



thermo scientific

Beverage resource guide for commercial testing laboratories

ThermoFisher
SCIENTIFIC




Fast, efficient analysis

For all beverages, the compositional quality and safety must be monitored to help track contamination, adulteration, and product consistency, as well as to ensure regulatory compliance from raw ingredients (water, additives, and fruits) to the final product.



Thermo Fisher Scientific is a recognized leader in providing analytical solutions. From carbohydrates and sugar substitutes, to vitamins and additives, we are unique in our commitment to provide fast, accurate testing for all applications performed in commercial testing laboratories.

Compendiums included in this directory

| | |
|---|----|
| Notebook of peer reviewed articles and papers | 4 |
| Additives in beverages | 5 |
| Beer | 6 |
| Bottled water | 7 |
| Composition analysis | 8 |
| Functional waters/drinks | 9 |
| Milk | 10 |
| Wine | 11 |



Key to the instrumentation acronyms

- Learn more
-  Beverage Testing Learning Center
 -  Beverage testing instrumentation

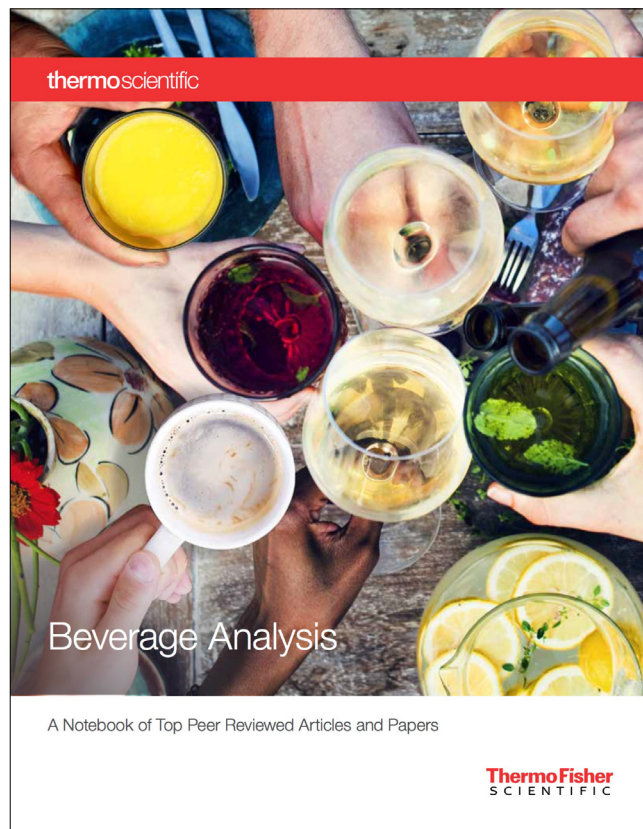


A key to the instrumentation acronyms

| Acronym | Description |
|-------------------|--|
| AA | Atomic Absorption Spectroscopy |
| GC-MS | Gas Chromatography with Single Quadrupole Mass Spectrometry Detection |
| GC-MS/MS | Gas Chromatography with Triple Quadrupole Mass Spectrometry Detection |
| HPLC-PAD | High Performance Liquid Chromatography with Pulsed Amperometry Detection |
| HPLC-CAD | High Performance Liquid Chromatography with Charged Aerosol Detection |
| HPLC-DAD | High Performance Liquid Chromatography with Diode Array Detection |
| HPLC-ECD | High Performance Liquid Chromatography with Electrochemical Detection |
| HPLC-Fluorescence | High Performance Liquid Chromatography with Fluorescence Detection |
| HPLC-UV | High Performance Liquid Chromatography with Ultraviolet Detection |

| Acronym | Description |
|----------------------------|---|
| IC-Suppressed Conductivity | Ion Chromatography with Suppressed Conductivity Detection |
| IC-UV | Ion Chromatography with Ultraviolet Detection |
| ICP-MS | Inductively Coupled Plasma Mass Spectrometry |
| ICP-OES | Inductively Coupled Plasma Optical Emission Spectrometry |
| Ion Exclusion-PAD | Ion Exclusion Chromatography with Pulsed Amperometry Detection |
| IRMS | Isotope Ratio Mass Spectrometry |
| LC-MS/MS | High Performance Liquid Chromatography with Triple Quadrupole Mass Spectrometry Detection |
| LC-HRAM | High Performance Liquid Chromatography with High Resolution Accurate Mass using Orbitrap Technology |
| UHPLC-MS/MS | Ultra High Performance Liquid Chromatography with Triple Quadrupole Mass Spectrometry Detection |

