

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



Chromatography for Foods and Beverages

Flavors, Colorants and Additives Analysis Applications Notebook

Analytical Methods for Food Ingredients

Table of Contents

Introduction

Analytical Technologies

Artificial Colorants

Blue Dyes

Caramel

General Dye Method

Global Method

Red Dyes

Sudan Dyes

Yellow Dyes

Natural Colorants

Turmeric

Flavorants

Amadori Compounds

Capsaicin

Quinine

Vanilla

Miscellaneous

Caffeine

Choline

Glucosamine

Iodide

Polyphosphates

Preservatives

Benzoate

Nitrate and Nitrite

Sulfites

Soft Drinks

References



Additives Including Colorants, Flavorants, and Preservatives

Introduction

Food additives are substances added to food to preserve flavor or enhance its taste, texture, and appearance. With the advent of processed foods many more additives are being used, both natural and artificial. Additives can maintain product quality and freshness (to stop deterioration, rancidity, and spoilage); improve or maintain the nutritional quality (to prevent diseases such as goiter, pellagra and rickets); make foods more appealing (make them look and taste good); and aid in the processing and preparation of foods.

Table of Contents

Introduction

Analytical Technologies

Artificial Colorants

Blue Dyes

Caramel

General Dye Method

Global Method

Red Dyes

Sudan Dyes

Yellow Dyes

Natural Colorants

Turmeric

Flavorants

Amadori Compounds

Capsaicin

Quinine

Vanilla

Miscellaneous

Caffeine

Choline

Glucosamine

Iodide

Polyphosphates

Preservatives

Benzoate

Nitrate and Nitrite

Sulfites

Soft Drinks

References



Analytical Technologies



High-Performance Liquid Chromatography

Thermo Scientific™ Vanquish™ UHPLC System and Thermo Scientific™ Dionex™ UltiMate™ 3000 UHPLC+ systems offer excellent chromatographic performance, operational simplicity and unrivaled flexibility. Choose from a wide range of standard and unique specialty detectors to extend your laboratory's analytical capabilities.

Table of Contents

Introduction

Analytical Technologies

Artificial Colorants

Blue Dyes

Caramel

General Dye Method

Global Method

Red Dyes

Sudan Dyes

Yellow Dyes

Natural Colorants

Turmeric

Flavorants

Amadori Compounds

Capsaicin

Quinine

Vanilla

Miscellaneous

Caffeine

Choline

Glucosamine

Iodide

Polyphosphates

Preservatives

Benzoate

Nitrate and Nitrite

Sulfites

Soft Drinks

References

Thermo
SCIENTIFIC



The next generation in UHPLC innovations

The Vanquish system takes high-end UHPLC to a new level, offering more resolution while meeting the throughput demands of modern laboratories.

The system delivers better separations, more results and easier interaction, simultaneously, without compromise.



Analytical Technologies

The Vanquish UHPLC System

Delivering the new standard in UHPLC

- More powerful separations with 1500 bar of pump pressure at flow rates up to 5 mL/min
- Industry-leading flow and gradient precision
- Excellent injections up to 100 μ L in 0.01 μ L increments
- Automated workflows with barcode reading for simplified setup and tracking
- Maximum sample capacity with up to 23 well plates, or 8832 samples
- More confident separations with a wide temperature range of 5 $^{\circ}$ C to 120 $^{\circ}$ C for two thermostating modes and active column pre-heating for improved precision
- UV detection with linear response up to 3000 mAu and noise levels as low as 3 μ Au
- Thermo Scientific™ LightPipe™ technology assures lowest peak dispersion with UV detection
- Available Vanquish Charged Aerosol detector for quantification of non-chromophoric compounds



Vanquish Diode Array Detector with LightPipe technology

Learn more at www.thermoscientific.com/vanquish

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Analytical Technologies

UHPLC Portfolio

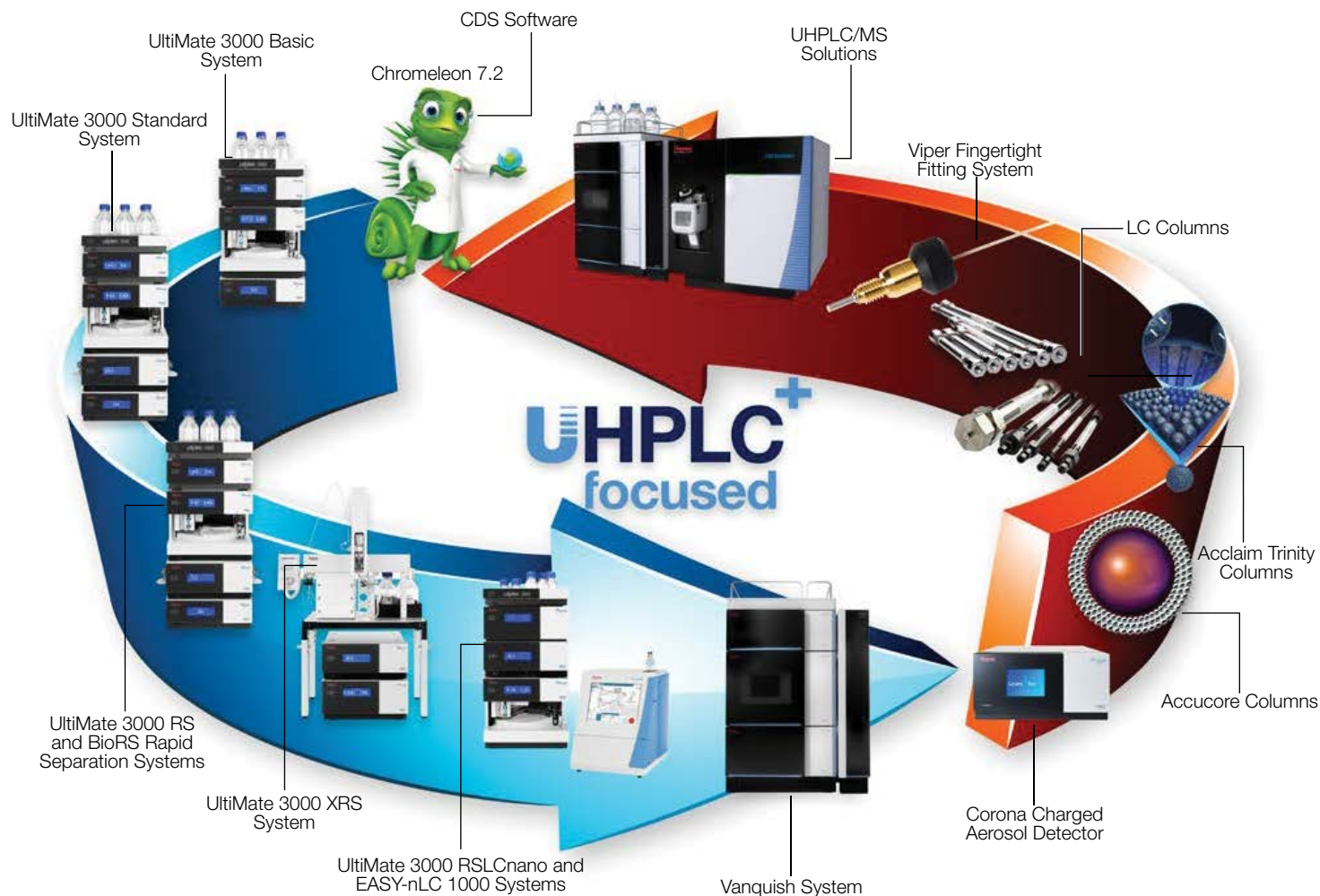


Table of Contents

Introduction

Analytical Technologies

Artificial Colorants

Blue Dyes

Caramel

General Dye Method

Global Method

Red Dyes

Sudan Dyes

Yellow Dyes

Natural Colorants

Turmeric

Flavorants

Amadori Compounds

Capsaicin

Quinine

Vanilla

Miscellaneous

Caffeine

Choline

Glucosamine

Iodide

Polyphosphates

Preservatives

Benzoate

Nitrate and Nitrite

Sulfites

Soft Drinks

References

Analytical Technologies



UltiMate 3000 UHPLC+ Systems

Best-in-class HPLC systems for all your chromatography needs

UltiMate 3000 UHPLC+ Systems provide excellent chromatographic performance while maintaining easy, reliable operation. The basic and standard analytical systems offer ultra HPLC (UHPLC) compatibility across all modules, ensuring maximum performance for all users and all laboratories.

Covering flow rates from 20 nL/min to 10 mL/min with an industry-leading range of pumping, sampling, and detection modules, UltiMate 3000 UHPLC+ Systems provide solutions from nano to semipreparative, from conventional LC to UHPLC.

Superior chromatographic performance

- UHPLC design philosophy throughout nano, standard analytical, and rapid separation liquid chromatography (RSLC)
- 620 bar (9,000 psi) and 100 Hz data rate set a new benchmark for basic and standard analytical systems
- RSLC systems go up to 1000 bar and data rates up to 200 Hz
- ×2 Dual System for increased productivity solutions in routine analysis
- Fully UHPLC compatible advanced chromatographic techniques
- Thermo Scientific™ Dionex™ Viper™ and nanoViper™ fingertight fittings—the first truly universal, fingertight fitting system even at UHPLC pressures

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

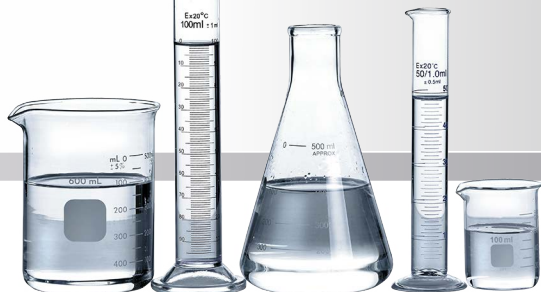
[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Analytical Technologies



UltiMate 3000 UHPLC+ Systems

We are uniquely focused on making UHPLC technology available to all users, all laboratories, and for all analytes.



Rapid Separation LC Systems

The extended flowpressure footprint of the RSLC system provides the performance for ultrafast high-resolution and conventional LC applications.



RSLCnano Systems

The Rapid Separation nano LC System (RSLCnano) provides the power for high resolution and fast chromatography in nano, capillary, and micro LC.



Standard LC Systems

Choose from a wide variety of standard LC systems for demanding LC applications at nano, capillary, micro, analytical, and semipreparative flow rates.



Basic LC Systems

UltiMate 3000 Basic LC Systems are UHPLC compatible and provide reliable, high performance solutions to fit your bench space and your budget.

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Thermo
SCIENTIFIC



Charged Aerosol Detection

Charged Aerosol Detection provides near universal detection independent of chemical structure for non- or semi-volatile analytes with HPLC and UHPLC. Thermo Scientific™ Dionex™ Corona™ Veo™ and Vanquish Charged Aerosol detectors are ideally suited as a primary detector for any laboratory, while providing complementary data to UV or MS methods. No other LC detector available today can match the performance of a Corona Veo detector.

- High sensitivity – single-digit nanogram on column
- Consistent response – independent of chemical structure
- Wide dynamic range – to four orders of magnitude or greater
- Simple to use – easy to integrate with any HPLC/UHPLC system

Charged aerosol detectors give the simplicity, reproducibility and performance required for a full range of applications from basic research to manufacturing QC/QA. With charged aerosol detection you get predictable responses to measure analytes in direct proportion to their relative amounts for quantitation without actual standards.

This detector offers the flexibility to use reversed-phase gradients, as well as normal phase and HILIC modes of separation on any LC system. And, in many cases eliminates the need for derivatization or sample pre-treatment to provide real dilute-and-shoot simplicity.

Analytical Technologies

Advanced Detection Capabilities



Corona Veo Charged Aerosol Detector



Vanquish system with Charged Aerosol Detector

Learn more at www.thermoscientific.com/CAD

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



CoulArray Multi-electrode Array Detector

The Thermo Scientific™ Dionex™ CoulArray™ Multi-electrode Array detector is the only practical multi-channel electrochemical detection system that allows you to measure multiple analytes simultaneously, including those that are chromatographically unresolved. The CoulArray detector delivers the widest dynamic range of any available electrochemical detector with unmatched selectivity for detection of trace components in complex matrixes, even when used with aggressive gradients.

- Measures analytes from femtomole to micromole levels
- Greatly simplify sample preparation and eliminate interferences
- Simultaneously analyze multiple analytes in very complex samples
- Easily produce qualitative information for compound identification

Multiple system configurations offer 4, 8, 12, or 16 channels that can be upgraded anytime. The unique data acquisition and processing software uses automatic signal ranging and a unique patented baseline correction algorithms to provide identification and quantitation of single or multiple analytes and powerful 3D data for quick sample fingerprint confirmation with integration to pattern recognition platforms.

With the power of coulometric array technology, the CoulArray detector can give you the qualitative data of a optical PDA with 1,000 fold greater sensitivity to profile the characteristic qualities of products, determine integrity, identify adulteration and even evaluate competitors' products.

Advanced Detection Capabilities



CoulArray Multi-electrode Array Detector

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



Advanced Detection Capabilities

RefractoMax 521 Refractive Index Detector

The Thermo Scientific RefractoMax 521 Refractive Index Detector from ERC Inc. This detector, in combination with the UltiMate 3000 system, is the right choice for the isocratic analysis of sugars, polymers, and fatty acids. It features fast baseline stabilization and excellent reproducibility, combined with high sensitivity. The RefractoMax 521 is fully controlled by Thermo Scientific™ Dionex™ Chromeleon™ Chromatography Data System Software (CDS), and can also operate in stand-alone mode.

- The detector is highly sensitive and applicable universally. It provides very stable baselines with a drift of 0.2 μ RIU/h and a noise specification of 2.5 nRIU or less
- The optical bench, thermostatically regulated from 30 °C to 55 °C, and the superior signal-to-noise ratio ensure highly precise measurement results

- The extended flow rate range from 1 mL/min up to 10 mL/min and the operating range of 1.00 to 1.75 RIU enable the use of this detector for a wide range of applications
- Applications include the analysis of all compounds with low UV-Vis activity, such as alcohols, mono- and polysaccharides, esters, fatty acids, or polymers
- An Auto Set-up function automates purging, equilibration, autozero, and the control baseline stability and noise
- Operation with Chromeleon CDS makes the detector easy to use and ensures maximum productivity in instrument control, data processing, and reporting of results



RefractoMax 521 Refractive Index Detector

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



UltiMate 3000 Diode Array and Multiple-Wavelength Detectors

The Thermo Scientific Dionex UltiMate DAD 3000 detector is a high-resolution, 1024-element diode array detector (DAD) available in Rapid Separation (200 Hz) and Standard (100 Hz) versions. It operates with Chromeleon CDS software to provide a variety of spectra views, including 3-D plotting and automated chromatogram handling. The high resolution and low-noise performance of the DAD-3000 family makes it ideal for the most sensitive and accurate library searches and peak purity analyses.

The detector is also available as a multiple wavelength detector (MWD) in Standard (100 Hz) and Rapid Separation (200 Hz) versions.

- Data collection at up to 200 Hz using a maximum of eight single-wavelength data channels and one 3-D field (3-D only with DAD-3000 (RS)) for best support of ultrafast separations
- Standard versions operate at up to 100 Hz data collection rate for optimum support of 62 MPa (9000 psi) UltiMate 3000 Standard systems
- Accurate compound confirmation with a 1024-element, high resolution photodiode array
- Flexibility in both UV and Vis applications with 190–800 nm wavelength range
- Low-noise over the full spectral range using deuterium and tungsten lamps
- Fast and accurate wavelength verification using a built-in holmium oxide filter

Advanced Detection Capabilities

- The detector can be upgraded with the UltiMate PCM 3000 for accurate monitoring pH gradients
- Excellent reliability and reproducibility with low baseline drift (typically < 500 μ AU/h)
- Simplified routine maintenance with front access to pre-aligned cells and lamps
- ID chips on flow cells and lamps for identification and life-span monitoring
- Chromeleon CDS software for full control and flexible data handling
- Front-panel display for easy monitoring of detector status to maximize uptime
- Flow cells for semi-micro, semi-analytical, analytical, and semi-preparative applications
- Flow cells available in stainless steel and biocompatible versions



UltiMate 3000 DAD-3000 Diode Array Detector

Table of Contents

Introduction

Analytical Technologies

Artificial Colorants

Blue Dyes

Caramel

General Dye Method

Global Method

Red Dyes

Sudan Dyes

Yellow Dyes

Natural Colorants

Turmeric

Flavorants

Amadori Compounds

Capsaicin

Quinine

Vanilla

Miscellaneous

Caffeine

Choline

Glucosamine

Iodide

Polyphosphates

Preservatives

Benzoate

Nitrate and Nitrite

Sulfites

Soft Drinks

References



Advanced Detection Capabilities

UltiMate 3000 Electrochemical Detector

Electrochemical detection delivers high sensitivity for neurotransmitter analysis, simplicity and robustness for pharmaceutical or clinical diagnostics, and the selectivity for the characterization of complex samples such as natural products, biological tissues and fluids. For today's researcher, there is a continuing need for detecting vanishingly small quantities of analyte and often in complex samples. Because electrochemical detection measures only compounds that can undergo oxidation or reduction it is both highly sensitive and very selective.

The Thermo Scientific Dionex UltiMate 3000 Electrochemical Detector, designed by the pioneers of coulometric electrochemical detection, delivers state-of-the-art sensor technologies complete with an entire range of high performance and ultra-high performance LC systems optimized for electrochemical detection. The UltiMate 3000 ECD-3000RS takes electrochemical detection to the next level with UHPLC compatibility, total system integration, and selection of detection mode, all with unprecedented operational simplicity.

Features include:

- Detection Modes – choose from DC and PAD for optimum analyte response
- Choice of sensors – both coulometric and amperometric sensors to meet the demands of any application
- UHPLC compatibility – ultralow peak dispersion and high data acquisition rates for conventional or fast, high resolution chromatography
- Modularity – easily expandable to multiple independent sensors for unrivaled flexibility
- Autoranging – simultaneously measure both low and high levels of analytes without losing data
- SmartChip™ technology – easy operation with automatic sensor recognition, event logging and electrode protection



UltiMate 3000 Electrochemical Detector

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



Advanced Detection Capabilities

UltiMate 3000 Fluorescence Detector

The Thermo Scientific Dionex UltiMate 3000 FLD-3000 is a high-sensitivity fluorescence detector series for UltiMate 3000 HPLC systems. It is available in Rapid Separation (RS) and Standard (SD) versions. The optics of the FLD-3000 series provide maximum stray-light suppression for best detection sensitivity. Operated with the Chromeleon CDS software, the detector provides automated qualification, various tools for method development, and instrument wellness monitoring for ease of use, maximum uptime, and the highest degree of regulatory compliance.

- Data collection at up to 200 Hz for optimal support of even the fastest UHPLC separations (FLD-3400RS)
- Standard detectors operate at up to 100 Hz data rate for optimum support of 62 MPa (9,000 psi) UltiMate 3000 standard systems
- Lowest limits of detection with a Raman signal-to-noise ratio (S/N): > 550 ASTM (> 2100 using dark signal as noise reference)

- Unsurpassed reproducibility with active flow cell temperature control for stable fluorophore activity independent of changes in ambient temperature
- Long-life xenon flash lamp for highest sensitivity and long-term operation without the need for frequent lamp changing
- Optional second photomultiplier (PMT) for unique Dual-PMT operation, offering an extended wavelength range up to 900 nm without sacrificing sensitivity in the standard wavelength range
- Two-dimensional (2D) or three dimensional (3D) excitation, emission, or synchro scans to provide the highest degree of flexibility for method development or routine sample characterization
- Innovative Variable Emission Filter for real-time compound-related sensitivity optimization (FLD-3400RS only)
- Large front-panel display for easy monitoring of the detector status
- Two flow-cell sizes for easy optimization to application requirements: the 8 μ L flow cell is ideal for trace analysis, and the 2 μ L flow cell offers best peak resolution with narrow-bore HPLC and UHPLC columns



Ultimate 3000 Fluorescence Detector

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



UltiMate 3000 Variable Wavelength Detectors

The Thermo Scientific Dionex UltiMate 3000 VWD-3000 is a variable wavelength detector (VWD) series for industry leading UV-Vis detection. The forward optics design and wide range of available flow cells ensure optimal performance over a flow rate range of five orders of magnitude. Automated qualification, performance optimization, and instrument wellness monitoring deliver maximum uptime, simplify work-flow, and give you full confidence in your analytical results. The detector is available in a standard 100 Hz (VWD-3100) and a 200 Hz Rapid Separation version (VWD-3400RS) for the most challenging UHPLC applications.

High-Performance UV-Vis Detection

- The VWD-3400RS variant provides data collection rates of up to 200 Hz for optimal support of today's and tomorrow's UHPLC separations
- The VWD-3100 standard detector operates at up to 100 Hz data rate for optimum support of 62 MPa (9000 psi) UltiMate 3000 Standard systems
- Superior detection of trace analytes with low noise ($< -2.0 \mu\text{AU}$) and drift ($< 100 \mu\text{AU/h}$)
- The detector's large linearity range of up to 2.5 AU is ideal for applications with widely varying analyte concentrations
- Up to four absorption channels (VWD-3400RS) and spectral scans support effective method development
- Active temperature control of optics and electronics for data acquisition independent of ambient conditions

Advanced Detection Capabilities

- Front panel access for quick and easy lamps and flow cells changes
- Automated qualification monitoring for full regulatory compliance
- Large front panel display for monitoring the detector status even from a distance
- Maximize uptime using predictive performance-based on monitoring the life cycle of detector lamps
- The detector can be upgraded with the Thermo Scientific Dionex pH/Conductivity Monitor (PCM-3000) for accurate and precise pH- and conductivity monitoring
- Unique 45 nL ultra-low dispersion UV monitor for dispersion-free UV detection in LC/MS



UltiMate 3000 VWD-3400 Variable Wavelength Detector.

Table of Contents

Introduction

Analytical Technologies

Artificial Colorants

Blue Dyes

Caramel

General Dye Method

Global Method

Red Dyes

Sudan Dyes

Yellow Dyes

Natural Colorants

Turmeric

Flavorants

Amadori Compounds

Capsaicin

Quinine

Vanilla

Miscellaneous

Caffeine

Choline

Glucosamine

Iodide

Polyphosphates

Preservatives

Benzoate

Nitrate and Nitrite

Sulfites

Soft Drinks

References



Analytical Technologies



Ion Chromatography

Thermo Scientific Dionex IC systems have led the analytical instrument industry for over 30 years with solutions that represent state-of-the-art technological advancements and patented technologies.

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



Innovative Ion Chromatography Solutions

Our High-Pressure™ Ion Chromatography (HPIC™) systems include the Thermo Scientific Dionex ICS-5000+ HPIC system, which is optimized for flexibility, modularity, and ease-of-use, combining the highest chromatographic resolution with convenience. In addition, the Thermo Scientific Dionex ICS-4000 Capillary HPIC system is the world's first commercially available dedicated capillary high-pressure Reagent-Free™ (RFIC™) IC system. The Dionex ICS-4000 system is always ready for the next analysis, delivering high-pressure IC on demand.

Reagent-Free IC systems eliminate daily tasks of eluent and regenerant preparation in turn saving time, preventing errors, and increasing convenience. RFIC-EG systems use electrolytic technologies to generate eluent on demand from deionized water, and to suppress the eluent back to

pure water to deliver unmatched sensitivity. RFIC-ER systems are designed to use carbonate, carbonate/ bicarbonate, or MSA eluents for isocratic separations.

