

FreeStyle 1.8 SP2 Release Notes

This release notes briefly list the changes in the Thermo FreeStyle™ 1.8 SP2 software, fixes to the previously reported issues, and known issues that exist in the FreeStyle 1.8 SP2 software release. Additionally, system requirements for the software installation are included in these release notes.

Contents

- [New Features](#)
- [System Requirements](#)
- [Important Information](#)
- [Resolved Issues](#)
- [Known Issues](#)
- [Trademarks](#)

For information on using the FreeStyle 1.8 SP2 application, refer to the user guide available as PDF file or the embedded Help.

New Features

This release incorporates the following new and enhanced features compared to version 1.7 SP2.

- Display peak labels produced by the Advanced Peak Determination (APD) algorithm that runs on the embedded instrument computer during data acquisition
 - Thermo Scientific™ Orbitrap Exploris™ MS Series: ICSW 4.0 or later
 - Thermo Scientific Orbitrap Tribrid™ MS Series: Tune 3.5 or later
- Elemental composition calculation based on mass only
- Thermo Scientific FAIMS Pro interface CV optimization data processing and export for TSQ™ ICSW version 3.4 or later
- Apply Xtract All deconvolution to chromatograms
- Peak area label with full numerical display
- MSn browser mass tolerance and precursor masses support 4 decimals
- Improved usability by reorganizing the ribbon menu and right-click menu options
- Backward compatibility with Thermo Scientific Foundation™ software 3.1 SP6 or later

System Requirements

These are the minimum required and recommended hardware and software configurations for operation of the FreeStyle 1.8 SP2 application.

System	Required or <i>recommended</i> configuration
Hardware	<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 Acrobat™ Reader™ 10.1 or later• Microsoft™ .NET Framework 4.7.2• Microsoft Office 2013 (for exported data)• Microsoft Windows™ 7 SP1 (64-bit) or Windows 10 (64-bit)• Thermo Foundation™ Platform 3.1 SP6 or later

Important Information

General

The FreeStyle 1.8 SP2 application does not require licensing in any form. The software may be installed on Windows 7 SP1 64-bit or Windows 10 64-bit operating systems.

Installation

You must have administrator rights to install the FreeStyle 1.8 SP2 software on your computer.

Compatibility

The FreeStyle 1.8 SP2 application requires Foundation 3.1 SP6 or later.

Install the FreeStyle 1.8 SP2 Application

Before upgrading to FreeStyle 1.8 SP2, verify the Foundation version installed on your system by going to Version Info.

Follow these instructions:

- [To upgrade to the FreeStyle 1.8 SP2 application for a system with the Foundation Platform 3.1 SP6 or later](#)
- [To upgrade to the FreeStyle 1.8 SP2 application for a system with the Foundation Platform 3.1 SP5 or earlier](#)

❖ To upgrade to the FreeStyle 1.8 SP2 application for a system with the Foundation Platform 3.1 SP6 or later

1. Download the FreeStyle 1.8 SP2 software from Flexera (<https://thermo.flexnetoperations.com/control/thmo/login>) to your local processing computer.
2. Open the Windows Control Panel and click **Programs and Features**.
3. Uninstall the previous version of the FreeStyle application.
4. Install the FreeStyle 1.8 SP2 application.

❖ To upgrade to the FreeStyle 1.8 SP2 application for a system with the Foundation Platform 3.1 SP5 or earlier

1. Download the FreeStyle 1.8 SP2 software from Flexera (<https://thermo.flexnetoperations.com/control/thmo/login>) to your local processing computer.
2. Open the Windows Control Panel and click **Programs and Features**.
3. Uninstall the FreeStyle application.
4. Uninstall any installed version of the Xcalibur application.
5. Uninstall any installed version of the Foundation Platform.
6. (Optional) Install the Xcalibur 4.2 SP1 application (or later.)
7. Install the FreeStyle 1.8 SP2 application.

Windows 10 Tip

In Windows 10, follow these steps to specify the FreeStyle application as the default application used to open all .RAW files.

1. In a File Explorer window, right-click a .RAW file and choose **Properties**.
2. On the General page of the Properties dialog box, click **Change**.
3. Select FreeStyle 1.8(C:\Program Files\Thermo\FreeStyle) and click **OK**.
4. In the Properties dialog box, click **OK**.

Supported Local Language for the Data System Computer

The FreeStyle 1.8 SP2 application was tested and supported for US-English Only locale settings.

To change the format on your computer, go to Control Panel > Region and specify the language format.

Table 1 lists the defects that were resolved between the FreeStyle 1.8 SP1 and the FreeStyle 1.8 SP2 release. The table excludes Help issues and any cosmetic fixes. In some cases, the Title has been amended or extended from the original to better describe the reported issue. The ID is the internal number assigned to each issue. Both an engineering fix and follow-up testing (verified by our product evaluation department) have resolved each of these issues.