Notebook of peer reviewed articles and papers



[View the full compendium](#)

[Learn more:](#) [Beverage Testing Learning Center](#)

Overview

The power of peer reviewed articles cannot be underestimated. Years of research within the beverage industry have led to the discovery of anti-oxidants in coffee and wine, the origin of raw ingredients when analyzing for beverage authenticity, and pesticides in fruit juices.

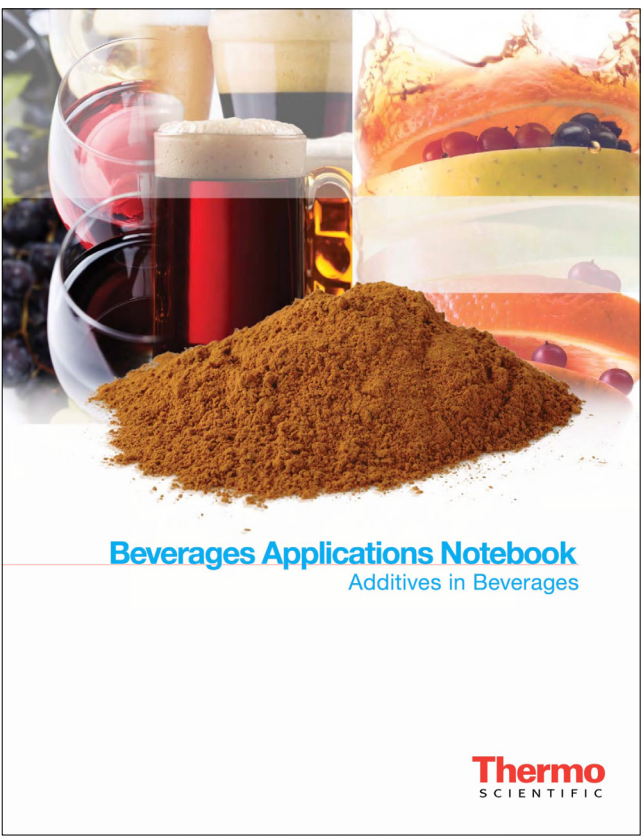
This comprehensive notebook represents a broad spectrum of publications that can be accessed easily online, specifically for the beverage industry. A collection of beverage analyses is presented, providing an insight into analytical chromatography research that is taking place today within the community. Techniques described include gas chromatography-mass spectrometry, liquid chromatography-mass spectrometry, ion chromatography, automated photometry, and trace elemental analyses.

Compendium contents

| Analysis | Instrumentation type |
|-------------------------------------|---|
| Beer and cider | Liquid Chromatography (LC) |
| Bottled water and functional drinks | Gas Chromatography (GC) |
| Coffee and cocoa | Ion Chromatography (IC) |
| Fruit juice | Triple Quadrupole Mass Spectrometry (LC, GC and IC) |
| Milk | High Resolution Accurate Mass |
| Soft drinks | ICP-OES, ICP-MS |
| Spirits | Atomic Absorption |
| Tea | Isotope Ratio Mass Spectrometer |
| Wine | Automated Discrete Analyzers |
| | Fourier Transform infrared spectroscopy |
| | UV-Vis Spectroscopy |



Additives in beverages



Overview

For all beverages, the compositional quality and safety must be monitored to help track contamination, adulteration, and product consistency, and to ensure regulatory compliance from raw ingredients (water, additives, and fruits) to the final product.