At the heart of our ion chromatography portfolio is a unique set of column chemistries that provide high selectivities and efficiencies with excellent peak shape and resolution. Thermo Scientific™ Dionex™ IonPac™ chromatography columns address a variety of chromatographic separation modes including ion exchange, ion exclusion, reversed-phase ion pairing, and ion suppression. Our column chemistries are designed to solve specific applications, and we offer a variety of selectivities and capacities for simple and complex samples. Additionally, our Dionex IonPac column line is available in standard bore, microbore and capillary formats for the ultimate application flexibility.



Thermo Scientific Dionex IC instrument family

Analytical Technologies

IC and RFIC Systems

Table of Contents

Introduction

Analytical Technologies

Artificial Colorants

Blue Dyes

Caramel

General Dye Method

Global Method

Red Dyes

Sudan Dyes

Yellow Dyes

Natural Colorants

Turmeric

Flavorants

Amadori Compounds

Capsaicin

Quinine

Vanilla

Miscellaneous

Caffeine

Choline

Glucosamine

Iodide

Polyphosphates

Preservatives

Benzoate

Nitrate and Nitrite

Sulfites

Soft Drinks

References



Analytical Technologies



Mass Spectrometry

We provide advanced integrated IC/MS and LC/MS solutions with superior ease-of-use and modest price and space requirements. UltiMate 3000 System Wellness technology and automatic MS calibration allow continuous operation with minimal maintenance. The Dionex ion chromatography family automatically removes mobile phase ions for effort-free transition to MS detection.

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



Single-Point Control and Automation

We provide advanced integrated IC/MS and LC/MS solutions with superior ease-of-use and modest price and space requirements. UltiMate 3000 System Wellness technology and automatic MS calibration allow continuous operation with minimal maintenance. The Dionex ion chromatography family automatically remove mobile phase ions for effort-free transition to MS detection.

- Thermo Scientific™ MSQ Plus™ mass spectrometer, the smallest and most sensitive single quadrupole on the market for LC and IC
- Self-cleaning ion source for low maintenance operation

Mass Spectrometry Instruments

- Chromeleon CDS software for single-point method setup, instrument control, and data management compatible with existing IC and LC methods
- The complete system includes the MSQ Plus mass spectrometer, PC data system, electrospray ionization (ESI) and atmospheric pressure chemical ionization (APCI) probe inlets, and vacuum system

Now, you no longer need two software packages to operate your LC/MS system. Chromeleon CDS software provides single-software method setup and instrument control; powerful UV, conductivity, and MS data analysis; and fully integrated reporting.



MSQ Plus Mass Spectrometer

Table of Contents

Introduction

Analytical Technologies

Artificial Colorants

Blue Dyes

Caramel

General Dye Method

Global Method

Red Dyes

Sudan Dyes

Yellow Dyes

Natural Colorants

Turmeric

Flavorants

Amadori Compounds

Capsaicin

Quinine

Vanilla

Miscellaneous

Caffeine

Choline

Glucosamine

Iodide

Polyphosphates

Preservatives

Benzoate

Nitrate and Nitrite

Sulfites

Soft Drinks

References



Analytical Technologies



Chromatography Data Systems

Tackle chromatography management challenges with the world's most complete chromatography software. Whether your needs are simple or complex or your scope is a single instrument, a global enterprise, or anything in between – the combination of Chromeleon CDS' scalable architecture and unparalleled ease-of use, makes your job easy and enjoyable with one Chromatography Data System for the entire lab.

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



The Fastest Way from Samples to Results

The 7.2 release of Chromeleon Chromatography Data System software is the first CDS that combines separation (GC/IC/LC) and Mass Spectrometry (MS) in an enterprise (client/server) environment. By extending Chromeleon 7.2 CDS beyond chromatography into MS, lab technicians can now streamline their chromatography and MS quantitation workflows with a single software package. MS support in Chromeleon 7.2 CDS is focused on routine and quantitative workflows, which provides access to rich quantitative data processing and automation capabilities — ultimately boosting your overall lab productivity and increasing the quality of your analytical results.



Chromeleon CDS Software

- Enjoy a modern, intuitive user interface designed around the principle of operational simplicity
- Streamline laboratory processes and eliminate errors with eWorkflows™, which enable anyone to perform a complete analysis perfectly with just a few clicks
- Access your instruments, data, and eWorkflows instantly in the Chromeleon Console
- Locate and collate results quickly and easily using powerful built-in database query features
- Interpret multiple chromatograms at a glance using MiniPlots
- Find everything you need to view, analyze, and report data in the Chromatography Studio
- Accelerate analyses and learn more from your data through dynamic, interactive displays
- Deliver customized reports using the built-in Excel® compatible spreadsheet

Excel is a registered trademark of Microsoft Corporation.

Table of Contents

Introduction

Analytical Technologies

Artificial Colorants

Blue Dyes

Caramel

General Dye Method

Global Method

Red Dyes

Sudan Dyes

Yellow Dyes

Natural Colorants

Turmeric

Flavorants

Amadori Compounds

Capsaicin

Quinine

Vanilla

Miscellaneous

Caffeine

Choline

Glucosamine

Iodide

Polyphosphates

Preservatives

Benzoate

Nitrate and Nitrite

Sulfites

Soft Drinks

References



Analytical Technologies



Process Analytical Systems

Thermo Scientific Dionex process analytical systems provide timely results by moving chromatography-based measurements on-line.

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

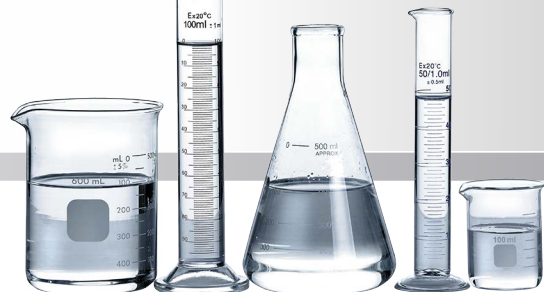
[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



Process Analytical Systems and Software

Improved Process Monitoring with On-line Chromatography IC and LC Systems

Information from the Thermo Scientific Dionex Integral process analyzer can help reduce process variability, improve efficiency, and reduce downtime. These systems provide comprehensive, precise, accurate information faster than is possible with laboratory-based results. From the lab to the factory floor, your plant's performance will benefit from the information provided by on-line LC.

- Characterize your samples completely with multicomponent analysis
- Reduce sample collection time and resources with automated multipoint sampling
- Improve your process control with more timely results
- See more analytes with unique detection capabilities
- The Thermo Scientific Integral Migration Path approach lets you choose the systems that best meets your needs



Integral process analyzer

Table of Contents

Introduction

Analytical Technologies

Artificial Colorants

Blue Dyes

Caramel

General Dye Method

Global Method

Red Dyes

Sudan Dyes

Yellow Dyes

Natural Colorants

Turmeric

Flavorants

Amadori Compounds

Capsaicin

Quinine

Vanilla

Miscellaneous

Caffeine

Choline

Glucosamine

Iodide

Polyphosphates

Preservatives

Benzoate

Nitrate and Nitrite

Sulfites

Soft Drinks

References



Analytical Technologies



Automated Sample Preparation

Solvent extractions that normally require labor-intensive steps are automated or performed in minutes, with reduced solvent consumption and reduced sample handling using the Thermo Scientific™ Dionex™ ASE™ Accelerated Solvent Extractor system or Thermo Scientific™ Dionex™ AutoTrace™ 280 Solid-Phase Extraction instrument.

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

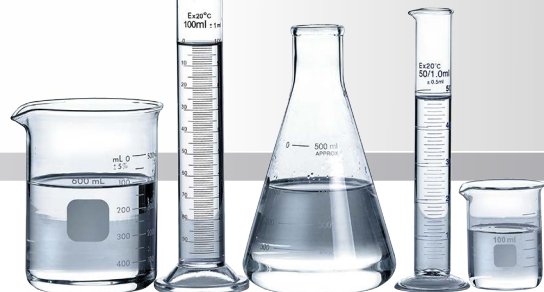
[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



Accelerated Solvent Extractor System

Complete Extractions in Less Time Using Less Solvent

Thermo Scientific Dionex ASE systems extract of solid and semisolid samples using common solvents at elevated temperature and pressure. The Dionex ASE 150 and 350 systems feature pH-hardened pathways with Dionium™ components to support extraction of acidic or alkaline matrices, and combine pretreatment, solvent extraction, and cleanup into one step. Dionium is zirconium that has undergone a proprietary

hardening process that makes it inert to chemical attack by acids and bases at elevated temperatures.

Dionex ASE systems are dramatically faster than Soxhlet, sonication, and other extraction methods, and require significantly less solvent and labor. Accelerated solvent extraction methods are accepted and established in the environmental, pharmaceutical, foods, polymers and consumer product industries. Accelerated solvent extraction methods are accepted and used by government agencies worldwide.



Dionex ASE 150/350 and Dionex AutoTrace 280 SPE instruments

Table of Contents

Introduction

Analytical Technologies

Artificial Colorants

Blue Dyes

Caramel

General Dye Method

Global Method

Red Dyes

Sudan Dyes

Yellow Dyes

Natural Colorants

Turmeric

Flavorants

Amadori Compounds

Capsaicin

Quinine

Vanilla

Miscellaneous

Caffeine

Choline

Glucosamine

Iodide

Polyphosphates

Preservatives

Benzoate

Nitrate and Nitrite

Sulfites

Soft Drinks

References



Additives Including Colorants, Flavorants, and Preservatives



Colorants

Food colorings (also called color additives), are any dyes, pigments, or substances that are used to impart color when they are added to food or drink. Colorants are used to enhance colors that occur naturally, correct natural variation in color, and to make food more attractive and appetizing. Colorants are also used to offset the loss of color due to processing and storage.



Artificial Colorants: Blue Dyes in Cereals

The FDA lists Food, Drugs, and Cosmetics (FD&C) certified color additives for use in foods in the United States. FD&C dyes are added to many popular breakfast cereal products to make them more visually appealing. This convenient and rapid assay uses ion-pair chromatography on a Thermo Scientific™ Acclaim™ C18 column to identify and quantify the dyes.

Pump: UltiMate LPG-3400
 Column: Acclaim 120 C18 3 μm, 3.0 × 75 mm
 Flow: 1.0 mL/min
 Temperature: UltiMate TCC-3100 at 30 °C
 Injection: UltiMate WPS-3000 SL sampler at 8 μL
 Mobile Phase: 677 g water, 0.97 g Na₂SO₄, 2.24 g KH₂PO₄, 3.20 g 55% Bu₄N OH (pH = 6.75), 250 g acetonitrile
 Detector: UltiMate VWD-3400, VIS at 427, 508, 625 nm, 2.5 Hz, 0.6 s time const.
 Sample Prep.: (1) Sort cereal by color
 (2) Grind to coarse powder
 (3) Extract 0.2 g of cereal with 5 mL mobile phase
 (4) Filter through 0.1 μm membrane
 Peaks: 1. FD&C Blue #2
 2. FD&C Blue #1

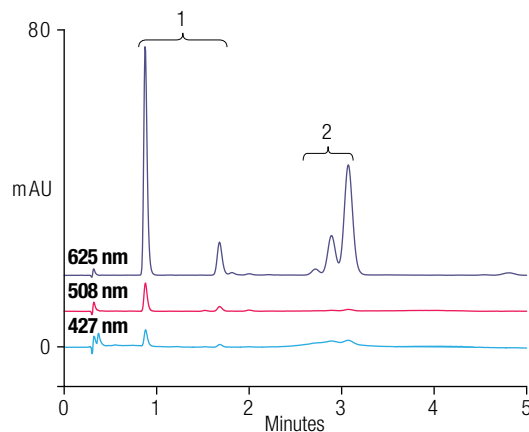
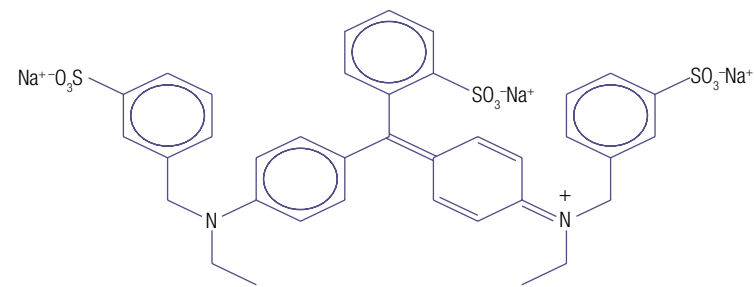
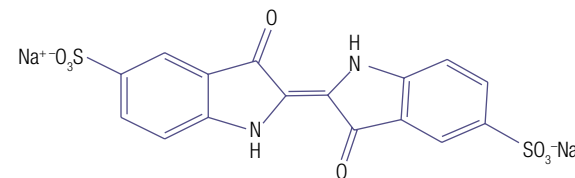


Figure 4-1. FD&C blue #1 and blue #2 in a blueberry-flavored cereal.



FDC Blue #1, Brilliant Blue, E133



FDC Blue #2, Indigocarmine, E132

Figure 4-2. FDC Blue dye structures.



Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

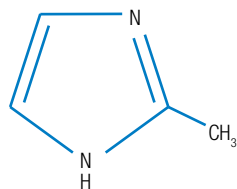
[References](#)



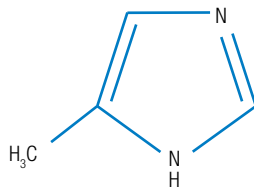
Table of Contents

- [Introduction](#)
- [Analytical Technologies](#)
- [Artificial Colorants](#)
 - [Blue Dyes](#)
 - [Caramel](#)
 - [General Dye Method](#)
 - [Global Method](#)
 - [Red Dyes](#)
 - [Sudan Dyes](#)
 - [Yellow Dyes](#)
- [Natural Colorants](#)
 - [Turmeric](#)
- [Flavorants](#)
 - [Amadori Compounds](#)
 - [Capsaicin](#)
 - [Quinine](#)
 - [Vanilla](#)
- [Miscellaneous](#)
 - [Caffeine](#)
 - [Choline](#)
 - [Glucosamine](#)
 - [Iodide](#)
 - [Polyphosphates](#)
- [Preservatives](#)
 - [Benzoate](#)
 - [Nitrate and Nitrite](#)
 - [Sulfites](#)
 - [Soft Drinks](#)
- [References](#)

The 2- and 4-methylimidazoles (2-MI and 4-MI) are undesired byproducts from the manufacturing of the caramel color ingredient used to darken food products such as carbonated beverages and soy sauces. Studies from the National Toxicology Program (NTP) and other researchers have concluded that there is clear evidence of both chemicals' carcinogenicity. California's Office of Environmental Health Hazard Assessment (OEHHA) listed 4-MI as a carcinogen in January 2011 with a proposed No Significant Risk Level (NSRL) of 16 µg per person per day. In February 2011, a group of scientists from the Center for Science in the Public Interests filed a petition with the U.S. FDA to bar the use of caramel coloring containing 2-MI and 4-MI, which is used for "purely cosmetic purposes." On the other hand, the European Food Safety Authority concluded that exposure to 4-MI from caramel coloring is not of concern. Although the issue is controversial, it seems prudent to take precautions, especially for sensitive populations susceptible to excessive exposure. Therefore, a quantitative analytical measurement method for both compounds is needed for exposure assessment.



2-methylimidazole
 Chemical Formula: C₄H₆N₂
 Exact Mass: 82.05
 Molecular Weight: 82.10



4-methylimidazole
 Chemical Formula: C₄H₆N₂
 Exact Mass: 82.05
 Molecular Weight: 82.10

Figure 4-3. Chemical structures for caramel colorants.

Artificial Colorants: Caramel

Column: Dionex IonPac CS19 (2 × 250 mm)
 Flow: 0.25 mL/min
 Temperature: 40 °C
 Injection Volume: 5 µL
 Eluent Source: Dionex EGC II MSA cartridge
 Eluent: Isocratic 1.7 mM MSA, gradient to 7 mM MSA
 From 9 to 13 min, isocratic to 25 min, back to
 1.7 mM MSA at 25.1 min
 Detection: Suppressed Conductivity, Thermo Scientific™ Dionex™ CSRS 300,
 AutoSuppression, recycle mode

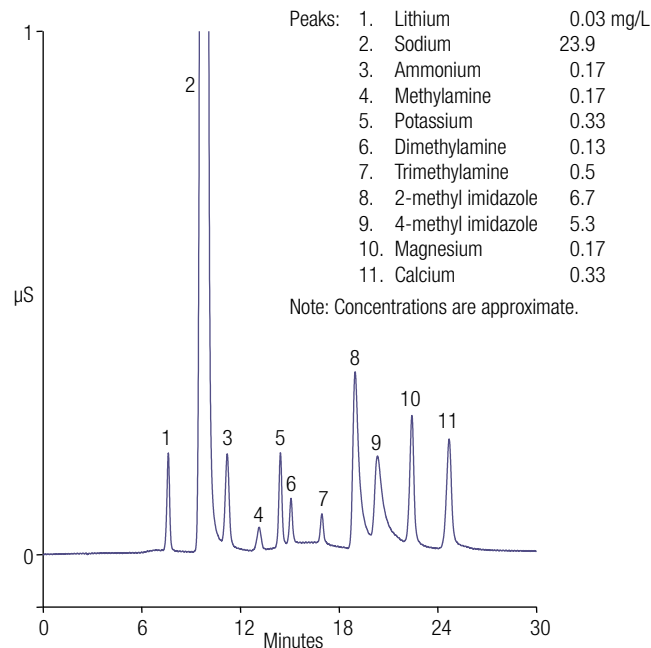


Figure 4-4. Gradient separation of six common cations, methylamines and imidazoles using the Dionex IonPac CS19 2 mm column.



Artificial Colorants: General Dye Method

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

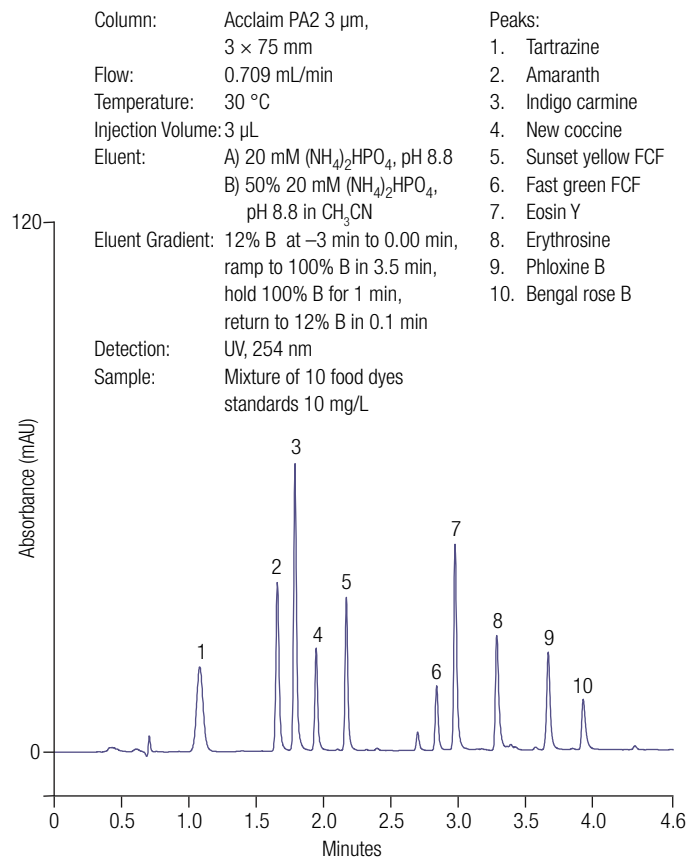


Figure 4-5. Chromatogram of the standard mixture of 10 dyes.

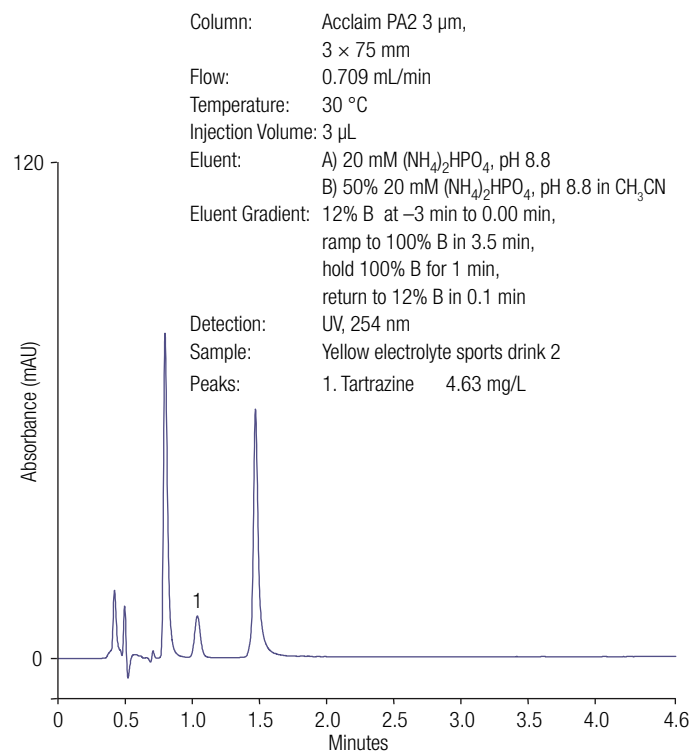


Figure 4-6. Chromatogram of an electrolyte sports drink.



Artificial Colorants: General Dye Method

Table of Contents

[Introduction](#)
[Analytical Technologies](#)

Artificial Colorants

[Blue Dyes](#)
[Caramel](#)
[General Dye Method](#)
[Global Method](#)
[Red Dyes](#)
[Sudan Dyes](#)
[Yellow Dyes](#)

Natural Colorants

[Turmeric](#)

Flavorants

[Amadori Compounds](#)
[Capsaicin](#)
[Quinine](#)
[Vanilla](#)

Miscellaneous

[Caffeine](#)
[Choline](#)
[Glucosamine](#)
[Iodide](#)
[Polyphosphates](#)

Preservatives

[Benzoate](#)
[Nitrate and Nitrite](#)
[Sulfites](#)
[Soft Drinks](#)

References

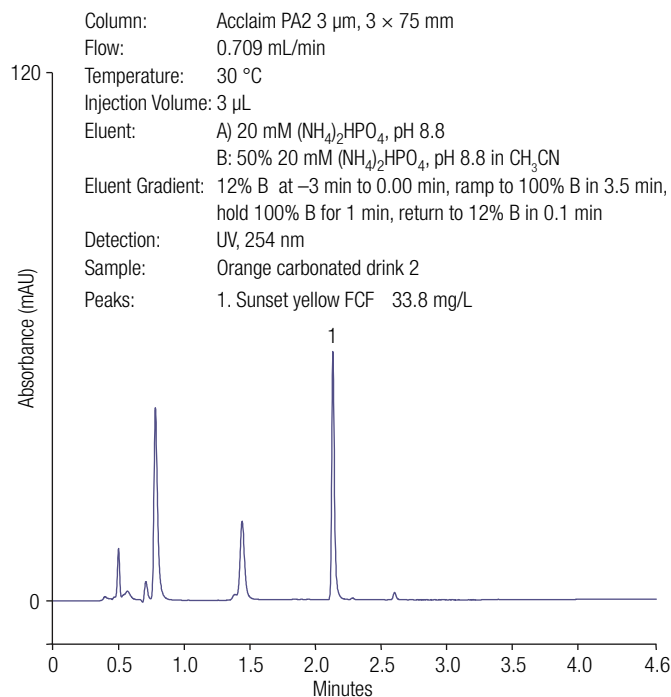


Figure 4-7. Chromatogram of an orange carbonated drink.