Table 1. Resolved issues in FreeStyle 1.8 SP2

ID	Title	Severity
216691	FreeStyle application hangs on clicking the refresh button when the LC run time is longer than the MS run time.	1 - Critical
196370	In the Chromatogram, the average spectrum stops after a few averaging when the spectrum view is closed.	2 - High
197607	When multiple workspace are open, the chromatogram ranges and the chromatogram view are not in sync after applying the auto filter to one of the workspaces.	2 - High
199404	Manual peak integration does not work on the Chromatogram after applying the layout.	2 - High
200608	The Elemental composition scan mode for negative scans displays incorrect elemental composition and delta PPM.	2 - High
206576	Chromatogram Ranges - unable to display the analog signals from all detectors when the detector channel names are the same.	2 - High
199603	Manual elemental composition parameters are reverted to the default on clicking different masses on the spectrum view.	3 - Medium
196394	In the Instrument method, CSV file export displays incorrect resolution.	3 - Medium
196359	A wrong message is displayed while exporting the spectrum list when the spectrum view is closed.	3 - Medium
200071	Wrong elemental composition calculated for scan mode when relative mass labeling is applied.	3 - Medium
207517	Chromatogram - wrong y-axis label displayed for the UV channel after unchecking the MS trace.	3 - Medium

Suggested Recovery Actions

- For some issues, restarting the application is the appropriate recovery action.
- In some cases (particularly issues that arise during data acquisition), restarting applications such as the Home Page might not ensure complete recovery. Typically, restarting the data system computer resolves these issues, but some devices with error conditions might require power cycling.
- As a fix, Thermo Fisher Scientific generally does not recommend reinstalling the software or the operating system, which more commonly occurs after you install a new hard drive.

Resolved Issues

Known Issues

Feature Requests and Other Removed Items

- We do not include issues where there is insufficient information logged to successfully reproduce the reported problem.
- We do not list feature requests as software issues, regardless of the reported significance or severity of the request. Product managers evaluate logged feature requests for future releases.
- We report only discrepancies in the documented software as known issues.

Terminology

Risk	Interpretation
1 - Critical	A problem that renders the system unusable because either an entire function is unusable and no workaround exists, or use of the current system compromises data integrity or results in data loss. Catastrophic problems also include significant and non-obvious quantitative errors, and all human and instrument safety issues.
2 - High	A serious issue that does not affect data integrity (meaning data loss, corruption of data, or the wrong answer), but affects the customer's ability to use the product as designed. It can be a failure, design issue, or documentation error or omission. A workaround might or might not exist.
3 - Medium	A minor error or poor behavior of a product feature. There is probably a workaround.
4 - Low	An issue that has a limited effect on customer usage of the product; for defects with visibility so low that a customer might never see it; or for ease of use issues or other items not causing any performance degradation.

Known Defects

Table 2 table contains known defects in the software, categorized by severity with a brief abstract. The item ID is the internal number assigned to each issue. Product management assesses risk, which can differ significantly from the reported severity.

Table 2. Known defects (Sheet 1 of 2)

ID	Title	Severity
29988	Product displays blank screen during remote access.	3 - Medium
30251	Blank spectrum is displayed when background subtraction 2 range is performed in the new chromatogram view.	3 - Medium
30467	Application does not accept decimal values as expected in mass range entry.	3 - Medium
44777	Unchecking and rechecking the trace from chromatogram ranges makes to the annotated text to disappear.	3 - Medium
63283	HRAM Data not Exported if Quick, Normal Search Types Selected.	3 - Medium
82778	Chromatogram becomes blank for a particular trace type.	3 - Medium
86938	Reports - Configured width and height is not used in "To Clipboard in EMF".	3 - Medium
107758	NIST similarity search type Neutral Loss results mismatch in FreeStyle and NIST browser.	3 - Medium
113915	Elemental results table does not track selected spectrum.	3 - Medium
163242	Undo zoom does not perform as expected upon toggling and un-toggling EIC mass and EIC range.	3 - Medium
172824	Xtract - Incorrect monoisotopic mass for large protein.	3 - Medium
178143	Unexpected message After Changing Something in Default Options When NIST Info Tab is Open.	3 - Medium
178155	Communicator Bar "No Results" Message Is Not Cleared When There Is Results.	3 - Medium
179721	Refresh issue in info bar Xtract <i>m/z</i> range fields.	3 - Medium
180146	Structure annotations are getting lost upon performing Undo Zoom.	3 - Medium

Table 2. Known defects (Sheet 2 of 2)

ID	Title	Severity
181647	Xtract - Monoisotopic Mass Mode is Off by 1amu for CAH Protein.	3 - Medium
187534	FreeStyle crashes when the layout is saved after moving Chromatogram as tabbed view with Spectrum View.	3 - Medium

Trademarks

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

Microsoft, Windows, and Excel are registered trademarks of Microsoft Corporation in the United States and other countries.

Adobe, Acrobat, and Reader are registered trademarks of Adobe Systems Inc. in the United States and other countries.

NIST is a registered trademark of the National Institute of Standards and Technology in the United States.

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