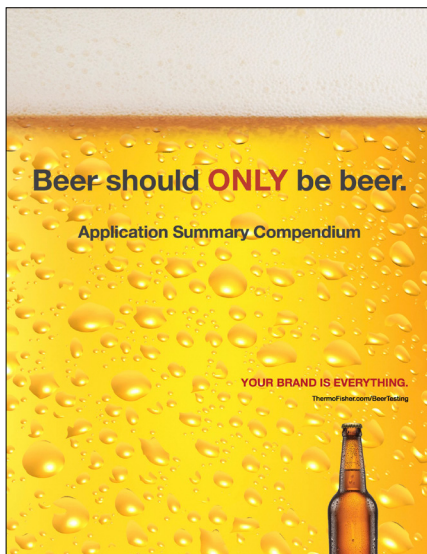
Compendium contents

| Analysis | Instrumentation type |
|--------------------|----------------------------|
| Benzoate | IC-Suppressed Conductivity |
| Dyes | HPLC-DAD |
| Mogroside V | HPLC-UV, HPLC-CAD |
| Steviol glycosides | HPLC-UV, HPLC-CAD |
| Sulfite | Ion Exclusion-PAD |
| Vitamins | UHPLC-MS/MS |

[View the full compendium](#)

Learn more: [Beverage Testing Learning Center](#)

Beer



[View the full compendium](#)



[View the full compendium](#)

Overview

Beer is the most widely consumed alcoholic beverage in the world. Beer is typically brewed from four basic ingredients: water, a starch source, brewer's yeast, and a flavoring agent such as hops. Many varieties of beer result from differences in these ingredients, the additives used, and the brewing process. Thermo Scientific instruments help in monitoring the brewing process and also testing the quality, consistency, and purity of the final product.

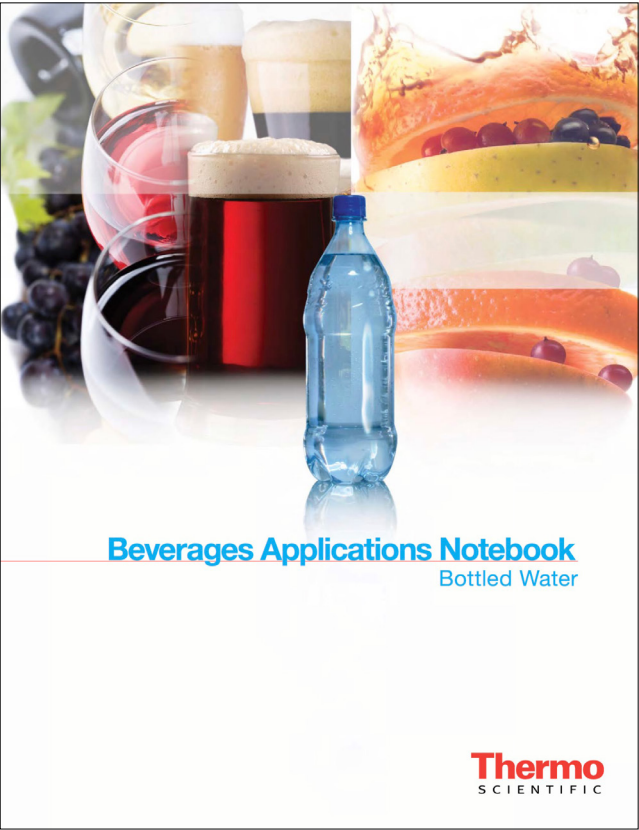
Compendium contents

| Analysis | Instrumentation type |
|---|--|
| IC analysis (anions, cations) | IC-Suppressed conductivity |
| Carbohydrates | HPAE-PAD |
| Organic acids | IC-Suppressed Conductivity |
| Acetaldehyde | Automated Discrete Photometry Analyzer |
| Beta-glucan | Automated Discrete Photometry Analyzer |
| Free Amino nitrogen | Automated Discrete Photometry Analyzer |
| Isohumulones (Iso- α -acids) | HPLC-UV |
| Sulfur dioxide | Automated Discrete Photometry Analyzer |
| Nitrosamines | GC-MS/MS |
| Chalconoids and bitter acids | HPLC-UV, HPLC-ECD |
| Multi-mycotoxin screening | LC-HRAM |
| Beer polyphenols, proanthocyanidins, and bitter acids | HPLC-UV, HPLC-ECD |
| Elemental contaminants | ICP-OES |