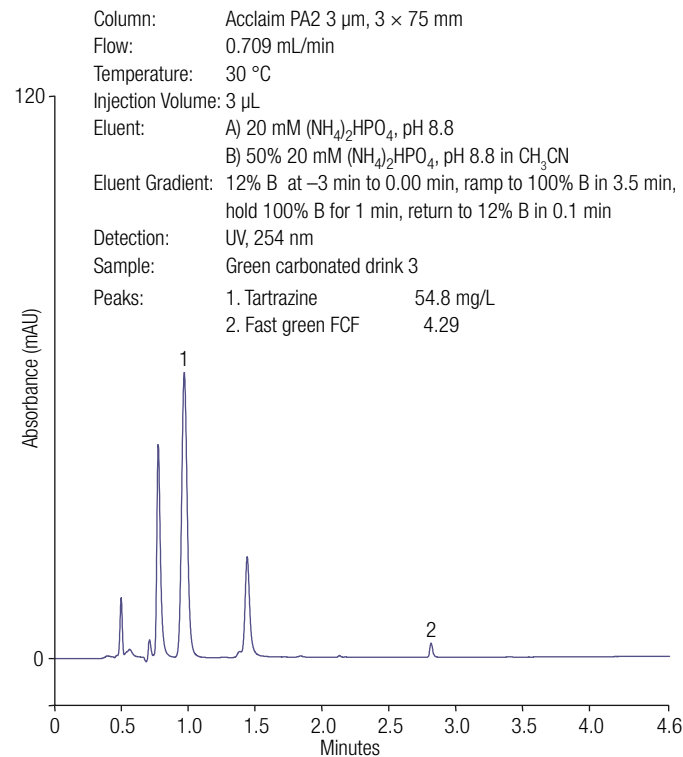


Figure 4-8. Chromatogram of a green carbonated drink.





Artificial Colorants: General Dye Method

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

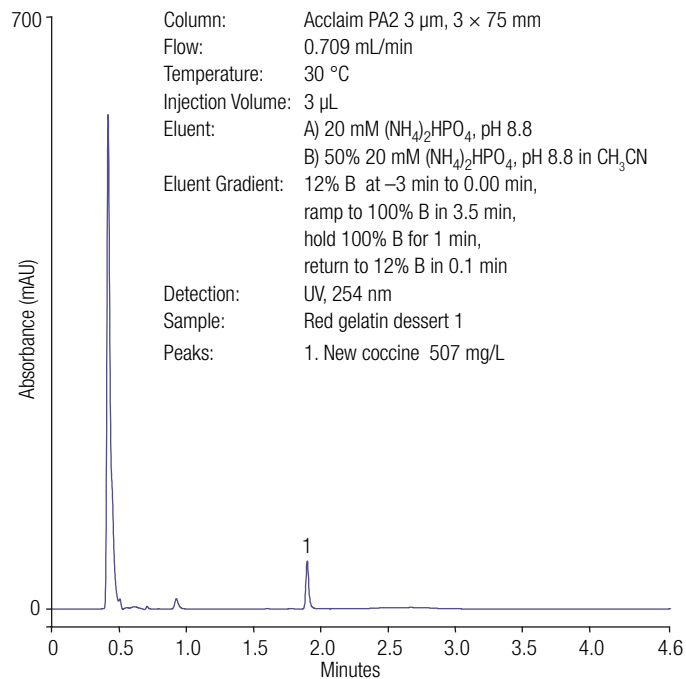


Figure 4-9. Chromatogram of a red gelatin dessert sample.





Table of Contents

- [Introduction](#)
- [Analytical Technologies](#)
- [Artificial Colorants](#)
 - [Blue Dyes](#)
 - [Caramel](#)
 - [General Dye Method](#)
 - [Global Method](#)
 - [Red Dyes](#)
 - [Sudan Dyes](#)
 - [Yellow Dyes](#)
- [Natural Colorants](#)
 - [Turmeric](#)
- [Flavorants](#)
 - [Amadori Compounds](#)
 - [Capsaicin](#)
 - [Quinine](#)
 - [Vanilla](#)
- [Miscellaneous](#)
 - [Caffeine](#)
 - [Choline](#)
 - [Glucosamine](#)
 - [Iodide](#)
 - [Polyphosphates](#)
- [Preservatives](#)
 - [Benzoate](#)
 - [Nitrate and Nitrite](#)
 - [Sulfites](#)
 - [Soft Drinks](#)
- [References](#)

Artificial Colorants: Global Method for Permitted and Banned Food Dyes

Colors are food additives that are regulated in the EU, USA, and China. Only approved food colors can be used and others are banned such as some azo compounds, because of their potential carcinogenicity. The so-called Sudan colors are usually used for coloring oils or as additives in shoe polish, but have been used illegally in foods. They are classified as class 3 carcinogens by the International Agency for Research on Cancer (IARC). Despite considerable publicity these banned colors are still found in foodstuffs, e.g. chili and chili containing products, according to recent RASFF alerts. Directive 2005/402/EC introduced emergency measures that has required certification of each product exported to the EU to be certified to be free of Sudan colors.

Method 52212 enables the measurement of both regulated water-soluble and banned fat-soluble colors and can be applied to liquid, semi-liquid and solid food samples. The method combines cloud point extraction and

LC-MS to quantify banned azo colors (Sudan I-IV, Sudan red G, Sudan red B, Sudan red 7B, Sudan black B, metanil yellow and rhodamine B), at an LOD of 0.5 mg/kg. Simultaneously, regulated food colors (sunset yellow, allura red, tartrazine and erythrosine) can be analyzed in liquid, semi-liquid and solid food and beverages.

Table 4-1. Limits of detection and limits of quantification (LODs and LOQs) of the LC-MS method for different matrices

Analyte	Liquid Matrix (wine)		Semi-liquid Matrix (sauce)		Solid Matrix (chili)	
	LOD [mg/kg]	LOQ [mg/kg]	LOD [mg/kg]	LOQ [mg/kg]	LOD [mg/kg]	LOQ [mg/kg]
Allura Red	47.0	140.0	5.0	15.0	13.0	38.0
Sunset Yellow	50.0	150.0	12.0	35.0	8.3	25.0
Tartrazine	**	**	40.0	120.0	**	**
Erythrosine	30.0	90.0	30.0	90.0	30.0	90.0
Metanil Yellow	*	*	0.5	1.5	0.3	0.9
Rhodamine B	*	*	0.03	0.1	0.03	0.1
Sudan I	*	*	0.5	1.5	0.4	1.3
Sudan II	*	*	0.3	0.9	0.3	0.9
Sudan III	*	*	0.2	0.6	0.2	0.6*
Sudan IV	*	*	0.1	0.3	0.1	0.3
Sudan Red G	*	*	0.5	1.5	0.4	1.2
Sudan Red B	*	*	0.5	1.5	0.5	1.5
Sudan Red 7B	*	*	0.5	1.5	0.5	1.5
Sudan Black B 1 st	*	*	0.03	0.1	0.03	0.1
Sudan Black B 2 nd	*	*	0.3	0.9	0.3	0.9

* Colors were not spiked because not relevant to food
 ** Compound not applicable for this matrix

Column: Thermo Scientific™ Hypersil GOLD™, 150 x 2.1 mm, 1.9 μm
 Tray Temperature Control: 22 °C
 Column Oven Temperature: 22 °C
 Injection Volume: 10 μL
 Injection Mode: Partial loop
 Mobile Phase: A; water (0.1 % formic acid); B: acetonitrile (0.1% formic acid)

Gradient:	Time	A[%]	B[%]
	0	99	1
	5	99	1
	10	30	70
	15	25	75
	35	25	75
	35.01	99	1
	42	99	1



Table of Contents

- [Introduction](#)
- [Analytical Technologies](#)
- [Artificial Colorants](#)
 - [Blue Dyes](#)
 - [Caramel](#)
 - [General Dye Method](#)
 - [Global Method](#)
 - [Red Dyes](#)
 - [Sudan Dyes](#)
 - [Yellow Dyes](#)
- [Natural Colorants](#)
 - [Turmeric](#)
- [Flavorants](#)
 - [Amadori Compounds](#)
 - [Capsaicin](#)
 - [Quinine](#)
 - [Vanilla](#)
- [Miscellaneous](#)
 - [Caffeine](#)
 - [Choline](#)
 - [Glucosamine](#)
 - [Iodide](#)
 - [Polyphosphates](#)
- [Preservatives](#)
 - [Benzoate](#)
 - [Nitrate and Nitrite](#)
 - [Sulfites](#)
 - [Soft Drinks](#)
- [References](#)

Artificial Colorants: Global Method for Permitted and Banned Food Dyes

MS analysis is carried out using a MSQ Plus LC single quadrupole mass spectrometer. Data acquisition and processing was performed using Thermo Scientific™ Xcalibur™ software. The MS conditions were as follows:

Ionization: Electrospray
Polarity: Positive negative switching
Probe Temperature: 450 °C
Cone Voltage: 150 V
Scan Mode: Multi ion monitoring (MIM)

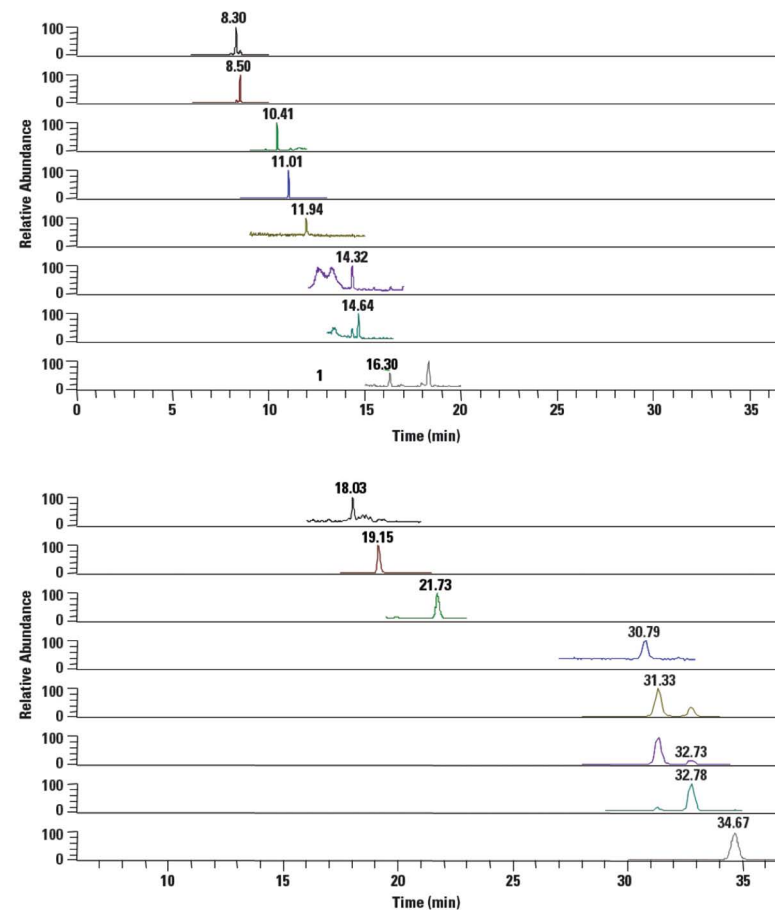


Figure 4-10. LC-MS chromatogram: separation and detection of all colors in sauce matrix.

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



Artificial Colorants: Red Dyes in Cereals

Pump: UltiMate LPG-3400
Column: Acclaim 120 C18 3 μm , 3.0 \times 75 mm
Flow: 1.0 mL/min
Temperature: UltiMate TCC-3100 at 30 $^{\circ}\text{C}$
Injection: UltiMate WPS-3000 SL sampler at 8 μL
Mobile Phase: 677 g water, 0.97 g Na_2SO_4 , 2.24 g KH_2PO_4 , 3.20 g 55% Bu_4NOH (pH = 6.75), 250 g acetonitrile
Detector: UltiMate VWD-3400, VIS at 427, 508, 625 nm, 2.5 Hz, 0.6 s time const.
Sample Prep.: (1) Sort cereal by color
(2) Grind to coarse powder
(3) Extract 0.2 g of cereal with 5 mL mobile phase
(4) Filter through 0.1 μm membrane
Peaks: 1. FD&C Red #40

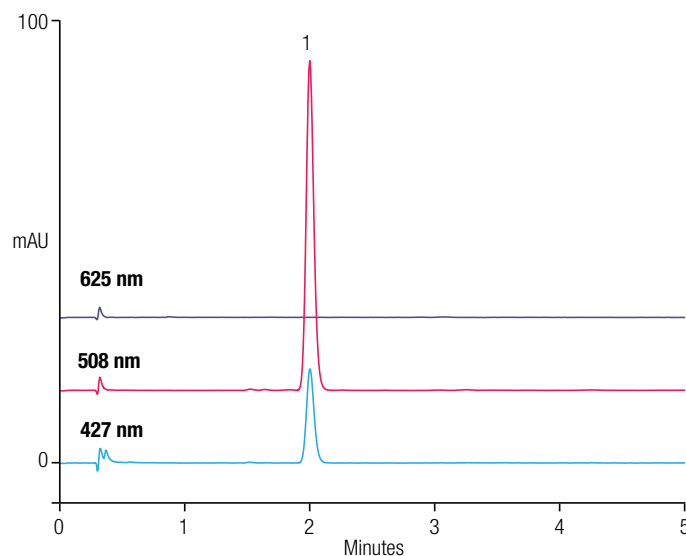
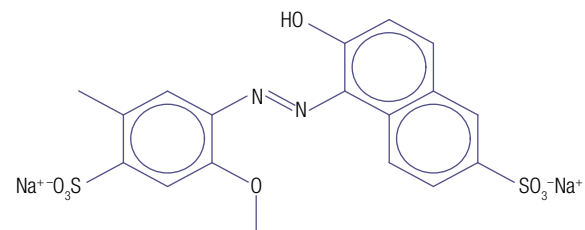


Figure 4-11. FD&C red #40 used in a cherry-flavored cereal.



FD&C Alura Red AC, E129

Figure 4-12. FD&C Red #40 dye structure.





Artificial Colorants: Sudan Dyes in Chili Oil

Table of Contents

- [Introduction](#)
- [Analytical Technologies](#)
- [Artificial Colorants](#)
 - [Blue Dyes](#)
 - [Caramel](#)
 - [General Dye Method](#)
 - [Global Method](#)
 - [Red Dyes](#)
 - [Sudan Dyes](#)
 - [Yellow Dyes](#)
- [Natural Colorants](#)
 - [Turmeric](#)
- [Flavorants](#)
 - [Amadori Compounds](#)
 - [Capsaicin](#)
 - [Quinine](#)
 - [Vanilla](#)
- [Miscellaneous](#)
 - [Caffeine](#)
 - [Choline](#)
 - [Glucosamine](#)
 - [Iodide](#)
 - [Polyphosphates](#)
- [Preservatives](#)
 - [Benzoate](#)
 - [Nitrate and Nitrite](#)
 - [Sulfites](#)
 - [Soft Drinks](#)
- [References](#)

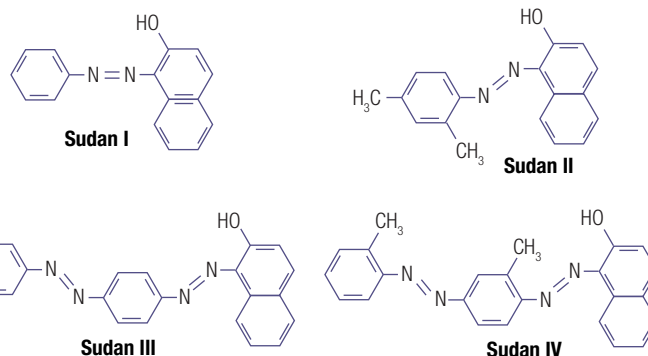


Figure 4-13. Structures of Sudan dyes I, II, III, and IV.

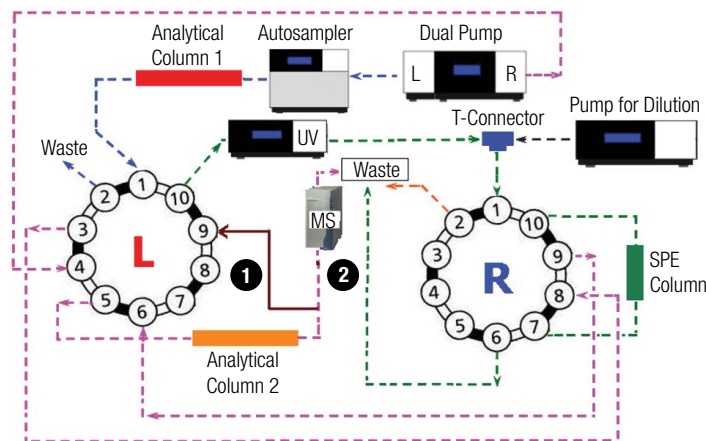


Figure 4-14. Flow schematics for a two-dimensional HPLC with online SPE intercolumn trapping system. Flow path 1 is configured for method development using a UV-vis detector; flow path 2 is configured for the determination of Sudan dyes in chili oil using an MS detector.





Artificial Colorants: Sudan Dyes in Chili Oil

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Column: Analytical column 1 (for the first dimension)
Acclaim PA2, 3 μm , 3.0 \times 150 mm (P/N 063705)
On-line SPE
Acclaim 120 C18, 5 μm , Guard Cartridge, 4.6 \times 10 mm (P/N 069580)
Analytical column 2 (for the second dimension)
Acclaim RSLC 120 C18, 2.2 μm , 2.1 \times 100 mm (P/N 068982)

Flow: For analytical column 1 (the first dimension), 0.6 mL/min
For on-line SPE, 1.0 mL/min
For analytical column 2 (the second dimension), 0.3 mL/min

Column Temp.: 30 $^{\circ}\text{C}$

Injection Volume: 20 μL on the analytical column 1

Mobile Phase: For analytical column 1 (the first dimension)
A: Water
B: CH_3CN
C: $\text{CH}_3\text{OH}/\text{THF}$, 1:1 (v/v)
In gradient: -3.5 ~ -0.5 min, B 50%, C 20%; 5.5 min, B 50%, C 50%; 11.0 ~ 14.0 min, B 20%, C 80%; 14.1 ~ 23 min, B 50%, C 20%

For on-line SPE
0.1% FA in water, isocratic
For analytical column 2 (the second dimension):
A: Water
B: CH_3CN
C: 0.1% FA in CH_3CN
In gradient: -3.5 ~ 7.0 min, B 50%; 16.0 ~ 22.4 min, B 90%; 22.5 ~ 23.0 min, B 50%.
A and C maintain 40% and 10%, respectively.

UV-vis Detection: Absorbance at 500 nm

Chromatograms: (A) Sudan dye standards (2 mg/L for each Sudan dye)
(B) Chili oil sample, spiked with 6 mg/L of each Sudan dye standard

Peaks:

1. Sudan dye I
2. Sudan dye II
3. Sudan dye III
4. Sudan dye IV

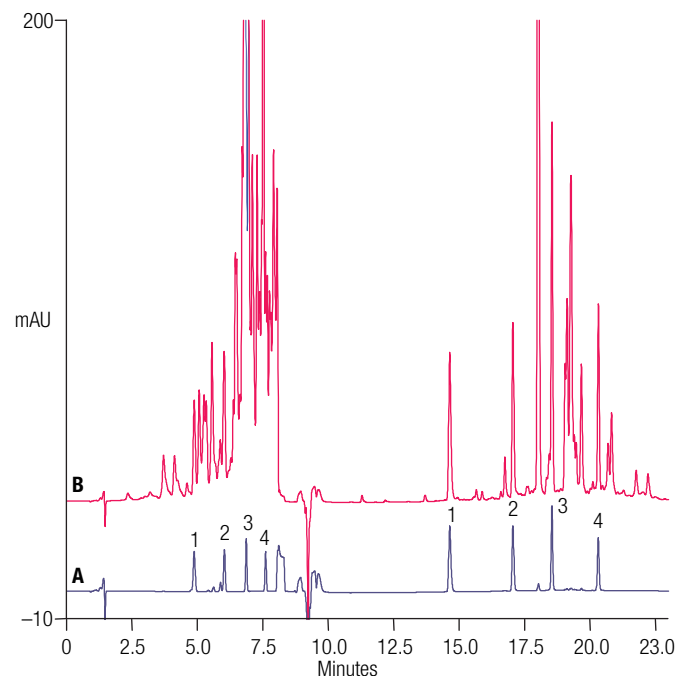


Figure 4-15. Chromatograms of (A) Sudan dye standards (2 mg/L for each Sudan dye), and (B) chili oil sample, spiked with 6 mg/L of each dye standard using the two-dimensional HPLC with on-line SPE intercolumn trapping system with UV-vis detection. The second dimension separation starts at 10 min.



Artificial Colorants: Sudan Dyes in Chili Oil

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Column: Analytical column 1 (for the first dimension)
Acclaim PA2, 3 μm , 3.0 \times 150 mm (P/N 063705)
On-line SPE
Acclaim 120 C18, 5 μm , Guard Cartridge, 4.6 \times 10 mm (P/N 069580)
Analytical column 2 (for the second dimension)
Acclaim RSLC 120 C18, 2.2 μm , 2.1 \times 100 mm (P/N 068982)

Flow: For analytical column 1 (the first dimension), 0.6 mL/min
For on-line SPE, 1.0 mL/min
For analytical column 2 (the second dimension), 0.3 mL/min

Column Temp.: 30 $^{\circ}\text{C}$
Injection Volume: 20 μL on the analytical column 1

Mobile Phase: For analytical column 1 (the first dimension)
A: Water
B: CH_3CN
C: $\text{CH}_3\text{OH}/\text{THF}$, 1:1 (v/v)
In gradient: -3.5 ~ -0.5 min, B 50%, C 20%; 5.5 min, B 50%, C 50%; 11.0 ~ 14.0 min, B 20%, C 80%; 14.1 ~ 23 min, B 50%, C 20%
For on-line SPE
0.1% FA in water, isocratic
For analytical column 2 (the second dimension):
A: Water
B: CH_3CN
C: 0.1% FA in CH_3CN
In gradient: -3.5 ~ 7.0 min, B 50%; 16.0 ~ 22.4 min, B 90%; 22.5 ~ 23.0 min, B 50%.
A and C maintain 40% and 10%, respectively.

