangular markers 152
 Tune Method view 303

U

undo
 background subtraction 149
 last deleted spectrum, MultiSpectrum view 182
 manual noise region 88
 manual peak 89
 spectrum averaging 147
 Undo Delete Chromatogram command 103
 Undo Delete Spectrum command, MultiSpectrum view 183
 units, y axis of the Chromatogram view 82
 unlinking a spectrum 55
 Use Fragments Matching check box 238

V

view arranger tool 23
 views
 arranging 23
 Chromatogram Ranges view 104
 Chromatogram view 102
 CV Plot view 281
 Deconvolved Spectrum view 317
 Elemental Composition Results view 231
 Error Log view 307
 File Header view 304
 Instrument Method view 301
 list of 46
 MultiSpectrum view 183

mzVault Search Results view 248
 NIST Search Result view 245
 Peaks List view 110
 Sample Information view 304
 Scan Header view 297
 Search Results view 245
 spectra-specific views 180
 Spectrum List view 298
 Spectrum view 180
 Status Log view 300
 Tune Method view 303
 Xtract Results view 317

W

warning messages 14
 window, FreeStyle
 description 11
 figure 18
 Info Bar pages 47
 toolbars 38
 views, list of 46
 Workspace As button 35
 Workspace Options toolbar
 buttons 39, 296
 description 38
 displaying 38, 295
 figure 39, 296
 Report area buttons 296
 workspace options, default settings for 333
 Workspace Processing toolbar
 buttons 100, 173, 266
 chromatograms 99
 description 173, 266
 displaying 99, 173, 266
 figure 99, 173, 266
 Peak Detection area (figure) 99
 Workspaces
 creating 39
 description 20
 opening 15
 Write to .RAW button
 for a chromatogram trace 91
 for exporting a spectrum to a new RAW file 151

X

XCALI-98082 i
 Xcalibur data system dependency x
 Xtract algorithm
 deconvolving spectra 309
 default settings 336
 deisotoping spectra 309

Index: Y

- description [309](#)
 - purpose [309](#)
 - results of deconvolution [318](#)
 - theoretical most abundant m/z markers [321](#)
- Xtract page [312](#)
- Xtract Results
- table parameters [318](#), [320](#)
 - view [317](#)

Y

- y-axis scaling, Chromatogram view [82](#), [98](#)
- y-axis scaling, Spectrum view [170](#)

Z

- Zoom In Y button [83](#), [172](#)
- zoom level, changing
 - chromatogram [82](#)
 - spectrum [171](#)
- Zoom Options toolbar [43–44](#)
- Zoom Out X button [83](#), [171](#)
- Zoom Out Y button [83](#), [172](#)