Learn more: [Beer Testing Learning Center](#)



Bottled water



Overview

Clean, safe, without any contamination—that is what bottled water should be. We help manufacturers of bottled water, commercial testing labs, and regulatory agencies with our instruments to test for all this and more. Our instruments and solutions help in identifying disinfection by products, water contamination, electrolytes, and phthalates.

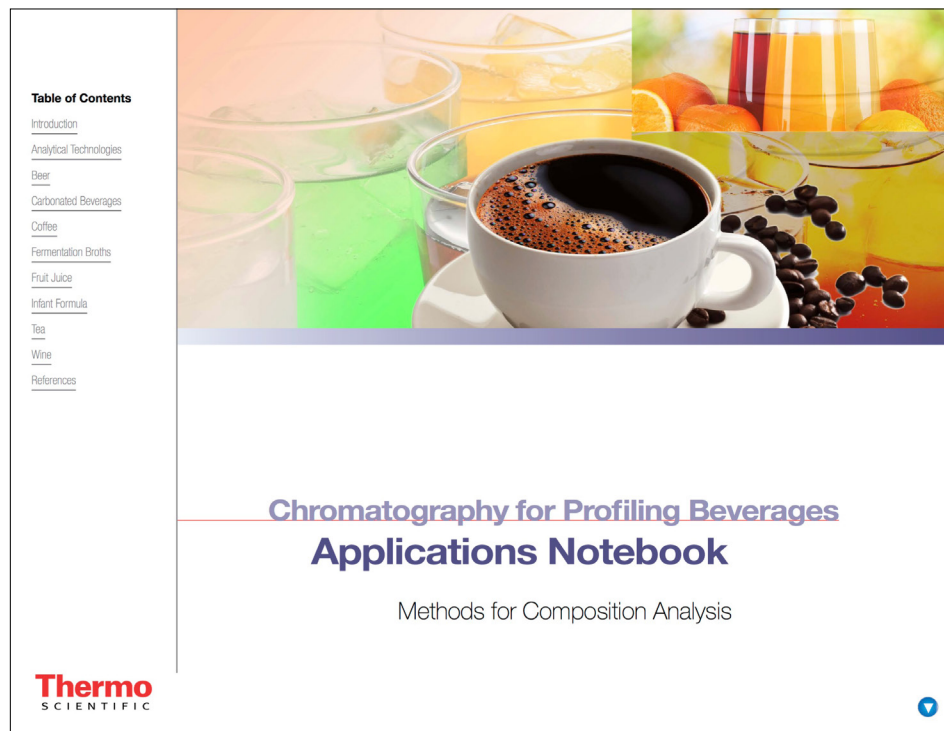
Compendium contents

| Analysis | Instrumentation type |
|---------------------------------|----------------------------|
| Oxyhalides and bromide | IC-Suppressed Conductivity |
| Chlorite, bromate, and chlorate | IC-Suppressed Conductivity |
| Phenols | HPLC-UV |

[View the full compendium](#)

Learn more: [Bottled Water Testing Learning Center](#)

Composition analysis



[View the full compendium](#)

Learn more: [Beverage Testing Learning Center](#)

Overview

The measurement of the different compounds found in a wide range of non-alcoholic and alcoholic drinks.