MS Detection: ESI mode, positive scan, probe temp at 450 $^{\circ}\text{C}$,
needle voltage at 4000 V, SIM at 249 m/z for Sudan dye I,
277 m/z for Sudan dye II, 353 m/z for Sudan dye III,
and 381 m/z for Sudan dye IV;
cone voltage at 35 V for Sudan dye I and II,
50 V for Sudan dye III and IV; nebulizer gas, Nitrogen at 75 psi

Chromatograms: (A) blank
(B) Sudan dye standards (5 $\mu\text{g}/\text{L}$ for Sudan dyes II and III, and 15 $\mu\text{g}/\text{L}$ for Sudan dyes I and IV)
(C) chili oil sample
(D) Chili oil sample, spiked with 10 $\mu\text{g}/\text{L}$ of Sudan dye II and III, and 30 $\mu\text{g}/\text{L}$ of Sudan dye I and IV standards

Peaks:
1. Sudan dye I
2. Sudan dye II
3. Sudan dye III
4. Sudan dye IV

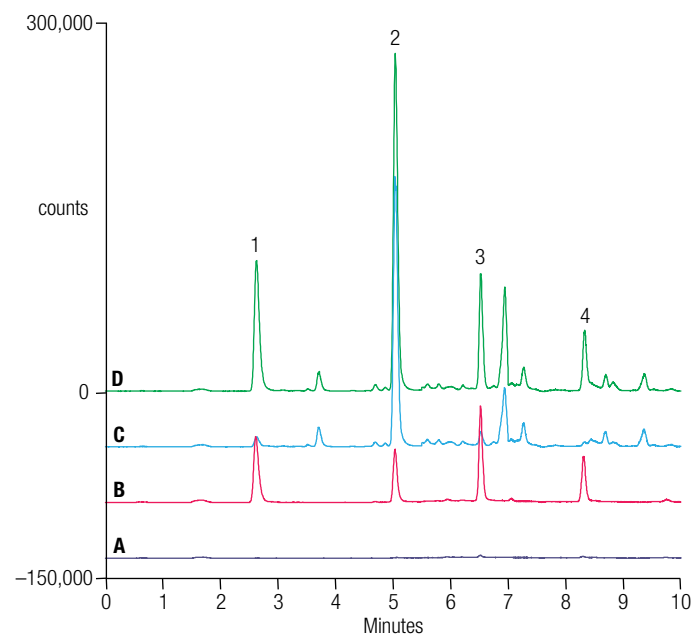


Figure 4-16. MS TIC chromatograms of (A) blank (CH_3CN), (B) mixed Sudan dye standards (5 $\mu\text{g}/\text{L}$ for Sudan dye II and III, and 15 $\mu\text{g}/\text{L}$ for Sudan dye I and IV), (C) chili oil sample, and (D) the same sample spiked with 10 $\mu\text{g}/\text{L}$ of Sudan dyes II and III, and 30 $\mu\text{g}/\text{L}$ of Sudan dyes I and IV using the two-dimensional HPLC with on-line SPE intercolumn trapping system with MS detection.



Artificial Colorants: Sudan Dyes in Curry Paste

Table of Contents

- [Introduction](#)
- [Analytical Technologies](#)
- [Artificial Colorants](#)
 - [Blue Dyes](#)
 - [Caramel](#)
 - [General Dye Method](#)
 - [Global Method](#)
 - [Red Dyes](#)
 - [Sudan Dyes](#)
 - [Yellow Dyes](#)
- [Natural Colorants](#)
 - [Turmeric](#)
- [Flavorants](#)
 - [Amadori Compounds](#)
 - [Capsaicin](#)
 - [Quinine](#)
 - [Vanilla](#)
- [Miscellaneous](#)
 - [Caffeine](#)
 - [Choline](#)
 - [Glucosamine](#)
 - [Iodide](#)
 - [Polyphosphates](#)
- [Preservatives](#)
 - [Benzoate](#)
 - [Nitrate and Nitrite](#)
 - [Sulfites](#)
 - [Soft Drinks](#)
- [References](#)

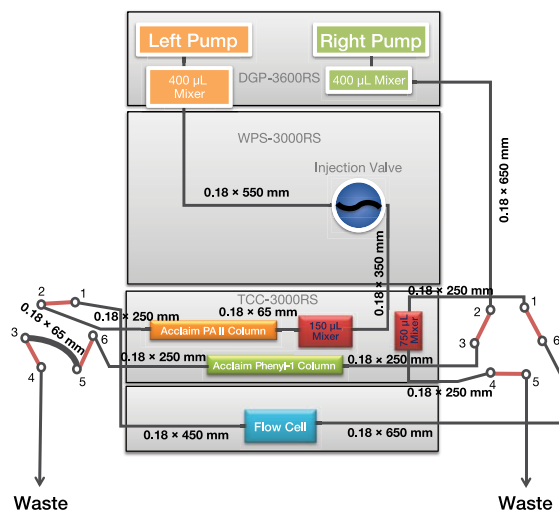


Figure 4-17. System configuration.

First Dimension

Column: Acclaim PA2 (3 μ m, 4.6 \times 150 mm)
 Mobile Phase: A: Water B: Acetonitrile C: 2-Propanol
 Gradient: See Table 2
 Flow: See Table 2
 Injection Volume: 300 μ L
 Temperature: 30 $^{\circ}$ C
 Detection: UV, 478 nm

Second Dimension

Column: Acclaim Phenyl-1 (3 μ m, 4.6 \times 150 mm)
 Mobile Phase: A: Water B: Acetonitrile C: Methanol
 Gradient: See Table 2
 Flow Rate: See Table 2
 Temperature: 30 $^{\circ}$ C
 Detection: UV, 478 nm

Samples: 1. Sudan dyes standard mixture with 150 μ L mixer
 2. Sudan dyes standard mixture without 150 μ L mixer

Table 4-2. Gradient program and valve switching.

First Dimension (Left Pump)					Valve Switching			Second Dimension (Right Pump)				
Time (min)	Flow Rate (mL/min)	% A (Water)	% B (CH ₃ CN)	% C (C ₃ H ₇ OH)	Time (min)	Right Valve Position	Left Valve Position	Time (min)	Flow Rate (mL/min)	% A (Water)	% B (CH ₃ CN)	% C (CH ₃ OH)
0.00	1.0	80	20	0	0.00	1-2	1-2	0.00	1.0	85	0	15
2.00	1.0	5	95	0	6.773	6-1	1-2	13.0	1.0	85	0	15
13.5	1.0	5	95	0	7.042	1-2	1-2	14.0	1.0	0	0	100
14.00	0.7	0	5	95	7.947	6-1	1-2	15.0	1.0	0	0	100
19.00	0.7	0	5	95	8.164	1-2	1-2	16.0	1.0	0	45	55
20.0	0.7	5	95	0	9.583	6-1	1-2	25.0	1.0	0	45	55
22.0	0.7	5	95	0	9.811	1-2	1-2	—	—	—	—	—
22.5	1.0	80	20	0	11.990	6-1	1-2	—	—	—	—	—
25.0	1.0	80	20	0	12.290	1-2	1-2	—	—	—	—	—
—	—	—	—	—	13.500	6-1	6-1	—	—	—	—	—



Artificial Colorants: Sudan Dyes in Curry Paste

Table of Contents

[Introduction](#)
[Analytical Technologies](#)

Artificial Colorants

[Blue Dyes](#)
[Caramel](#)
[General Dye Method](#)
[Global Method](#)
[Red Dyes](#)
[Sudan Dyes](#)
[Yellow Dyes](#)

Natural Colorants

[Turmeric](#)

Flavorants

[Amadori Compounds](#)
[Capsaicin](#)
[Quinine](#)
[Vanilla](#)

Miscellaneous

[Caffeine](#)
[Choline](#)
[Glucosamine](#)
[Iodide](#)
[Polyphosphates](#)

Preservatives

[Benzoate](#)
[Nitrate and Nitrite](#)
[Sulfites](#)
[Soft Drinks](#)

References

First Dimension

Column: Acclaim PA2
 (3 μ m, 4.6 \times 150 mm)
 Flow: See Table 2
 Temperature: 30 $^{\circ}$ C
 Injection Volume: 300 μ L
 Mobile Phase: A: Water
 B: Acetonitrile
 C: 2-Propanol
 Gradient: See Table 2
 Detection: UV, 478 nm

Second Dimension

Column: Acclaim Phenyl-1
 (3 μ m, 4.6 \times 150 mm)
 Flow: See Table 2
 Temperature: 30 $^{\circ}$ C
 Mobile Phase: A: Water
 B: Acetonitrile
 C: Methanol
 Gradient: See Table 2
 Detection: UV, 478 nm
 Samples: Sudan I, II, III, and IV
 working standards

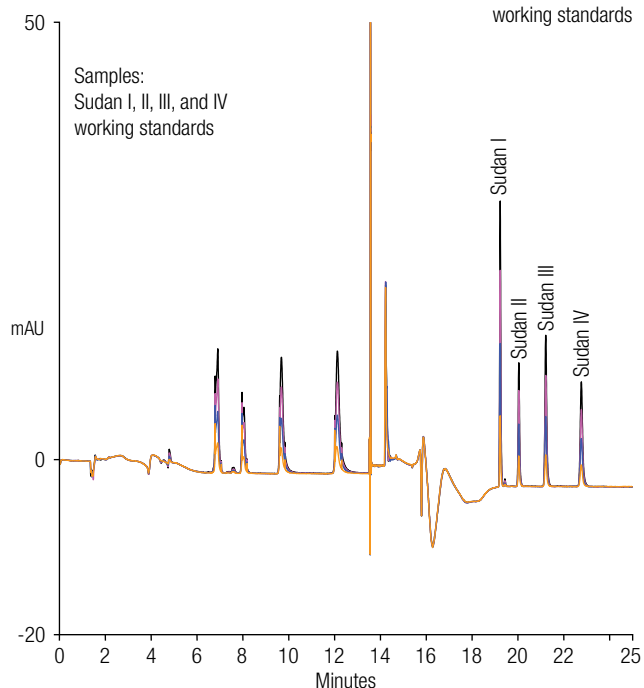


Figure 4-18. Overlay of chromatograms of working standards.

First Dimension

Column: Acclaim PA2
 (3 μ m, 4.6 \times 150 mm)
 Flow: See Table 2
 Temperature: 30 $^{\circ}$ C
 Injection Volume: 300 μ L
 Mobile Phase: A: Water
 B: Acetonitrile
 C: 2-Propanol
 Gradient: See Table 2
 Detection: UV, 478 nm

Second Dimension

Column: Acclaim Phenyl-1
 (3 μ m, 4.6 \times 150 mm)
 Flow: See Table 2
 Temperature: 30 $^{\circ}$ C
 Mobile Phase: A: Water
 B: Acetonitrile
 C: Methanol
 Gradient: See Table 2
 Detection: UV, 478 nm
 Samples: 1. Fresh curry paste
 2. Spiked fresh curry paste

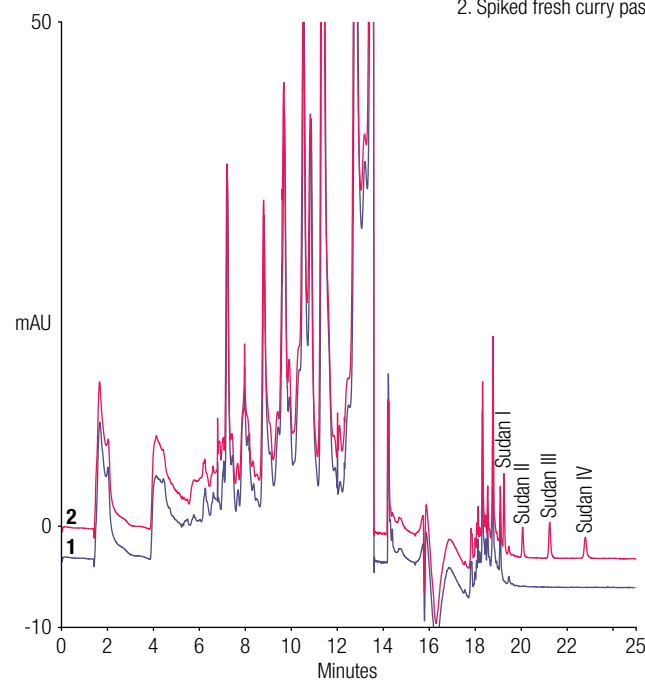


Figure 4-19. Overlay chromatograms of fresh curry paste and spiked fresh curry paste samples.



Artificial Colorants: Sudan Dyes in Paprika

Sudan dyes are a class of synthetic dyes that are mainly used for industrial applications such as the coloring of plastic. These dyes are banned as a food-coloring agent because they are classified as carcinogens. For economical reasons, however, Sudan dyes are sometimes illegally used to color food to improve its appearance. Therefore, methods are needed to determine if food products have been adulterated with these dyes.

There have been cases of adulteration of capsicum products with Sudan Orange and other illegal dyes that resulted in large-scale product recalls. The chromatography shown here clearly resolves the Sudan dyes while compressing the natural pigments into a well-removed band. The example shown here has >85% recovery for all dyes.



Column: Acclaim 120 C18, 5 μ m, 4.6 \times 150 mm
Flow: 1.25 mL/min
Temperature: 30 $^{\circ}$ C
Injection: Summit ASI-100, 15 μ L
Pump: Summit P680 DGP
Mobile Phase: (A) Water
(B) 1:1 tetrahydrofuran:methanol
Gradient: Times 0 7.0 7.1 13
%A 20 8 0 0
%B 80 92 100 100
Detector: Diode array, 500 nm, and spectra 200–595 nm
Samples: (A) Control paprika without artificial coloring
(B) Paprika spiked with approx. 25 mg/kg of each dye
Peaks: 1. Sudan I
2. Sudan II
3. Sudan III
4. Sudan IV
5. Oil Red O (2 peaks)
(Natural pigments 8–12 min)

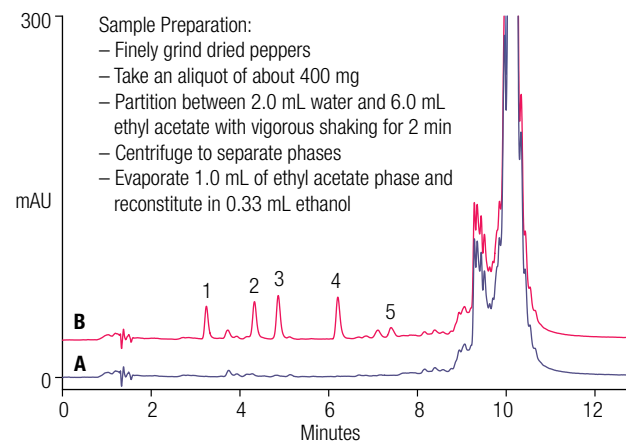


Figure 4-20. Overlay chromatograms of control and paprika spiked with sudan dyes.

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



Artificial Colorants: Yellow Dyes in Cereals

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Pump: UltiMate LPG-3400
Column: Acclaim 120 C18 3 μm , 3.0 \times 75 mm
Flow: 1.0 mL/min
Temperature: UltiMate TCC-3100 at 30 $^{\circ}\text{C}$
Injection: UltiMate WPS-3000 SL sampler at 8 μL
Mobile Phase: 677 g water, 0.97 g Na_2SO_4 , 2.24 g KH_2PO_4 ,
3.20 g 55% Bu_4NOH (pH = 6.75), 250 g acetonitrile
Detector: UltiMate VWD-3400, VIS at 427, 508,
625 nm, 2.5 Hz, 0.6 s time const.
Sample Prep.: (1) Sort cereal by color
(2) Grind to coarse powder
(3) Extract 0.2 g of cereal with 5 mL mobile phase
(4) Filter through 0.1 μm membrane
Peaks: 1. FD&C Yellow #6

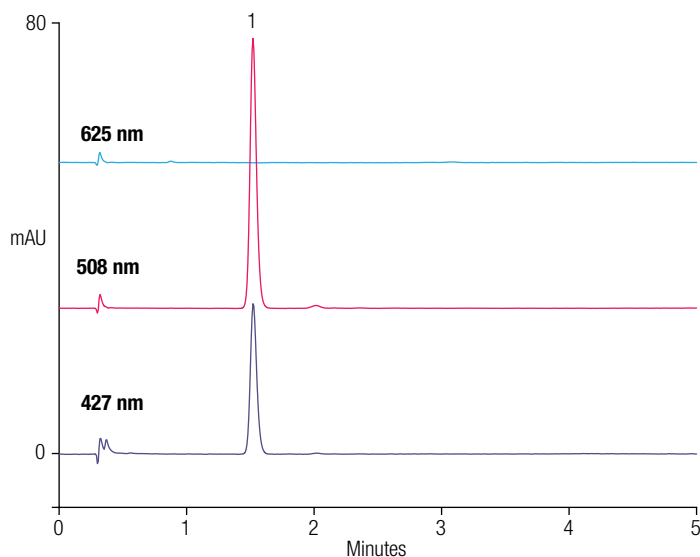


Figure 4-21. FD&C yellow #6 used in an orange-flavored cereal.

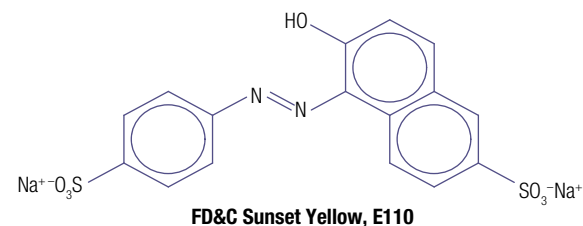


Figure 4-22. FD&C yellow #6 dye structure.





Natural Colorants: Turmeric by UV

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Pigments in Turmeric

Turmeric (*Curcuma longa*) is used as a spice, a natural yellow food color, an herbal medicine, and even a colorimetric reagent. This versatile plant has three major yellow pigments, but they are not well resolved on C18 columns. Using the alternate selectivity of the Acclaim PolarAdvantage II columns gives excellent separation.

Columns:	Acclaim RSLC 120 C18 Acclaim RSLC PA2	Mobile Phase: A: 15 mM H ₃ PO ₄ Mobile Phase: B: Methanol
Dimensions:	2 μm, 2.1 × 100 mm	Isocratic: C18: 70% B (v/v) PA2: 80% B (v/v)
Flow:	0.41 mL/min	Detector: UV, 428 nm
Temperature:	30 °C	Sample: Turmeric extract
		Peaks: 1. Curcumin 2. Demethoxycurcumin 3. Bis-demethoxycurcumin

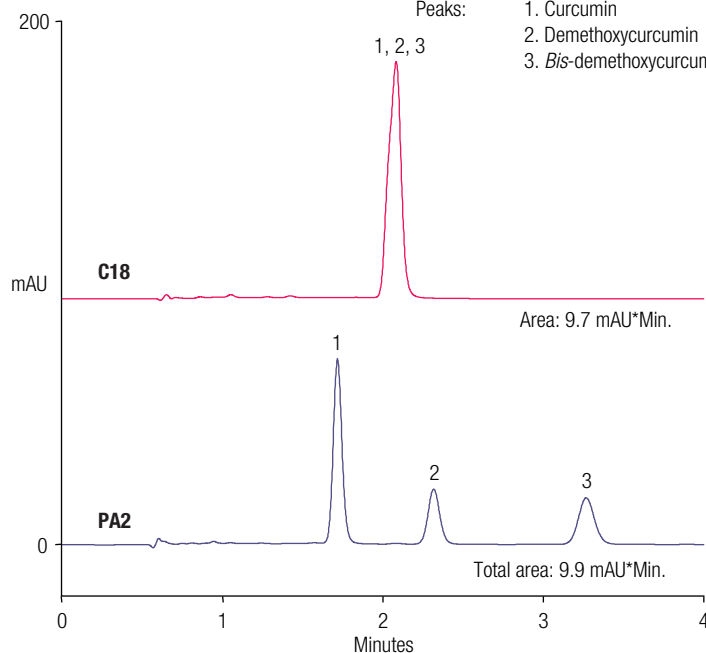


Figure 4-23. Improved resolution of curcuminoids using an Acclaim PA2 column.

Did You Know?

- Turmeric is a member of the ginger family.
- It is used to color foods yellow (designated E100) and as a fabric dye, although it is not very light stable.
- It is used in Ayurvedic medicine to treat stomach and liver ailments and is being investigated in Western medicine as a treatment to Alzheimer's, diabetes, and other diseases

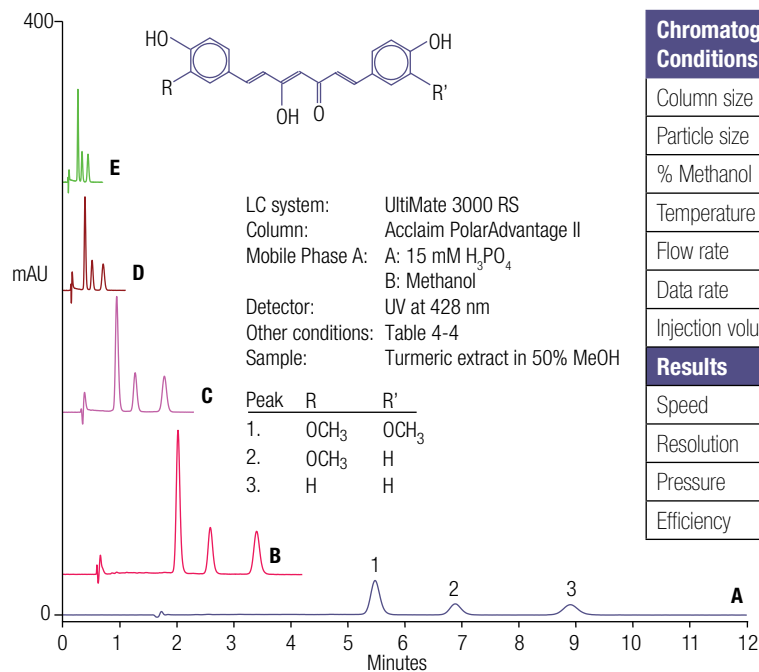




Table of Contents

- [Introduction](#)
- [Analytical Technologies](#)
- [Artificial Colorants](#)
 - [Blue Dyes](#)
 - [Caramel](#)
 - [General Dye Method](#)
 - [Global Method](#)
 - [Red Dyes](#)
 - [Sudan Dyes](#)
 - [Yellow Dyes](#)
- [Natural Colorants](#)
 - [Turmeric](#)
- [Flavorants](#)
 - [Amadori Compounds](#)
 - [Capsaicin](#)
 - [Quinine](#)
 - [Vanilla](#)
- [Miscellaneous](#)
 - [Caffeine](#)
 - [Choline](#)
 - [Glucosamine](#)
 - [Iodide](#)
 - [Polyphosphates](#)
- [Preservatives](#)
 - [Benzoate](#)
 - [Nitrate and Nitrite](#)
 - [Sulfites](#)
 - [Soft Drinks](#)
- [References](#)

Natural Colorants: Turmeric by UV



Chromatogram Conditions	A	B	C	D	E
Column size	4.6 × 150 mm	3.0 × 75 mm	2.1 × 50 mm	2.1 × 150 mm	2.1 × 150 mm
Particle size	4.5 μm	3.0 μm	2.2 μm	2.2 μm	2.2 μm
% Methanol	80%	80%	80%	80%	75%
Temperature	30 °C	30 °C	30 °C	30 °C	45 °C
Flow rate	1.00 mL/min	0.61 mL/min	0.41 mL/min	1.00 mL/min	1.50 mL
Data rate	2 hz	4 hz	4 hz	10 hz	20 hz
Injection volume	3.0 μL	3.0 μL	1.2 μL	1.2 μL	1.2 μL
Results					
Speed	1 time	3 times	4 times	11 times	20 times
Resolution	4.2 (R1,2)	4.2 (R1,2)	3.2 (R1,2)	2.3 (R1,2)	2.0 (R1,2)
Pressure	89 bar	140 bar	174 bar	417 bar	529 bar
Efficiency	5270 Th. Plates	4860 Th. Plates	2370 Th. Plates	1150 Th. Plates	1160 Th. Plates

Figure 4-24. Separation of pigments in turmeric. The table shows increasing degrees of acceleration by using smaller particle sizes, shorter columns, faster flow rates, and higher temperatures. With a speed-up factor as high as 20, and a solvent savings of 93%, a baseline-resolved chromatogram is still obtained.

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



Additives Including Colorants, Flavorants, and Preservatives

Natural Colorants: Turmeric by UV, FLD, and ECD

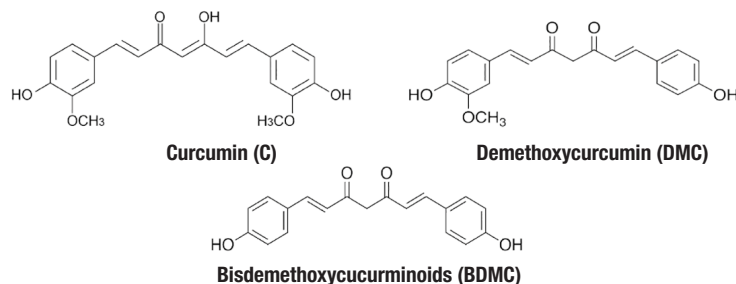


Figure 4-25. Chemical structure of curcuminoids.

Method for Figures 4-26, 4-27, and 4-28

Liquid Chromatography

An UltiMate 3000 HPLC system consisting of:

- LPG-3400BM Pump
- WPS-3000 TBSL Autosampler
- TCC-3000RS Thermostatted Column Compartment
- ECD-3000RS Electrochemical Detector with 6011RS ultra Coulometric Analytical cell, E1 = E2 = 700 mV
- FLD-3400RS Fluorescence Detector with Dual PMT, Excitation 426 nm, Emission 539 nm
- VWD-3100 Variable Wavelength Detector, 426 nm

Data Analysis and Processing

Chromatographic data were collected using Chromeleon CDS 6.8 (SR9)

Standard Preparation

Stock solution of C, DMC, and BDMC were prepared in methanol at 1 mg/mL. Working standards were prepared by dilution of stock solution in mobile phase.

Sample Preparation

Turmeric powder, yellow curry powder, and a pellet curry sauce sample were prepared at 1 mg/mL in methanol, then vortex-mixed and sonicated for 10 min. The solution was filtered through a 0.22 µm centrifuge tube filter and diluted with mobile phase prior to analysis.

Analytical Column: Acclaim RSLC PA2, 2.1 × 50 mm, 3 µm
Flow: 0.4 mL/min
Injection Volume: 2 µL
Mobile Phase: 25% 100 mM phosphate buffer, pH 3, 75% methanol

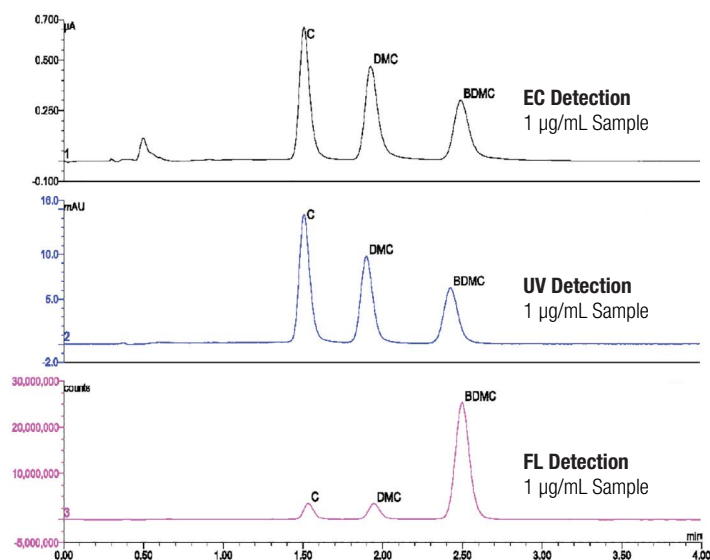


Figure 4-26. Chromatograms showing separation of curcuminoids in 1 µg/mL standard using EC, UV, or FL detection

Table 4-3. Limits of detection (LoD) of the three curcuminoids using UV, FL, and EC detection

Detector	LoD (pg on column)	
	C	DMC
UV	10	20
FL	20	20
EC	2	2

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



Additives Including Colorants, Flavorants, and Preservatives

Natural Colorants: Turmeric by UV, FLD, and ECD

Curcuminoids can be measured using HPLC with UV, FL, or EC detection. However, EC detection is the preferred approach for measurement of degradation and related products, and also has the sensitivity and selectivity for measurement of these analytes *in vivo*.

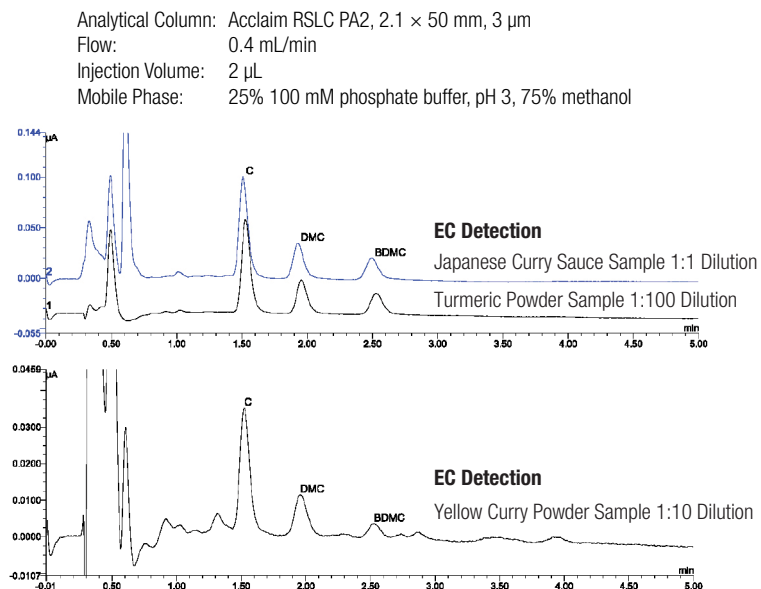


Figure 4-27. EC chromatogram of turmeric powder, yellow curry powder, and a Japanese curry sauce sample.

Table 4-4. Curcuminoid content in tested samples (ng/mg sample).

Sample Name	C	DMC
Japanese curry sauce	127	57.5
Turmeric powder	17,255	7,797
Yellow curry powder	1,804	813

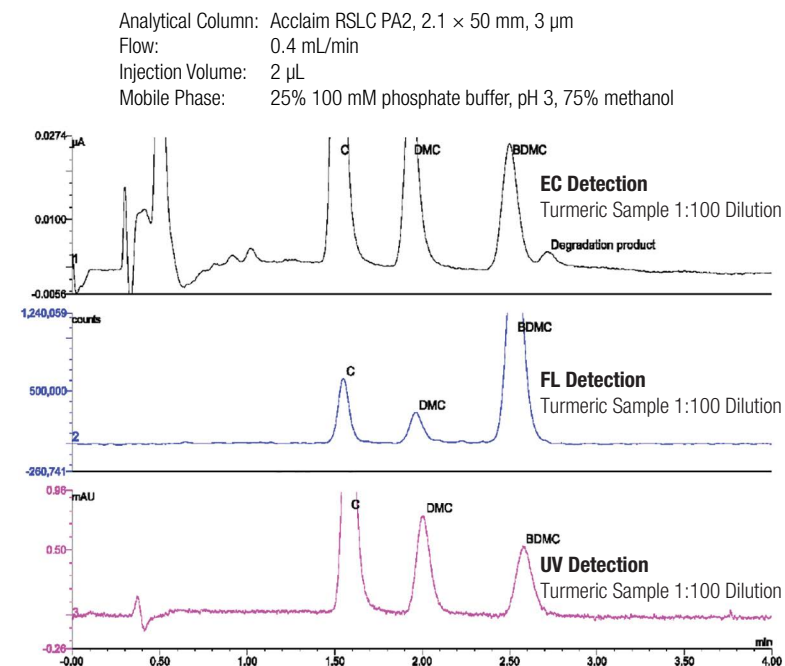


Figure 4-28. Chromatograms showing a degradation product of curcuminoids determined by EC, but not by UV or FL detection.

Table of Contents

Introduction

Analytical Technologies

Artificial Colorants

Blue Dyes

Caramel

General Dye Method

Global Method

Red Dyes

Sudan Dyes

Yellow Dyes

Natural Colorants

Turmeric

Flavorants

Amadori Compounds

Capsaicin

Quinine

Vanilla

Miscellaneous

Caffeine

Choline

Glucosamine

Iodide

Polyphosphates

Preservatives

Benzoate

Nitrate and Nitrite

Sulfites

Soft Drinks

References



Additives Including Colorants, Flavorants, and Preservatives



Flavorants

A flavorant is defined as an additive, either artificial or natural, that is used to give foods and beverages flavor. Flavorants include chemicals that can impart or enhance a particular taste e.g., sweet, sour, tangy, savory, or bitter.



Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Flavorants: Amadori Compounds

Although not an additive, *per se*, Amadori products add to flavor and appearance. Amadori products are reactive amino acid sugar conjugates formed by the Maillard reaction of an α -hydroxy carbonyl moiety of a reducing sugar and an amino group of an amino acid. This reaction can be initiated even under mild conditions; food products stored at room temperature may eventually turn brown due to the reaction of reducing sugars with amino acids. However, during high-temperature processing of food, further degradation of Amadori products, formed in the early phase of the Maillard reaction, results in the formation of complex mixtures of heterocyclic, polymeric, and other compounds. These products are directly responsible for the distinctive aroma and brown color of baked and roasted food products.



Column: Acclaim PA (3 μ m, 120 , 2.1 \times 150 mm)
Flow: 0.25 mL/min
Temperature: 30 $^{\circ}$ C
Injection Volume: 10 μ L
Eluent: 40 mmol/L formic acid in water
Detection: MSQ, ESI+, needle voltage: 3.0 kV;
probe temperature: 350 $^{\circ}$ C; cone voltage: 50 V;
dwell time: 1 μ s; SIM as indicated, time window: 0—15 min

Peaks:
1. Fru-Gly
2. Fru-Pro
3. Fru-Glu

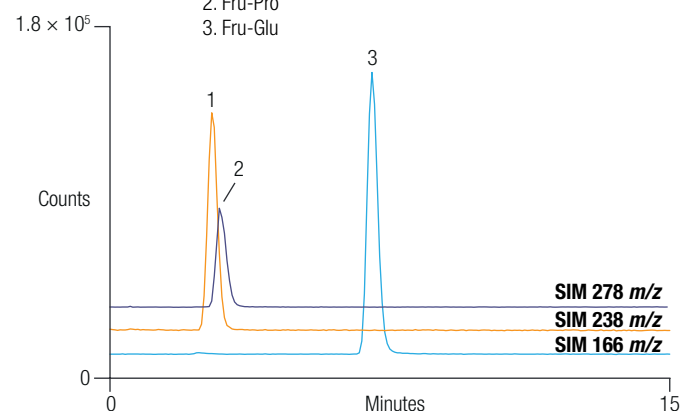


Figure 4-29. SIM chromatogram of different amadori compounds 1) fructosyl-glycine; 2) fructosyl proline and 3) fructosyl-glutamate.



Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Chili peppers (*Capsicum* sp.) are very popular ingredients in cooking worldwide. The active principles (mainly capsaicin and dihydrocapsaicin) are present in variable amounts depending on the growth conditions, so a reliable quality control method is needed. HPLC has displaced the older Scoville test based on sensory evaluation. Chilies also contain pigments, vitamins, flavinoids, and many other substances. Therefore, the ideal method should have high-resolution, high-throughput, and easy, sample cleanup. The use of the 2.2 μm Acclaim RSLC column in 2.1 mm i.d. format allows fast analysis time with reduced solvent consumptions. Shown here are two examples of separation of capsaicin and dihydrocapsaicin in dried pasilla chilies and habanero sauce. The Acclaim RSLC 120 C18 column provides excellent resolution and fast analysis at the same time, even with a very minimal sample cleanup.

Did You Know?

- The perceived heat of a chili is measured in Scoville units.
- Pure capsaicin is equivalent to 16M Scoville units.
- Resiniferatoxin, found in resin spurge is equivalent to 15B Scoville units and actually can inflict chemical burns.

Flavorants: Capsaicin by UV

LC System:	UltiMate 3000 RS
Column:	Acclaim RSLC 120 C18, 2.2 μm , 2.1 x 100 mm
Flow:	0.80 mL/min
Temperature:	60 $^{\circ}\text{C}$
Injection Volume:	2 μL ; bypass mode at 0.25 min, reset at -1.0 min.
Mobile Phases:	A: 2 mM NH_4OAc , pH 5.4 B: Acetonitrile
Time (min)	-3.0 0.0 0.25 3.0 3.01 5.0
%A	55 55 55 30 0 0
%B	45 45 45 70 100 100
Detection:	UV at 280 nm (shown); spectra 220-800 nm 10 Hz data rate, 0.3 sec time constant
Samples:	A. Natural capsaicin 50 $\mu\text{g}/\text{mL}$ in EtOH (denatured 3A) B. Dried pasilla chili - Extract 0.5 g ground chili with 20 mL EtOH and sonication. - To a 4.5 mL aliquot, add 1.0 mL heptane, 1.0 mL water, and shake. - Analyze lower (ethanol) layer C. Habanero sauce - Mix 4.5 g of sauce and 16 mL of EtOH with sonication. - To a 5 mL aliquot, add 1.0 mL heptane and shake. - Analyze lower layer.

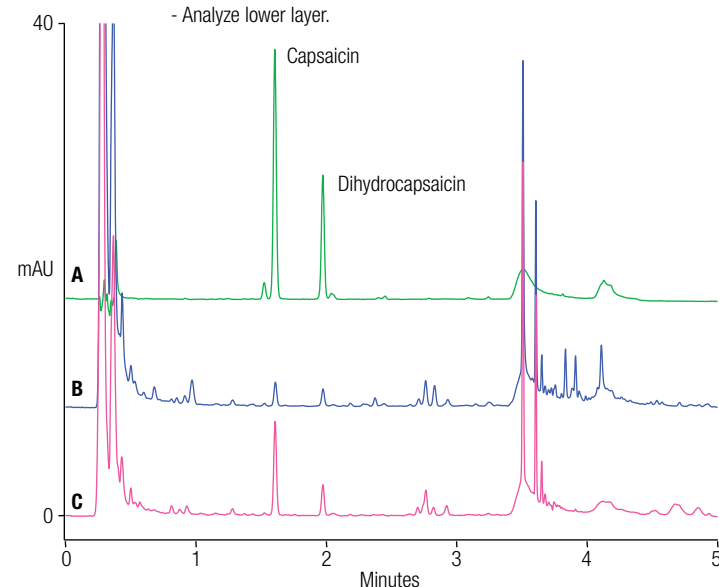


Figure 4-30. Capsaicin in chilies on Acclaim RSLC 120 C18 column.

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Additives Including Colorants, Flavorants, and Preservatives



Flavorants: Characteristics of Capsaicin

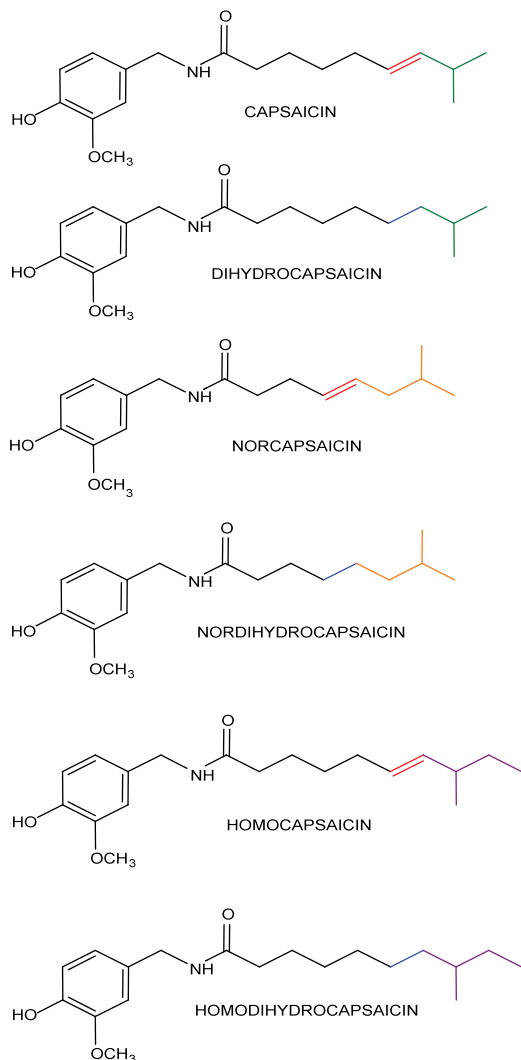


Figure 4-31. Chemical structures of different capsaicin compounds.

Table 4-5. Heat ranking of different chili peppers.

Scoville Heat Units	Examples
1,500,000–2,000,000	Trinidad Moruga Scorpion
855,000–1,463,700	Naga Viper pepper, Infinity Chilli, Bhut Jolokia chili pepper, Trinidad Scorpion Butch T pepper, Bedfordshire Super Naga, 7-Pot Chili
350,000–580,000	Red Savina habanero
100,000–350,000	Habanero chili, Scotch bonnet pepper, Datil pepper, Rocoto, Piri Piri Ndungu, Madame Jeanette, Peruvian White Habanero, Jamaican hot pepper, Guyana Wiri Wiri, Fatalii
50,000–100,000	Byadgi chilli, Bird's eye chili (aka. Thai Chili Pepper), Malagueta pepper, Chiltepin pepper, Piri piri (African bird's eye), Pequin pepper, Siling Labuyo (native chili cultivar from the Philippines)
30,000–50,000	Guntur chilli, Cayenne pepper, Aji pepper, Tabasco pepper, Cumari pepper (Capsicum Chinese)
10,000–23,000	Serrano pepper, Peter pepper, Aleppo pepper
3,500–8,000	Espelette pepper, Jalapeño pepper, Chipotle, Guajillo pepper, New Mexican peppers, Hungarian wax pepper, Tabasco sauce
1,000–2,500	Anaheim pepper (cultivar of New Mexican peppers), Poblano pepper, Rocotillo pepper, Peppadew, Sriracha sauce, Gochujang
100–900	Pimento, Peperoncini, Banana pepper, Cubanelle
No significant heat	Bell pepper, Aji dulce

Did You Know?

- Capsaicin gives chilies their burning taste
- The seeds themselves do not contain capsaicin. The highest concentration of capsaicin can be found in the white pith of the inner wall where the seeds are attached.

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Additives Including Colorants, Flavorants, and Preservatives



Flavorants: Capsaicin by ECD

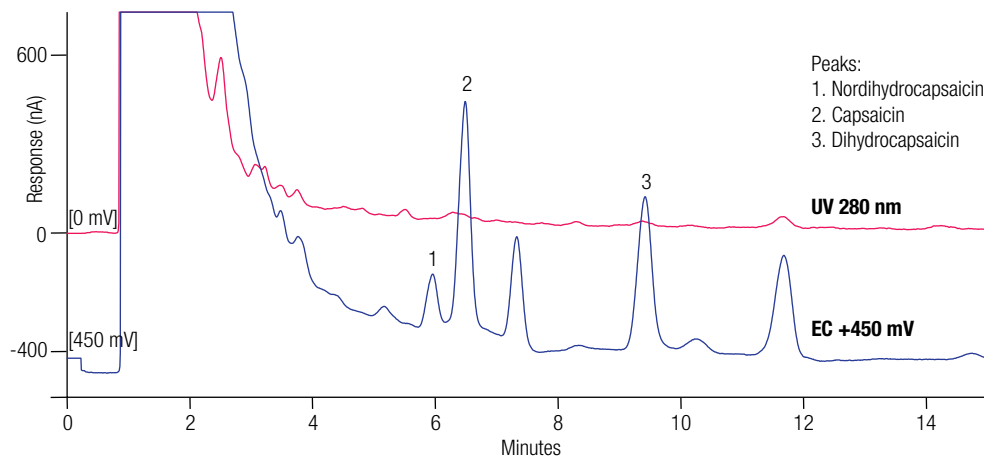


Figure 4-32. Targeted analysis of an American Spice Trade Association chili sample showing the resolution of capsaicinoids and the superior sensitivity of electrochemical detection.

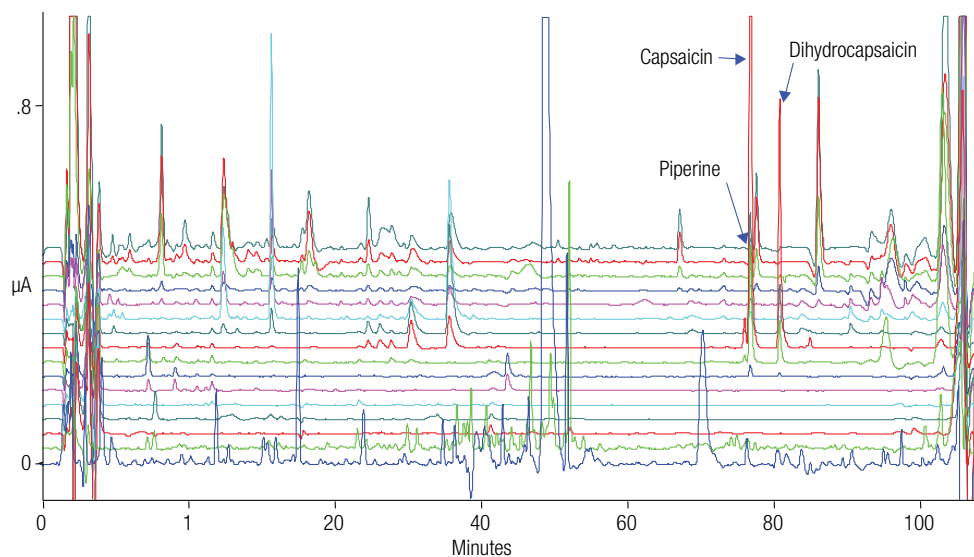


Figure 4-33. Gradient array chromatogram of chili extract.



Table of Contents

- [Introduction](#)
- [Analytical Technologies](#)
- [Artificial Colorants](#)
 - [Blue Dyes](#)
 - [Caramel](#)
 - [General Dye Method](#)
 - [Global Method](#)
 - [Red Dyes](#)
 - [Sudan Dyes](#)
 - [Yellow Dyes](#)
- [Natural Colorants](#)
 - [Turmeric](#)
- [Flavorants](#)
 - [Amadori Compounds](#)
 - [Capsaicin](#)
 - [Quinine](#)
 - [Vanilla](#)
- [Miscellaneous](#)
 - [Caffeine](#)
 - [Choline](#)
 - [Glucosamine](#)
 - [Iodide](#)
 - [Polyphosphates](#)
- [Preservatives](#)
 - [Benzoate](#)
 - [Nitrate and Nitrite](#)
 - [Sulfites](#)
 - [Soft Drinks](#)
- [References](#)

Flavorants: Quinine in Tonic Water

Quinine is a toxic bitter alkaloid extracted from the bark of the cinchona tree (*Cinchona* spp.). Quinine was the first medication to successfully treat malaria. Its addition to tonic water gave rise to the classic summer cocktail—gin and tonic—which proved an easy way for British colonists in India to take a mild dose of medicine. Today, quinine is still added to tonic water, but at a much lower level.

Did You Know?

Quinine is naturally fluorescent, causing tonic water to glow blue when exposed to UV light. Too much quinine causes quininism also known as cinchonism. Symptoms include dizziness, tinnitus and vision problems amongst others. Pilots in the armed forces are advised not to consume tonic water 3 days before flying and not to consume more than 36 oz of tonic water per day.