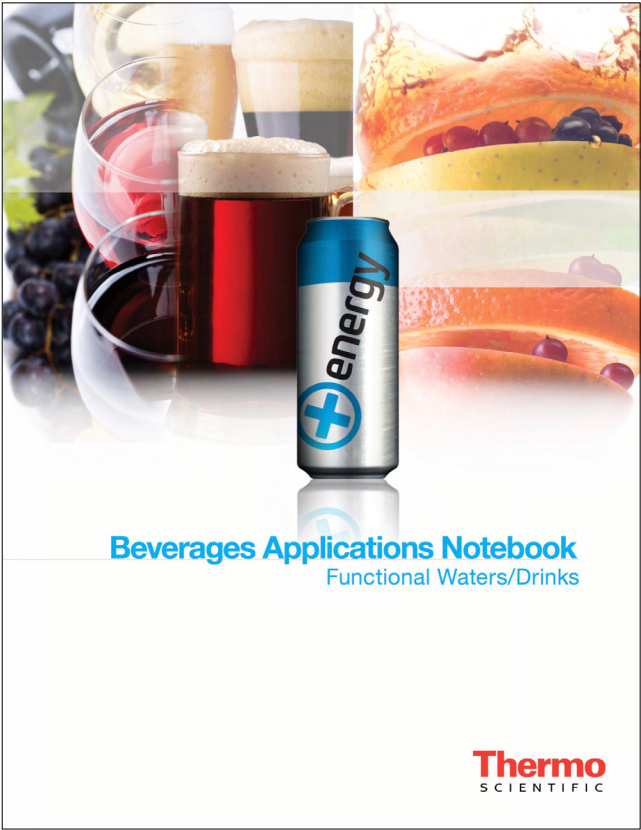
Included is an overview of instrumentation for the testing of beverages and a multitude of references from peer reviewed journals using our HPLC, ion chromatography, and sample preparation solutions.

Compendium contents

| Analysis | Instrumentation type |
|----------------------|---|
| Beer | HPAE-PAD, HPLC-UV, HPLC-ECD, IC-Suppressed Conductivity |
| Carbonated beverages | IC-Suppressed Conductivity |
| Coffee | HPAE-PAD, IC-Suppressed Conductivity, HPLC-UV |
| Fermentation broths | HPAE-PAD, IC-Suppressed Conductivity |
| Fruit juice | HPAE-PAD, IC-Suppressed Conductivity, HPLC-ECD |
| Infant formula | HPAE-PAD, IC-Suppressed Conductivity, HPLC-DAD |
| Tea | HPLC-UV, IC-Suppressed Conductivity |
| Wine | HPLC-UV, IC-Suppressed Conductivity, HPLC-ECD |



Functional waters/drinks



Overview

The beverage industry is growing each year with the introduction of new functional beverages, such as vitamin fortified water, energy drinks, anti-aging water, and sports drinks. With this growth come many more analytical challenges. Our instruments and applications help in overcoming these challenges.