Column: Acclaim PA C16, 5 μ m,
4.6 \times 150 mm

Mobile Phase: A. 896 mL water, 100 mL acetonitrile,
1.40 mL methanesulfonic acid,
0.80 mL acetic acid,
2.00 mL diethylamine
B. 70:30 v:v acetonitrile:water

Gradient:

Time	0.0	10.5	10.6	15.0
%A	100	100	50	50
%B	0	0	50	50

Flow: 1.0 mL/min
Temperature: 30 $^{\circ}$ C
Inj. Volume: 25 μ L
Detection: UV at 235 nm
Sample Prep.: Filter

Peaks:

1. Cinchonidine
2. Quinine
3. Dihydroquinine
4. Benzoic acid

Reference: U.S. Pharmacopeia,
22nd Ed., 1990, 1209.

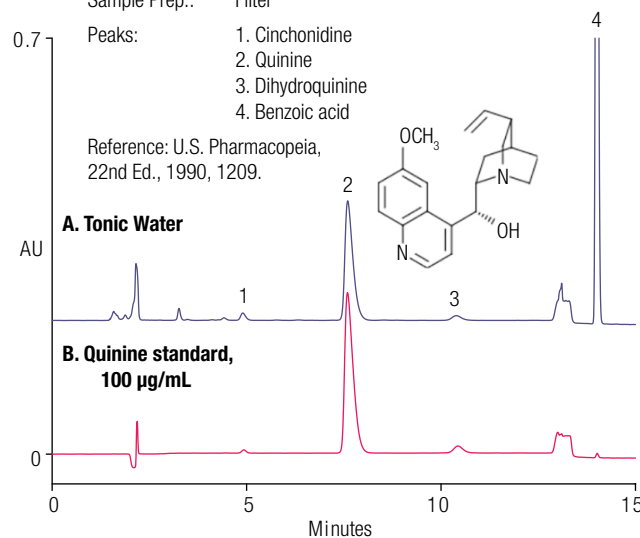


Figure 4-34. The Acclaim PolarAdvantage column can be used to rapidly determine the flavor additive quinine, as well as preservatives including benzoic acid.



Table of Contents

- [Introduction](#)
- [Analytical Technologies](#)
- [Artificial Colorants](#)
 - [Blue Dyes](#)
 - [Caramel](#)
 - [General Dye Method](#)
 - [Global Method](#)
 - [Red Dyes](#)
 - [Sudan Dyes](#)
 - [Yellow Dyes](#)
- [Natural Colorants](#)
 - [Turmeric](#)
- [Flavorants](#)
 - [Amadori Compounds](#)
 - [Capsaicin](#)
 - [Quinine](#)
 - [Vanilla](#)
- [Miscellaneous](#)
 - [Caffeine](#)
 - [Choline](#)
 - [Glucosamine](#)
 - [Iodide](#)
 - [Polyphosphates](#)
- [Preservatives](#)
 - [Benzoate](#)
 - [Nitrate and Nitrite](#)
 - [Sulfites](#)
 - [Soft Drinks](#)
- [References](#)

Flavorants: Quinine in Tonic Water

Quinine can be determined by HPLC with UV or fluorescence detection with high sensitivity. For peak confirmation, MS is used.

Column: Acclaim 120 C18, 3 μ , 120 Å, 2.1 \times 150 mm
Temperature: 35 °C
Flow: 0.25 mL/min
Injection Volume: 10 μ L
Eluent: 20% methanol, 10% acetonitrile, 70% water
Detection: 1. UV: 254 nm; 3-D: 200–800 nm
2. MSQ, ESI+, needle voltage: 3.0 kV; probe temperature: 350 °C; cone voltage: 50 V; dwell time: 0.2 s;

SIM: 324.7–325.7 m/z
Peaks: 1. Quinine 325 m/z

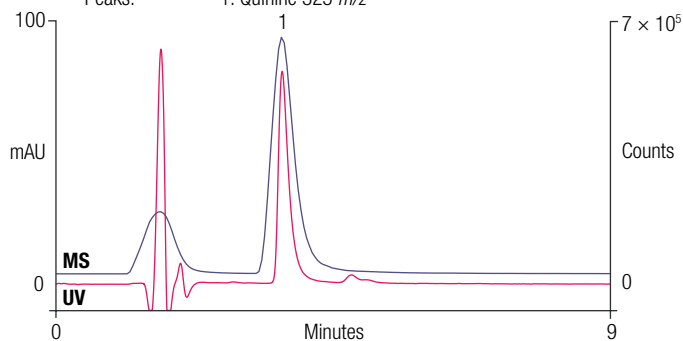


Figure 4-35. Comparison between UV and SIM chromatograms for quinine determination in tonic water.

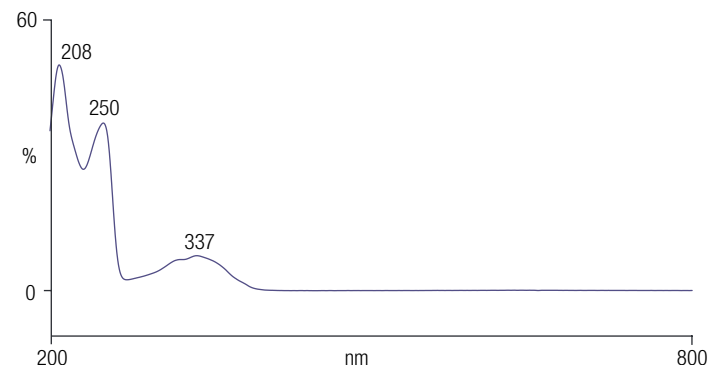


Figure 4-36. UV spectrum of quinine.

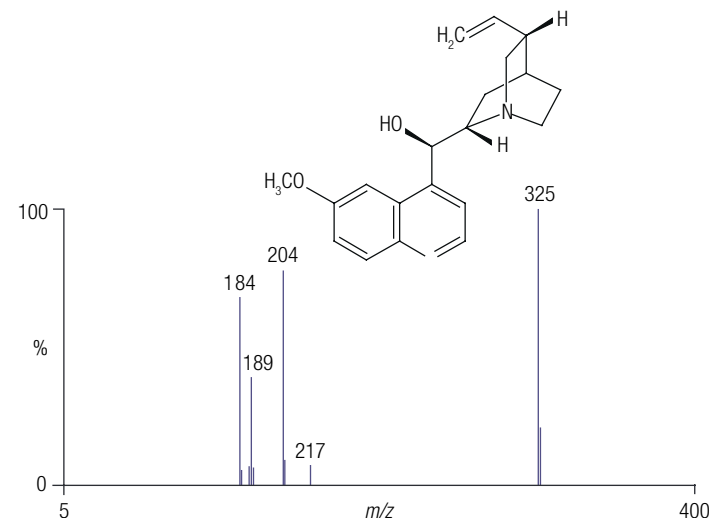


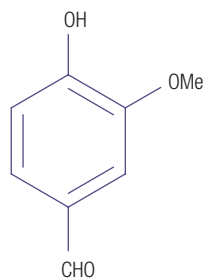
Figure 4-37. MS spectrum of quinine, ESI positive: cone voltage, 75 V; scan time, 1 s; mass range, 100–800 m/z .



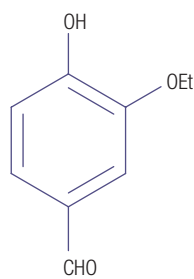
Table of Contents

- [Introduction](#)
- [Analytical Technologies](#)
- [Artificial Colorants](#)
 - [Blue Dyes](#)
 - [Caramel](#)
 - [General Dye Method](#)
 - [Global Method](#)
 - [Red Dyes](#)
 - [Sudan Dyes](#)
 - [Yellow Dyes](#)
- [Natural Colorants](#)
 - [Turmeric](#)
- [Flavorants](#)
 - [Amadori Compounds](#)
 - [Capsaicin](#)
 - [Quinine](#)
 - [Vanilla](#)
- [Miscellaneous](#)
 - [Caffeine](#)
 - [Choline](#)
 - [Glucosamine](#)
 - [Iodide](#)
 - [Polyphosphates](#)
- [Preservatives](#)
 - [Benzoate](#)
 - [Nitrate and Nitrite](#)
 - [Sulfites](#)
 - [Soft Drinks](#)
- [References](#)

Artificial vanilla flavor is significantly less expensive to manufacture than natural vanilla. Its main components are vanillin and ethyl vanillin. Coumarin is considered an adulterant in artificial vanilla, and its use is illegal in the United States. The Acclaim Polar Advantage (PA) II column demonstrates excellent selectivity for this matrix, with separation and detection of the components complete in approximately 3.5 minutes.



Vanillin
Mw = 152.15



Ethyl Vanillin
Mw = 166.17

Figure 4-38. Chemical structures for vanillin and ethyl vanillin.

Flavorants: Vanilla (Artificial)

Column: Acclaim PA2, 3 μ m 3.0 \times 33 mm
 Flow: 1.00 mL/min
 Temperature: TCC-100 at 30 $^{\circ}$ C
 Injection: ASI-100 sampler at 4 μ L
 Mobile Phase: 25:75:0.2 Methanol:water:phosphoric acid
 Detector: UVD 340U at 254 nm, spectra 220–350 nm
 Samples: A. Artificial vanilla flavor, diluted 5 \times
 B. Standards, 330 μ g/mL in 2% EtOH

Peaks:
 1. Vanillin
 2. Coumarin
 3. Ethyl vanillin
 4. Benzoic acid
 Notes:
 - Coumarin is an illegal additive
 - Benzoic acid is a preservative

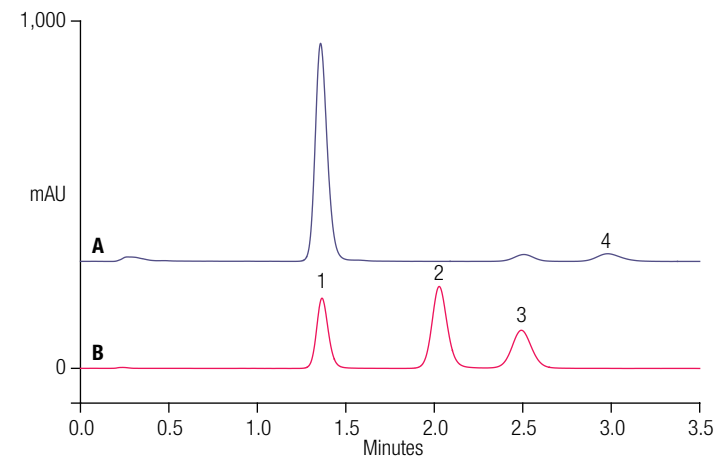


Figure 4-39. Components of artificial vanilla flavor on an Acclaim PA2 column.

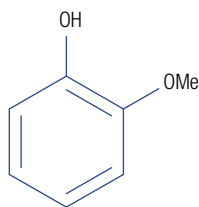


Table of Contents

- [Introduction](#)
- [Analytical Technologies](#)
- [Artificial Colorants](#)
 - [Blue Dyes](#)
 - [Caramel](#)
 - [General Dye Method](#)
 - [Global Method](#)
 - [Red Dyes](#)
 - [Sudan Dyes](#)
 - [Yellow Dyes](#)
- [Natural Colorants](#)
 - [Turmeric](#)
- [Flavorants](#)
 - [Amadori Compounds](#)
 - [Capsaicin](#)
 - [Quinine](#)
 - [Vanilla](#)
- [Miscellaneous](#)
 - [Caffeine](#)
 - [Choline](#)
 - [Glucosamine](#)
 - [Iodide](#)
 - [Polyphosphates](#)
- [Preservatives](#)
 - [Benzoate](#)
 - [Nitrate and Nitrite](#)
 - [Sulfites](#)
 - [Soft Drinks](#)
- [References](#)

Artificial Vanilla Precursors

Vanillin was first synthesized from eugenol (found in oil of clove) and later synthesized from lignin-containing sulfite liquor, a byproduct of wood pulp processing in paper manufacture. While some vanillin is still made from lignin wastes, today most synthetic vanillin is synthesized in a two-step process from the petrochemical precursors: guaiacol and glyoxylic acid.



Guaiacol
Mw = 124.14)

OCHCOOH

Glyoxylic Acid
Mw = 74.04)

Figure 4-40. Structures of guaiacol and glyoxylic acid.

Flavorants: Vanilla (Artificial)

Column:	Acclaim PA2 3 × 75 mm, 3 μm	Sample:	A. Solution of second step reaction (oxidation) B. Solution of first step reaction (condensation)
Temperature:	30 °C		
Flow:	0.5 mL/min	Peaks:	1. Impurity 2. Ethyl vanillin 3. Impurity 4. Byproduct 5. Ethyl vanillyl mandelic acid 6. Byproduct 7. Ethyl vanillin 8. Guaethol
Injection Volume:	1 μL		
Eluent:	A. Water B. Acetonitrile C. 10% Acetic acid		
Detection:	UV, at 280 nm		

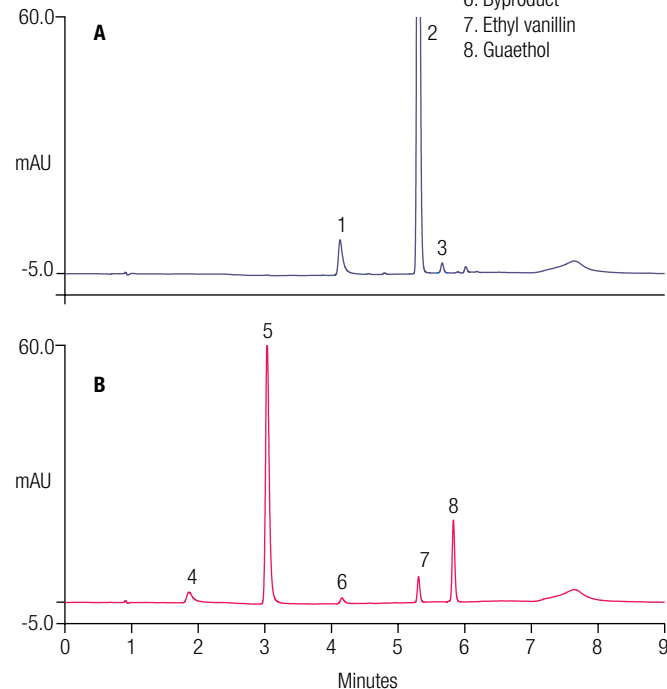


Figure 4-41. Chromatograms of solutions of ethyl vanillin synthesis process A) solution of the second step of the synthetic process. Peak 2 is ethyl vanillin, and peaks 1 and 3 are impurities, and B) solution of the first step of the synthetic process.



Table of Contents

- [Introduction](#)
- [Analytical Technologies](#)
- [Artificial Colorants](#)
 - [Blue Dyes](#)
 - [Caramel](#)
 - [General Dye Method](#)
 - [Global Method](#)
 - [Red Dyes](#)
 - [Sudan Dyes](#)
 - [Yellow Dyes](#)
- [Natural Colorants](#)
 - [Turmeric](#)
- [Flavorants](#)
 - [Amadori Compounds](#)
 - [Capsaicin](#)
 - [Quinine](#)
 - [Vanilla](#)
- [Miscellaneous](#)
 - [Caffeine](#)
 - [Choline](#)
 - [Glucosamine](#)
 - [Iodide](#)
 - [Polyphosphates](#)
- [Preservatives](#)
 - [Benzoate](#)
 - [Nitrate and Nitrite](#)
 - [Sulfites](#)
 - [Soft Drinks](#)
- [References](#)

Flavor Components of Vanilla

Natural vanilla (*Vanilla planifolia*, *V. tahitensis*, and *V. pompona*) contains many flavor components, but only four are used in official methods to assess the quality of the vanilla. The AOAC Official Method 990.25 specifies a 4.6 x 250 mm column packed with C8, 10 μ m irregular particles—a column that has only about 7000 plates, and takes over an hour to run. This figure shows the AOAC conditions implemented on Acclaim C18 columns. The RSLC column has the same number of plates and better selectivity than the originally specified column, and runs in minutes. Using well-known geometric scaling rules, the assay can be accelerated six-fold. Doubling the flow rate accelerates the analysis twelve-fold, with no sacrifice in the quality of the chromatogram.

Did You Know?

- Natural vanilla is obtained from the cured fruit of an orchid vine.
- Mexican *V. planifolia* flowers can only be pollinated by a specific *Melipone* bee found in Mexico (*abeja de monte* or mountain bee). Vines transplanted to Europe, flowered but failed to produce fruit unless hand pollinated.



Flavorants: Vanilla (Natural)

Column:	A: Acclaim 120 C18, 5 μ m, 4.6 x 150 mm B, C: Acclaim RSLC C18, 2 μ m, 2.1 x 50 mm	Detector: UV, 254 nm, A: 1 Hz data rate B: 5 Hz data rate C: 10 Hz data rate
Flow:	A: 1.00 mL/min B: 0.41 mL/min C: 0.82 mL/min	Peaks: 1. p-Hydroxybenzoic acid 2. p-Hydroxybenzaldehyde 3. Vanillic acid 4. Vanillin
Temperature:	20 °C	Sample: Commercial vanilla extract in 40% ethanol, filtered
Injection Volume:	A: 10 μ L B: 1.2 μ L C: 1.2 μ L	Reference: AOAC Official Method 990.25
Mobile Phase:	200 mM HOAc in 10% (v/v) MeOH	

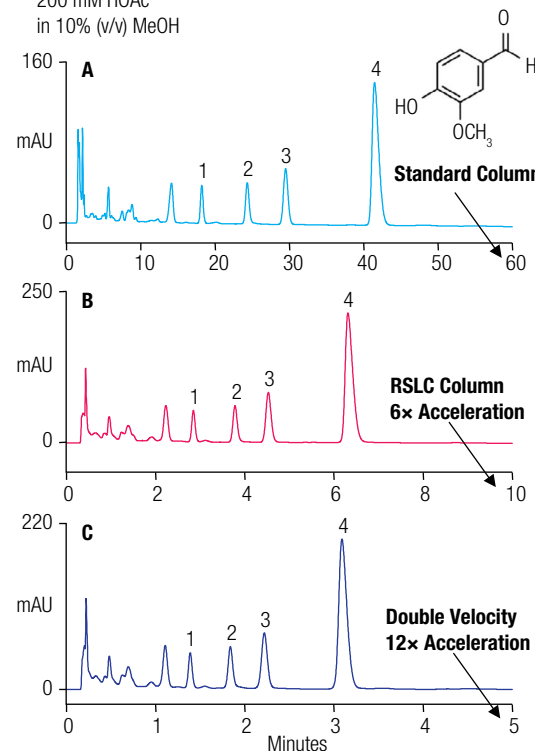


Figure 4-42. Major improvement in sample throughput when using UHPLC conditions.

Table of Contents

Introduction

Analytical Technologies

Artificial Colorants

Blue Dyes

Caramel

General Dye Method

Global Method

Red Dyes

Sudan Dyes

Yellow Dyes

Natural Colorants

Turmeric

Flavorants

Amadori Compounds

Capsaicin

Quinine

Vanilla

Miscellaneous

Caffeine

Choline

Glucosamine

Iodide

Polyphosphates

Preservatives

Benzoate

Nitrate and Nitrite

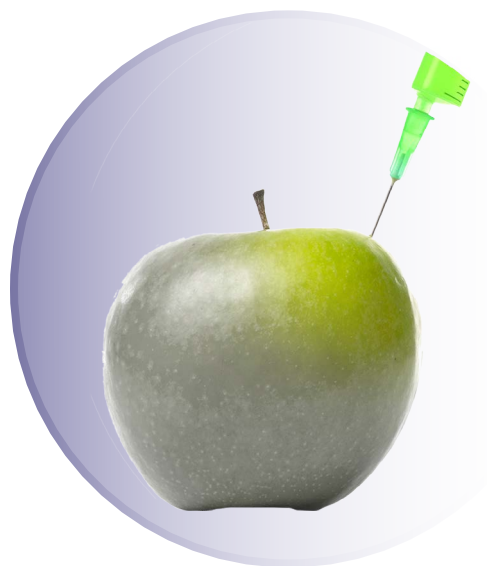
Sulfites

Soft Drinks

References



Additives Including Colorants, Flavorants, and Preservatives



Miscellaneous

There are a wide variety of chemicals that can be added to foods and beverages. These include the addition of stimulants (e.g., caffeine), fortification by inclusion of micronutrients (e.g., choline, iodide), and the use of additives to improve appearance and texture of the product (e.g., polyphosphates).



Caffeine

Table of Contents

- [Introduction](#)
- [Analytical Technologies](#)
- [Artificial Colorants](#)
 - [Blue Dyes](#)
 - [Caramel](#)
 - [General Dye Method](#)
 - [Global Method](#)
 - [Red Dyes](#)
 - [Sudan Dyes](#)
 - [Yellow Dyes](#)
- [Natural Colorants](#)
 - [Turmeric](#)
- [Flavorants](#)
 - [Amadori Compounds](#)
 - [Capsaicin](#)
 - [Quinine](#)
 - [Vanilla](#)
- [Miscellaneous](#)
 - [Caffeine](#)
 - [Choline](#)
 - [Glucosamine](#)
 - [Iodide](#)
 - [Polyphosphates](#)
- [Preservatives](#)
 - [Benzoate](#)
 - [Nitrate and Nitrite](#)
 - [Sulfites](#)
 - [Soft Drinks](#)
- [References](#)

Caffeine is a bitter xanthine alkaloid found in varying quantities in the seeds, leaves, and fruit of some plants, where it acts as a natural pesticide. It is most commonly consumed by humans in infusions extracted from the seed of the coffee plant, the leaves of the tea bush, and from products that are derived from the kola nut.

Caffeine acts as a central nervous system stimulant, temporarily warding off drowsiness and restoring alertness.

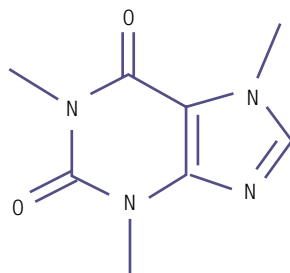


Figure 4-43. Structure of caffeine.

Did You Know?

- Unlike humans caffeine is considerably more toxic to various animals, such as dogs and birds. This is at least partly due to their poorer ability to metabolize the compound.
- Chocolate, derived from cocoa beans, typically contains only a small amount of caffeine, along with the other stimulants theobromine and theophylline. A typical 28-gram serving of a milk chocolate bar has about as much caffeine as a cup of decaffeinated coffee. Some dark chocolate contains as much as 160 mg per 100 g, about double the caffeine content by weight of the strongest caffeinated drip coffee.

Caffeine in Beverages

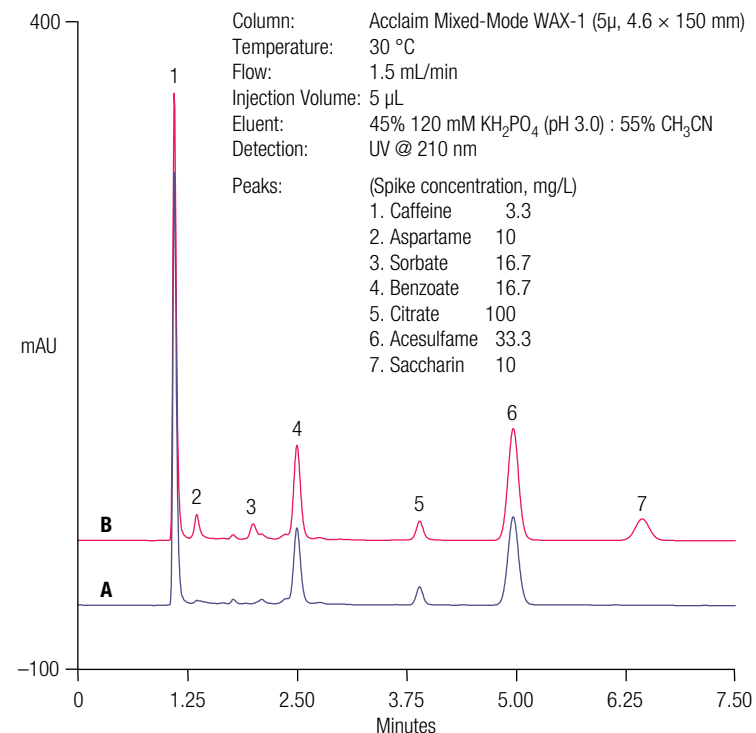


Figure 4-44. Overlay of chromatograms of a lemon-lime carbonated beverage, diluted ten-fold (A) and the same sample spiked (B).



Caffeine

Table of Contents

- [Introduction](#)
- [Analytical Technologies](#)
- [Artificial Colorants](#)
 - [Blue Dyes](#)
 - [Caramel](#)
 - [General Dye Method](#)
 - [Global Method](#)
 - [Red Dyes](#)
 - [Sudan Dyes](#)
 - [Yellow Dyes](#)
- [Natural Colorants](#)
 - [Turmeric](#)
- [Flavorants](#)
 - [Amadori Compounds](#)
 - [Capsaicin](#)
 - [Quinine](#)
 - [Vanilla](#)
- [Miscellaneous](#)
 - [Caffeine](#)
 - [Choline](#)
 - [Glucosamine](#)
 - [Iodide](#)
 - [Polyphosphates](#)
- [Preservatives](#)
 - [Benzoate](#)
 - [Nitrate and Nitrite](#)
 - [Sulfites](#)
 - [Soft Drinks](#)
- [References](#)

Caffeine in Beverages

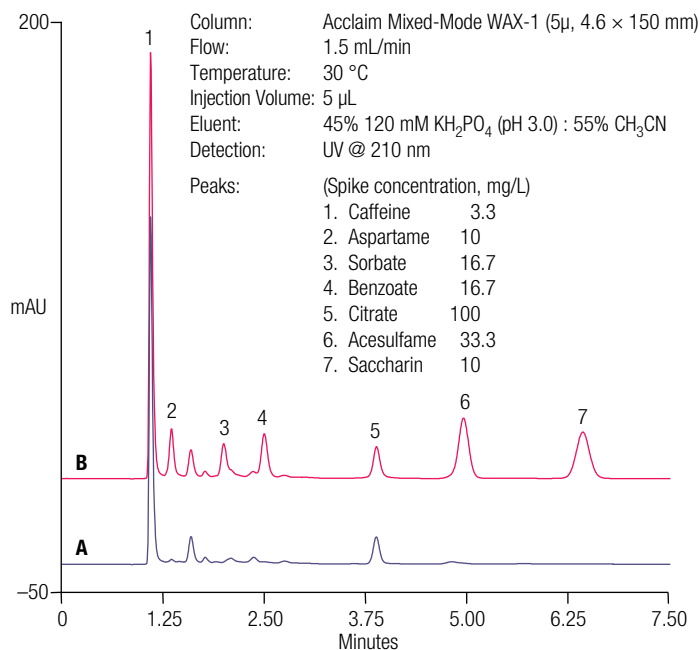


Figure 4-45. Overlay of chromatograms of cola #2, (A) diluted five-fold and (B) the same sample spiked.

Column: Acclaim RSLC 120, C18, 2.2 μ m, analytical (2.1 \times 150 mm)
 Flow: 0.45 mL/min
 Temp.: 25 $^{\circ}$ C
 Injection Volume: 1.0 μ L
 Eluent: A: 0.1% TFA, 5% acetonitrile
 B: 0.1% TFA in acetonitrile
 Gradient: 100% A, 0.0–1.2 min:
 0 to 28.5% B, 1.2–15.5 min:
 Hold at 28.5% B for 1.5 min
 Detection: Absorbance, UV 280 nm
 Sample: 1:20 Dilute white tea

Peaks:	mg/g*
1. Gallic Acid	—
2. Galocatechin	15.77
3. Epigallocatechin	16.51
4. Caffeine	—
5. Catechin	3.12
6. Epicatechin	2.73
7. Epigallocatechin Gallate	42.60
8. Galocatechin Gallate	8.83
9. Epicatechin Gallate	8.96

*Calculated concentration

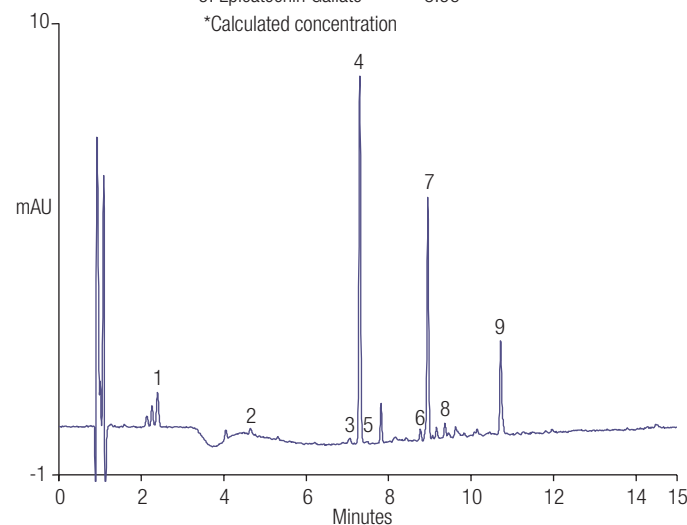


Figure 4-46. Overlay of chromatograms of cola #2, (A) diluted five-fold and (B) the same sample spiked. Separation of catechins in white tea using the Acclaim RSLC 120, C18, 2.2 μ m column.



Choline

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Choline is a water-soluble micronutrient vital to cell membrane integrity, support of methyl group metabolism, and nervous system activity. It is present as free choline in small quantities in a wide variety of foods and frequently found in its esterified forms. Choline can also be found in fortified foods and dietary supplements; e.g., choline is a required additive in many infant formulas. It is therefore important to determine the choline content of common foods.

Application Update 189 describes an improved IC method for determination of choline in infant formula and other food samples (e.g., egg powder and soy flour) using an RFIG system with suppressed conductivity detection.



Columns: Dionex IonPac CG19 and CS19, 2 mm
 Flow: 0.25 mL/min
 Injection Volume: 5 μ L
 Eluent Source: Dionex EGC III MSA cartridge with CR-CTC II trap column
 Eluent: 6.4 mM MSA
 Detection: Suppressed conductivity, Dionex CSRS 300 suppressor, 2 mm, 4 mA, recycle mode

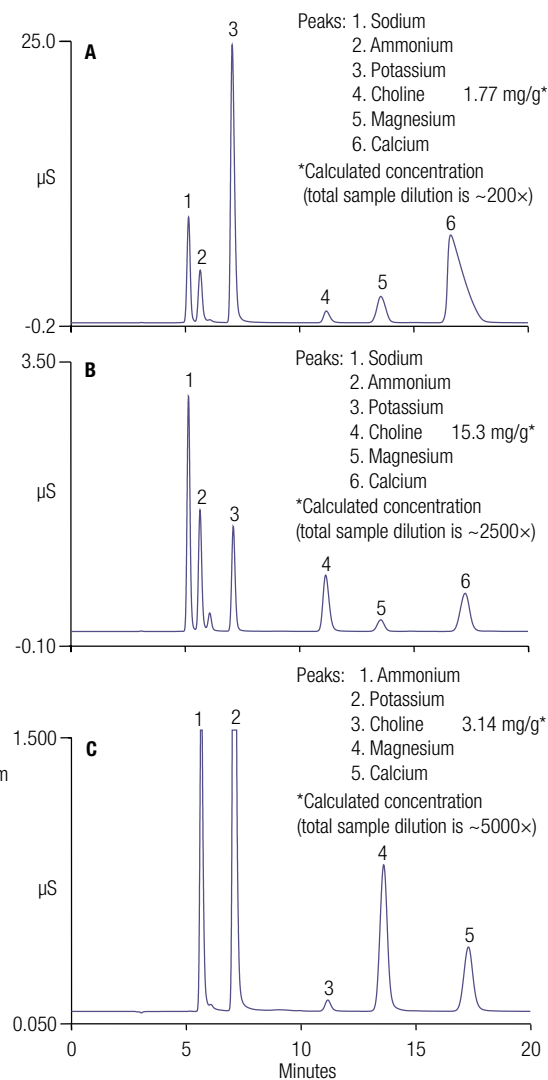


Figure 4-47. A) Determination of choline in infant formula, B) determination of choline in egg powder, and C) determination of choline in soy flour.

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



Additives Including Colorants, Flavorants, and Preservatives

Glucosamine

Glucosamine (GlcN), an amino sugar, occurs naturally in the human body. It is a major structural component in the biosynthesis of glycosaminoglycans, compounds involved in normal joint function. Use of GlcN as a dietary supplement in the management of osteoarthritis has attracted considerable attention. Results of the 2002 National Health Interview Survey showed that GlcN was one of the five nonvitamin, nonmineral herbal products/dietary supplements most frequently used by adults in the U.S.A. While the principal use for GlcN dietary supplements is for arthritis management, especially in older adults, its use as a preventive measure to maintain health and in veterinary medicine also has been reported.

The 1994 Dietary Supplement Health and Education Act granted the US FDA authority to prescribe good manufacturing practices for dietary supplements. The final rule, published in June, 2007, established regulations requiring current good manufacturing practices (cGMP) for dietary supplements. Using the cGMP regulation model for foods, the rule ensures that dietary supplements are produced in a quality manner, do not contain contaminants or impurities, and are accurately labeled.

Application Note 197 describes the use HPAE-PAD chromatography for the sensitive, selective and direct-detection of GlcN.

Column: Thermo Scientific™ Dionex™ CarboPac™ PA20, 3 mm
Flow: 0.5 mL/min
Temperature: 30 °C
Injection Volume: 10 µL
Eluent: 20 mM KOH
Detection: PAD, disposable Au electrode
Sample: Dietary supplements, about 10 µM (1.8 µg/mL)

Peaks:
1. Propylene glycol
2. Glycerol
3. myo-Inositol
4,5. Unknown
6. Sorbitol
7. Unknown
8. Mannitol
9. Glucosamine
10. Glucose
11. Fructose
12. Sucrose

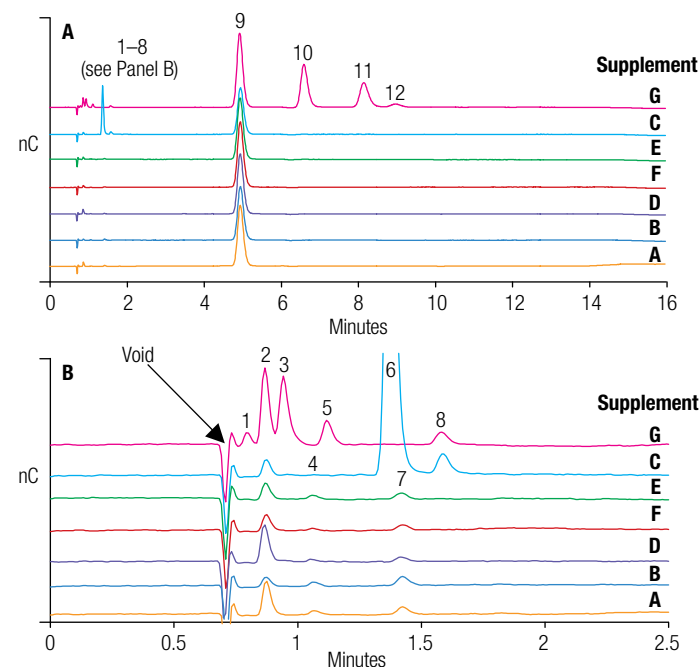


Figure 4-48. HPAE-PAD analysis of GlcN-containing dietary supplements. Seven dietary supplement samples diluted to approximately 10 µM (1.8 µg/mL) GlcN, 10-µL injection. A) Full chromatogram. B) Expanded early RT region of the chromatogram.

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



Additives Including Colorants, Flavorants, and Preservatives

Iodide

Iodide in Infant Formulas

Iodine is an important micronutrient essential for the production of thyroid hormones that are involved in the regulation of many key biochemical reactions. Iodine deficiency can lead to varying degrees of growth and developmental abnormalities in children and adults, including such illnesses as goiter and cretinism. However, an excess of iodine can also lead to thyroid disorders, especially in infants. As iodine is primarily absorbed from our diet, supplementation of iodine in food is a common practice.

The concentration of iodine in iodine-fortified foods is often regulated and monitored. Infant formula is the most highly regulated consumer food product on the market today. The Infant Formula Act of 1980 specifies minimum and maximum amounts of several nutrients, and authorizes the U.S. FDA to create and enforce standards for commercial infant formulas.

Application Note 37 includes an acetic acid digestion method for iodide extraction, coupled with an IC-PAD method for iodide detection. The IC method coupled with electrochemical detection enables selective and sensitive determination of iodide in complex matrices. An acid digestion procedure to extract iodide was optimized for milk- and soy-based infant formulas. In addition, sample preparation conditions to convert iodate to iodide for determining total iodine (i.e., iodide and iodate) are described. This method is shown to be fast, robust, and sensitive.

Column:	Dionex IonPac AG11/AS11, 4 mm	Samples:	Iodide (mg/L)
Flow:	1.5 mL/min	Standard	0.040
Temperature:	30 °C	Infant Formula Brand 1	0.027
Injectio Volume:	100 µL	Infant Formula Brand 2	0.042
Eluent:	50 mM Nitric acid	Infant Formula Brand 3	0.023
Detection:	PAD	Infant Formula Brand 4	0.035
Electrode:	Ag	Soy Infant Formula	0.046

Peaks: 1. Iodide
2. Thiocyanate

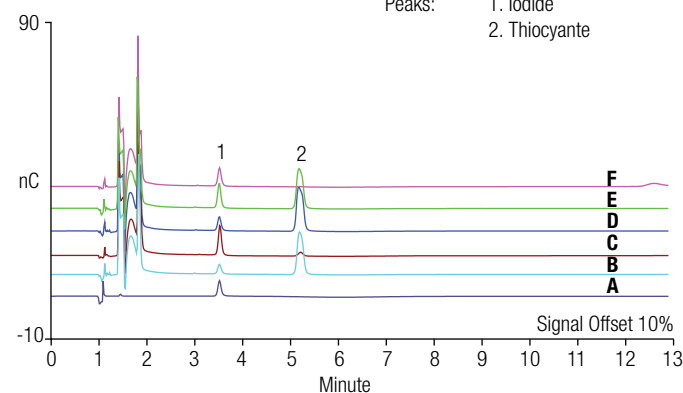


Figure 4-49. Determination of iodide in (A) DI water, (B-E) milk-based infant formulas, and (F) soy-based infant formula.

Table of Contents

- [Introduction](#)
- [Analytical Technologies](#)
- [Artificial Colorants](#)
 - [Blue Dyes](#)
 - [Caramel](#)
 - [General Dye Method](#)
 - [Global Method](#)
 - [Red Dyes](#)
 - [Sudan Dyes](#)
 - [Yellow Dyes](#)
- [Natural Colorants](#)
 - [Turmeric](#)
- [Flavorants](#)
 - [Amadori Compounds](#)
 - [Capsaicin](#)
 - [Quinine](#)
 - [Vanilla](#)
- [Miscellaneous](#)
 - [Caffeine](#)
 - [Choline](#)
 - [Glucosamine](#)
 - [Iodide](#)
 - [Polyphosphates](#)
- [Preservatives](#)
 - [Benzoate](#)
 - [Nitrate and Nitrite](#)
 - [Sulfites](#)
 - [Soft Drinks](#)
- [References](#)



Additives Including Colorants, Flavorants, and Preservatives

Iodide

Iodide in Iodized Table Salt

Iodine is an essential element naturally present in seawater and seafood. The most common forms of natural iodine in the diet are iodide and iodate, with additional iodized organic compounds providing a small fraction of bioavailable iodine. To combat disease, table salt is routinely iodized. Iodization levels vary greatly and can range from 5–100 µg/g (ppm) of iodine in salt depending on the country of manufacture and the storage conditions. Potassium iodide and potassium iodate are safe sources of dietary iodine used to iodize salt and prevent iodine deficiencies. In the United States and Canada, potassium iodide is frequently used to iodize salt whereas in many other countries, potassium iodate is preferred due to its greater stability.

Iodide can be oxidized to iodine under many conditions, including exposure to humidity, reaction with existing moisture present in the salt, exposure to sunlight, and exposure to heat. This conversion is also catalyzed by metal ions, particularly ferrous ions. Iodine readily sublimates, and is therefore easily lost from iodized salt. However, iodate is comparably more stable and is not lost by such pathways.

Application Note 236 describes a High Performance Liquid Chromatography-photodiode array UV detection method to determine iodide and iodate in, synthetic sea salt, and table salt. This method is specific, sensitive, and rapid.

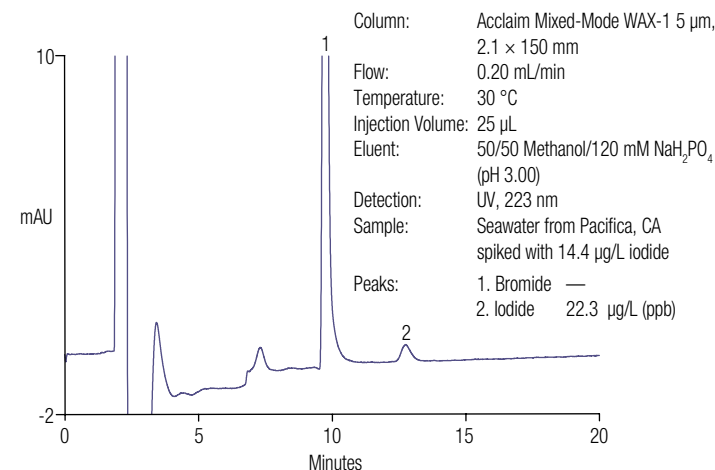


Figure 4-50. Separation of iodide and iodate in synthetic sea salt.

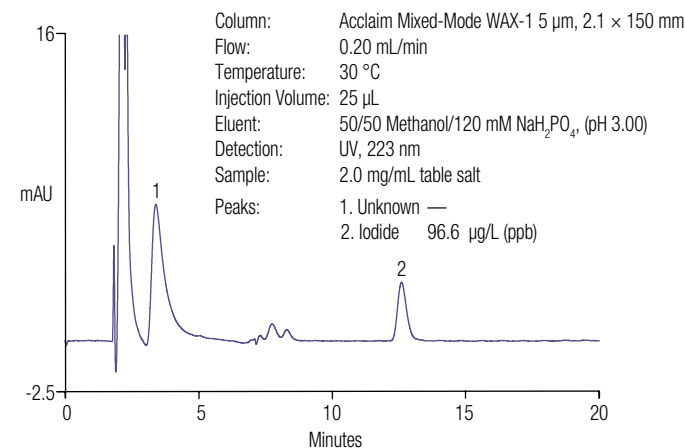


Figure 4-51. Separation of iodide in table salt on the Acclaim Mixed-Mode WAX-1 column.



Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Polyphosphates

Polyphosphates are legally added to some food products such as meat, fish, and seafood. In the seafood industry, polyphosphates are used in both fresh and frozen products to increase their water-binding capacity. This improves the appearance and texture of the product and also increases the weight of the seafood. Polyphosphates are legal, but some countries require that polyphosphate usage be declared and may also limit the amount that can be added. Citric acid is sometimes added to shrimp as a preservative to maintain an acidic environment.

Application Note 1007 describes the simultaneous determination of mono-, di-, and tri-phosphates and citrate in shrimp by ion chromatography.



Column: Dionex IonPac AS11 Analytical, 4 × 250 mm
Dionex IonPac AG11 Guard, 4 × 50 mm
Flow: 1.00 mL/min
Column Tem.: 35 °C
Tray Temp.: 10 °C
Injection Volume: 25 µL
Eluent Source: Dionex EGC III KOH, gradient mode
Gradient: 30 mM from -5 to 3 min, 45 mM from 3.1 to 7 min
50 mM from 7.5 to 13 min, (gradient curve 5)
Pressure: ~2000 psi
Detection: Suppressed conductivity, Dionex ASRS 300 suppressor,
4 mm, with external water mode, Thermo Scientific Dionex
SRS suppressor, current 130 mA
Sample: Single phosphate, polyphosphates, and citrate standard

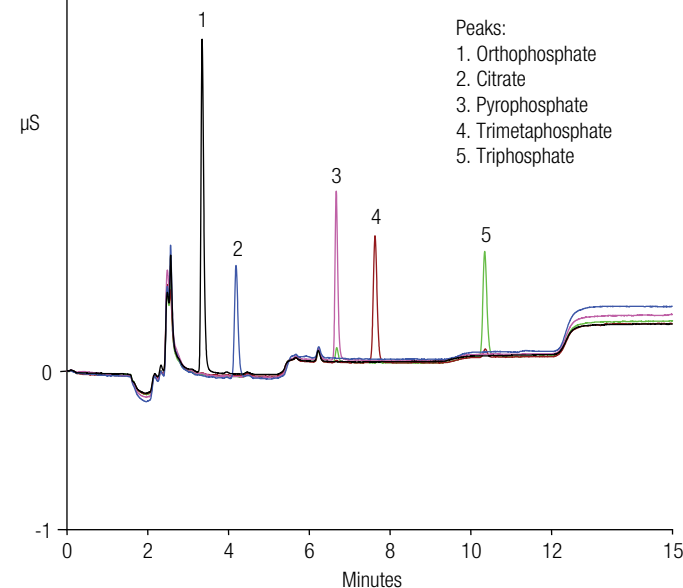


Figure 4-52. Overlay of chromatograms of single standard injections of a phosphate, three polyphosphates, and citrate.