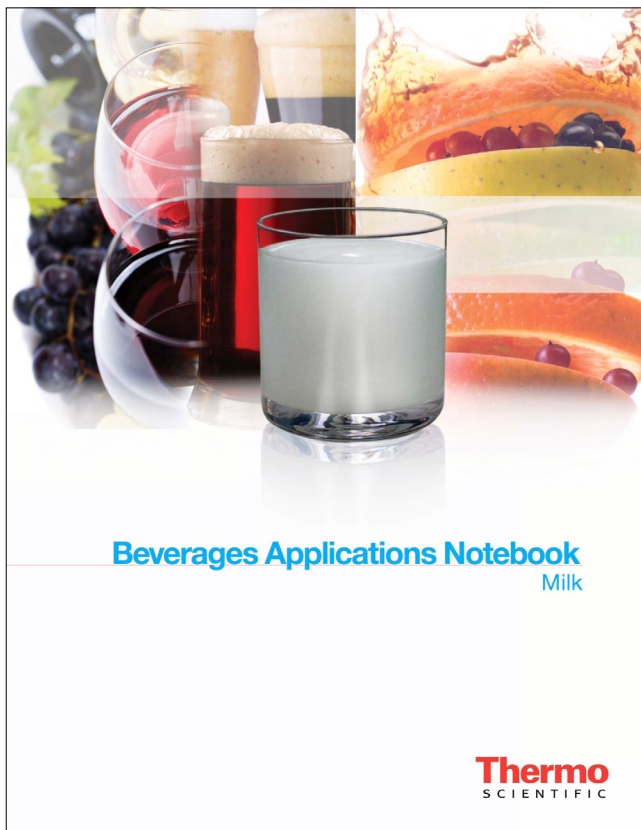
Compendium contents

| Analysis | Instrumentation type |
|--------------------------------|----------------------|
| Water and fat-soluble vitamins | HPLC-DAD |
| Glucosamine | HPAE-PAD |
| Vitamin B12 | HPLC-UV |

[View the full compendium](#)

Learn more: [Functional Beverage Testing Learning Center](#)

Milk



Overview

People trust their milk to be safe, consistent, and unadulterated. Our instruments help in adhering to food label claims, product consistency, and purity, as well as identifying any contaminants and adulterants. Discover our solutions from carbohydrate analysis to vitamin D analysis to choline analysis in infant formula.

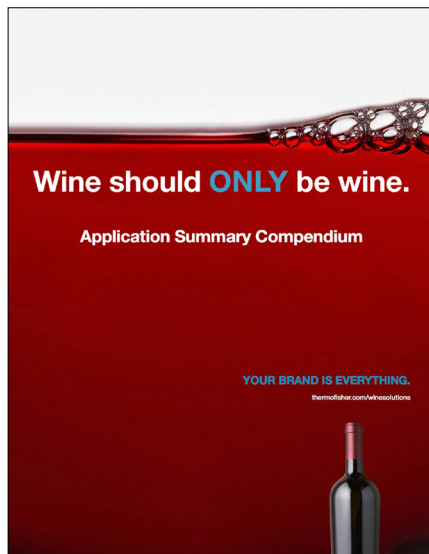
Compendium contents

| Analysis | Instrumentation type |
|-----------------|--------------------------------|
| Choline | IC-Suppressed Conductivity |
| Fat extraction | Accelerated Solvent Extraction |
| Iodide | HPAE-PAD |
| Lactose | HPAE-PAD |
| Melamine | HPLC-UV, IC-UV |
| Nitrate/Nitrite | IC-Suppressed Conductivity |
| Sialic acids | HPAE-PAD, HPLC-Fluorescence |

[View the full compendium](#)

Learn more: [Milk Testing Learning Center](#)

Wine



[View the full compendium](#)



[View the full compendium](#)

Overview

People trust their wine to be exactly as they expect: unadulterated, safe, and consistent. Our customers know how long it takes to build a brand's reputation and how quickly one can disappear. So, they come to us and our widest instrumentation portfolio—ion, liquid, and gas chromatography, metal analysis, mass spectrometry, discrete analyzers, and data management—for accurate, reliable answers for their wine testing needs.

Compendium contents

| Analysis | Instrumentation type |
|------------------------------|--|
| Catechins and phenolic acids | HPLC-UV |
| Pesticide residues | LC-MS/MS |
| Isotopes | GasBench II-IRMS |
| Phenolics, flavonoids | LC-HRAM |
| Iron, copper and zinc | AA |
| Nutrients and trace elements | ICP-OES |
| Impurities | GC-MS |
| Free sulfite | Automated Discrete Photometry Analyzer |
| Acetaldehyde | Automated Discrete Photometry Analyzer |
| Total acidity | Automated Discrete Photometry Analyzer |
| Glycerol | Automated Discrete Photometry Analyzer |

Learn more: [Wine Testing Learning Center](#)



Thermo Fisher Scientific is a recognized leader in providing analytical solutions. From carbohydrates and sugar substitutes, to vitamins and additives, we are unique in our commitment to provide fast, accurate testing and labeling information for all applications in this industry.

Find out more at thermofisher.com/powerofpartnership

ThermoFisher
SCIENTIFIC

© 2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. **A173834-EN 1120M**