Polyphosphates

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Column: Dionex IonPac AS11 Analytical, 4 × 250 mm
Dionex IonPac AG11 Guard, 4 × 50 mm
Flow: 1.00 mL/min
Column Temp: 35 °C
Injection Volume: 25 µL
Eluent Source: Dionex EGC III KOH, gradient mode
Gradient: 30 mM from -5 to 3 min, 45 mM from 3.1 to 7 min, 50 mM from 7.5 to 13 min, (gradient curve 5)
Pressure: ~2000 psi
Tray Temperature: 10 °C
Detection: Suppressed conductivity, Dionex ASRS 300 suppressor, 4 mm, with external water mode, Dionex SRS suppressor, current 130 mA

Sample: A. Shrimp sample at the beginning without using Dionex InGuard HRP cartridge
B. Shrimp sample at 300th injection without using Dionex InGuard HRP cartridge

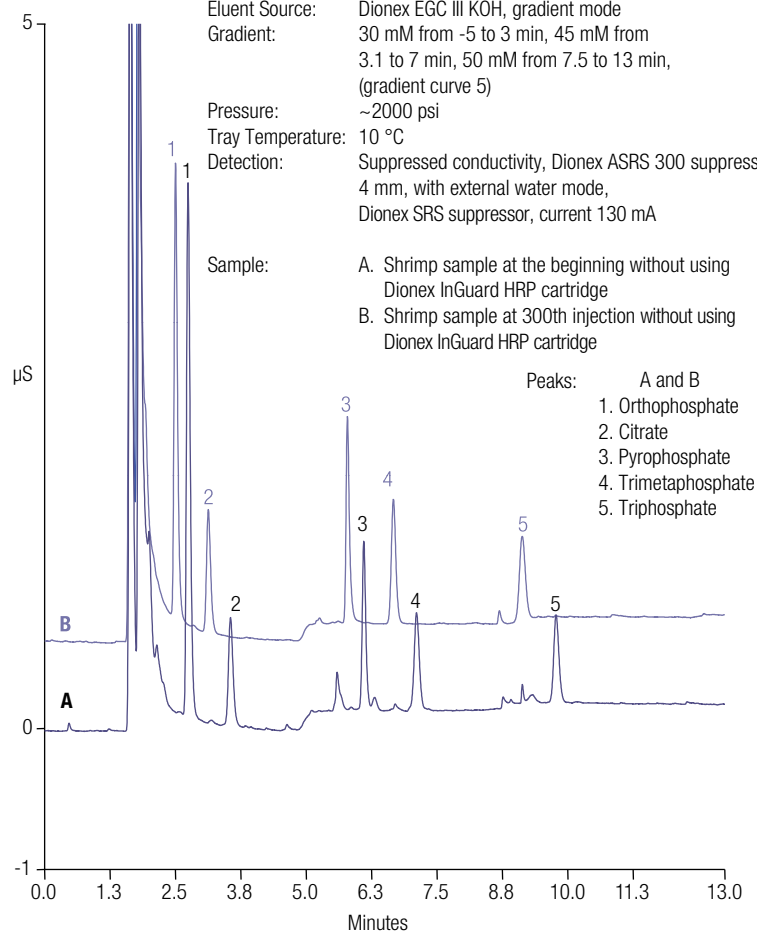


Figure 4-53. Overlay of chromatograms of a shrimp sample at the first and 300th injection using a Dionex InGuard HRP cartridge showing the value of sample preparation prior to IC analysis.

Table of Contents

Introduction

Analytical Technologies

Artificial Colorants

Blue Dyes

Caramel

General Dye Method

Global Method

Red Dyes

Sudan Dyes

Yellow Dyes

Natural Colorants

Turmeric

Flavorants

Amadori Compounds

Capsaicin

Quinine

Vanilla

Miscellaneous

Caffeine

Choline

Glucosamine

Iodide

Polyphosphates

Preservatives

Benzoate

Nitrate and Nitrite

Sulfites

Soft Drinks

References



Additives Including Colorants, Flavorants, and Preservatives



Preservatives

Artificial preservatives are man-made. They are used in foods to prevent undesirable chemical changes, reduce the risk of foodborne infections, decrease microbial spoilage, and preserve fresh attributes and nutritional quality.



Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Preservatives: Structures

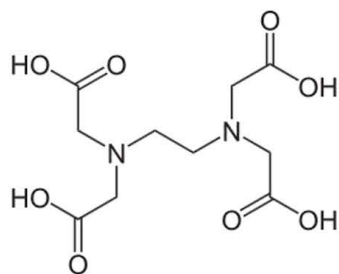


Figure 4-54A. **EDTA**. Ethylenediaminetetraacetic acid (EDTA) is an antioxidant that acts by sequestering transition metal ions (e.g., iron). EDTA is added to some food as a preservative or stabilizer to prevent catalytic oxidative decoloration. In soft drinks containing ascorbic acid and sodium benzoate, EDTA helps to prevent the formation of benzene, a carcinogen.

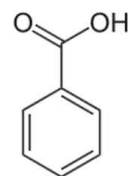


Figure 4-54B. **Benzoic acid** (E210) and its salts are preservatives. Sodium benzoate (E211) is bacteriostatic and fungistatic under acidic conditions. It is commonly added to acidic foods such as salad dressings, carbonated drinks, jams, pickles, and fruit juices.

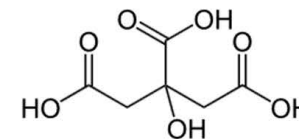


Figure 4-54C. **Citric acid**, like EDTA, acts as an antioxidant through sequestration of transition metal ions. However, due to its chemical structure, it is far less effective in this role than EDTA. Citric acid is added to caramel to prevent re-crystallization of sucrose, to ice cream as an emulsifying agent to keep fats from separating, and is generally used as a substitute for fresh lemon juice.

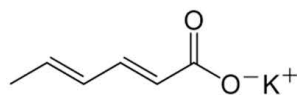


Figure 4-54D. **Potassium sorbate** (E202) is a preservative used to inhibit molds and yeasts in many foods, such as cheese, wine, yogurt, dried meats, soft drinks, fruit drinks, and baked goods. It is also added to herbal dietary supplement products to increase shelf life.

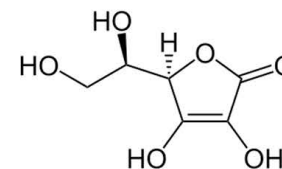


Figure 4-54E. **Erythorbic acid** (E315), previously referred to as isoascorbic acid, is a stereoisomer of ascorbic acid (vitamin C). It is a vegetable-derived food additive produced from sucrose and is widely used as an antioxidant in processed foods.



Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Since the early 1900s, benzoate has been widely used worldwide as a preservative due to its antimicrobial properties combined with its low toxicity and taste. The soft drink industry is the largest user of benzoate as a preservative due to the amount of high fructose corn syrup in many carbonated beverages.

Application Note 165 describes a simple and reliable ion chromatography method for the direct determination of benzoate in liquid food products. This method can be used to ensure that the concentration of benzoate in a sample is within product and regulatory specifications.

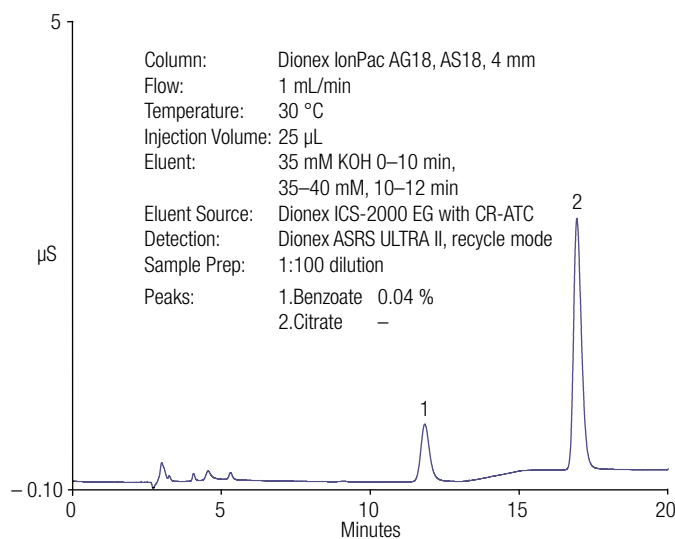


Figure 4-55. Determination of benzoate in flavored soda.

Preservatives: Benzoate

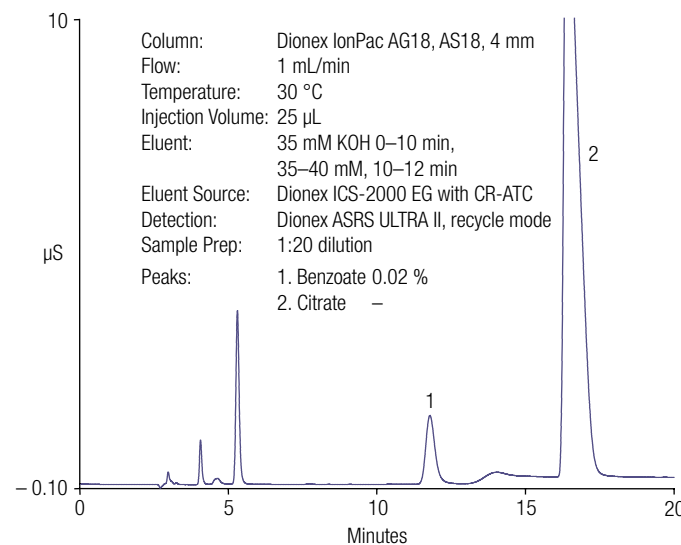


Figure 4-56. Determination of benzoate in diet soda.



Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Preservatives: Benzoate

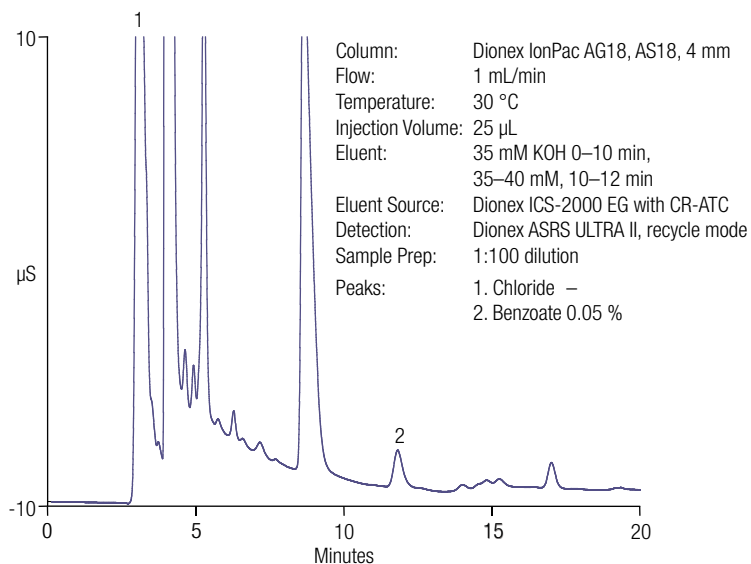


Figure 4-57. Determination of benzoate in soy sauce.

Column: Dionex IonPac AG18, AS18, 4 mm
Flow: 1 mL/min
Temperature: 30 °C
Injection Volume: 25 µL
Eluent: 35 mM KOH 0–10 min,
35–40 mM, 10–12 min
Eluent Source: Dionex ICS-2000 EG with CR-ATC
Detection: Dionex ASRS ULTRA II, recycle mode
Sample Prep: 1:100 dilution
Peaks: 1. Benzoate 0.05 %
2. Citrate –

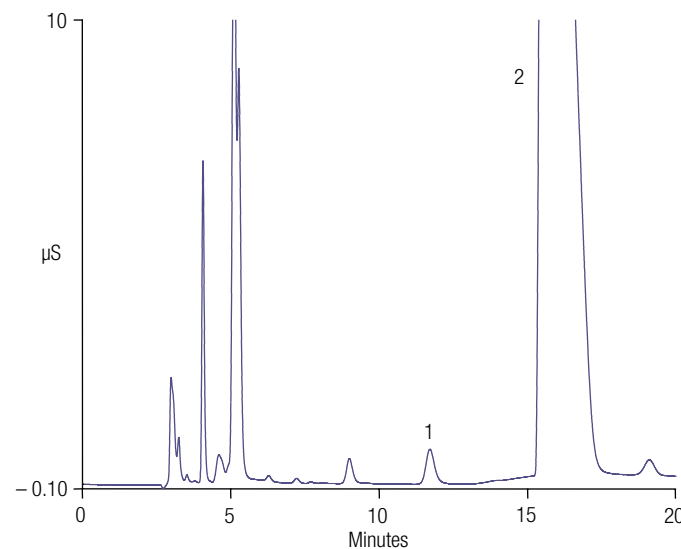


Figure 4-58. Determination of benzoate in lemon juice.



Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



Additives Including Colorants, Flavorants, and Preservatives

Preservatives: Nitrate and Nitrite

Nitrate and nitrite are usually added to processed meat products to protect against microorganisms that can cause food poisoning, such as *Clostridium botulinum*. However, nitrite can react with secondary amines to form nitrosoamines, a class of carcinogenic compounds, in food products or in the digestive system. Nitrate, although more stable than nitrite, can act as a reservoir for nitrite. Also, nitrate can readily be converted into nitrite by microbial reduction. Thus, both nitrate and nitrite must be monitored to ensure the quality and safety of meat products.

Application Note 112 describes an accurate and sensitive method in which nitrate and nitrite are extracted from meat products and then determined directly using anion exchange chromatography with UV detection. Commercially available ham and salami were used as model samples.

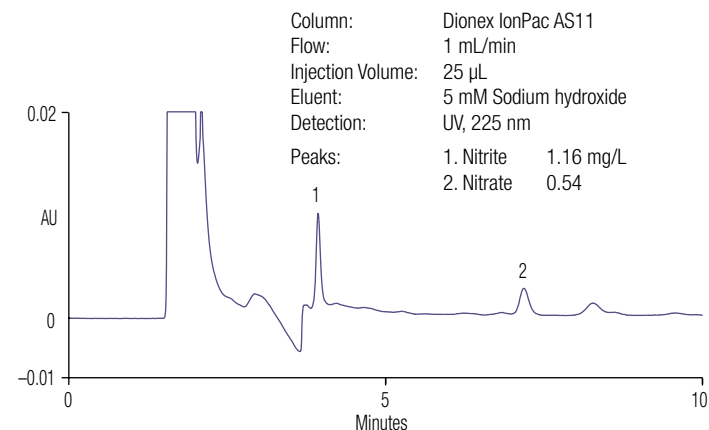


Figure 4-59. Separation of nitrate and nitrite from ham.

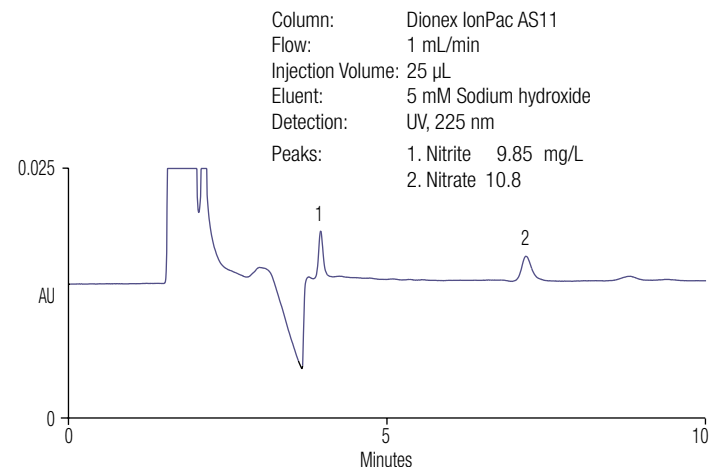


Figure 4-60. Separation of nitrate and nitrite from salami.



Table of Contents

- [Introduction](#)
- [Analytical Technologies](#)
- [Artificial Colorants](#)
 - [Blue Dyes](#)
 - [Caramel](#)
 - [General Dye Method](#)
 - [Global Method](#)
 - [Red Dyes](#)
 - [Sudan Dyes](#)
 - [Yellow Dyes](#)
- [Natural Colorants](#)
 - [Turmeric](#)
- [Flavorants](#)
 - [Amadori Compounds](#)
 - [Capsaicin](#)
 - [Quinine](#)
 - [Vanilla](#)
- [Miscellaneous](#)
 - [Caffeine](#)
 - [Choline](#)
 - [Glucosamine](#)
 - [Iodide](#)
 - [Polyphosphates](#)
- [Preservatives](#)
 - [Benzoate](#)
 - [Nitrate and Nitrite](#)
 - [Sulfites](#)
 - [Soft Drinks](#)
- [References](#)

In the food and beverage industries, sulfites are a group of compounds that includes sulfur dioxide and sulfite salts. Sulfites can occur naturally in some foods and beverages due to fermentation. For centuries, sulfiting agents—such as sodium sulfite, sodium bisulfite, and sodium metabisulfite—have been used as preservatives to prevent microbial spoiling and browning reactions in a wide variety of food and beverage products.

Conditions for Figures 4-61–4-64

Columns: Dionex IonPac ICE-AS1 Guard, 4 × 50 mm (P/N 067842)
 Dionex IonPac ICE-AS1 Analytical, 4 × 250 mm (P/N 064198)

Temperature: 25 °C (upper compartment, detector)
 30 °C (lower compartment, column)

Eluent: 20 mM MSA

Tray Temp: 4 °C

Detection: PAD, disposable Pt working electrode

System Backpressure: ~1760 psi

Background: 50–60 nC

Noise: 0.04 nC peak-to-peak

Run Time: 25 min

Waveform for the ED:

Time (s)	Potential (V)	Last Step*	Ramp*	Gain Region*	Integration
0.00	0.8	Off	On	Off	Off
0.40	0.8	Off	On	On	On
0.60	0.8	Off	On	On	Off
0.61	1.2	Off	On	Off	Off
0.70	1.2	Off	On	Off	Off
0.71	0.1	Off	On	Off	Off
1.00	0.1	On	On	Off	Off

*These settings are required in the Dionex ICS-3000/5000 system but not used in older Dionex systems; the reference electrode is in AgCl mode (Ag/AgCl reference electrode).

Preservatives: Sulfites

Sulfites have been implicated as the cause of allergic reactions that range in severity from minor to life threatening. Sulfites also have been reported as the cause of some asthmatic responses in certain people. Therefore, since 1986, the U.S. Food and Drug Administration (FDA) has required labeling of any food or beverage containing a sulfite concentration >10 ppm.

Application Note 54 describes an improved ion-exclusion chromatography method for the determination of total and free sulfite in foods and beverages, that does not have the analytical problems associated with AOAC Methods 990.28 and 990.31.

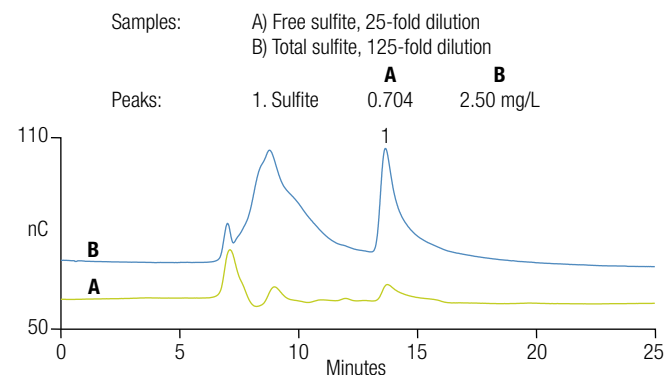


Figure 4-61. Chromatograms of A) total and B) free sulfite in red wine. A 15% signal offset has been applied.



Table of Contents

- [Introduction](#)
- [Analytical Technologies](#)
- [Artificial Colorants](#)
 - [Blue Dyes](#)
 - [Caramel](#)
 - [General Dye Method](#)
 - [Global Method](#)
 - [Red Dyes](#)
 - [Sudan Dyes](#)
 - [Yellow Dyes](#)
- [Natural Colorants](#)
 - [Turmeric](#)
- [Flavorants](#)
 - [Amadori Compounds](#)
 - [Capsaicin](#)
 - [Quinine](#)
 - [Vanilla](#)
- [Miscellaneous](#)
 - [Caffeine](#)
 - [Choline](#)
 - [Glucosamine](#)
 - [Iodide](#)
 - [Polyphosphates](#)
- [Preservatives](#)
 - [Benzoate](#)
 - [Nitrate and Nitrite](#)
 - [Sulfites](#)
 - [Soft Drinks](#)
- [References](#)

Preservatives: Sulfites

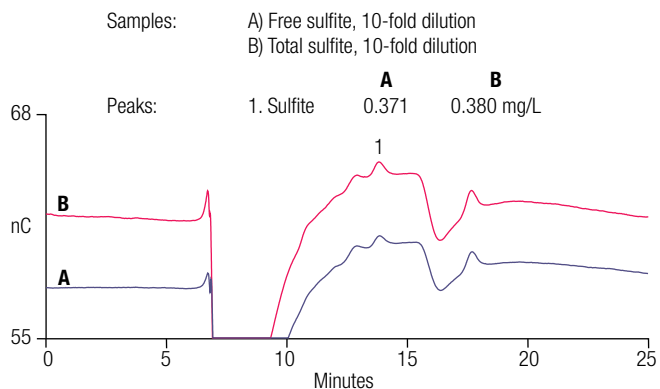


Figure 4-62. Chromatograms of A) free and B) total sulfite in coconut water.

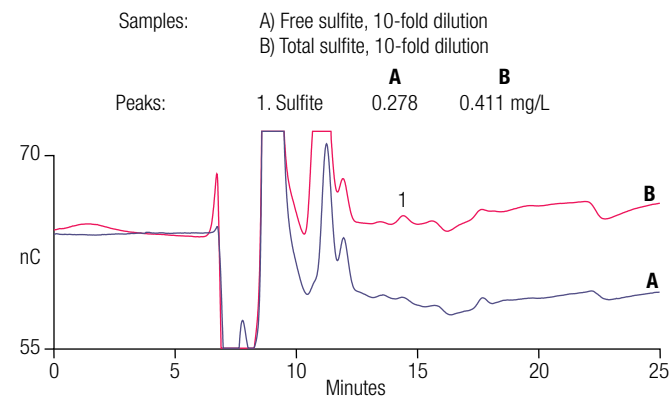


Figure 4-64. Chromatograms of A) free and B) total sulfite in coconut water.

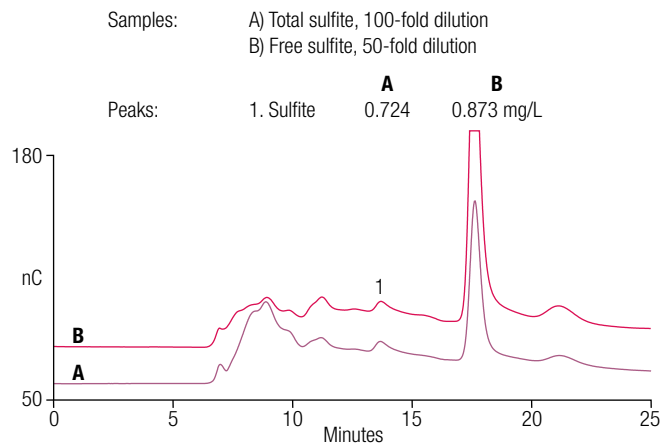


Figure 4-63. Chromatograms of A) total and B) free sulfite in red wine. A 15% signal offset has been applied.





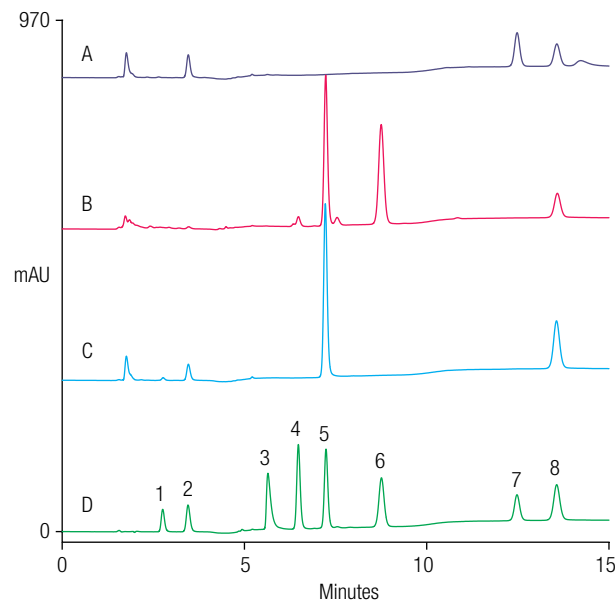
Preservatives: Soft Drinks

Soft drink formulations often contain additives such as artificial sweeteners, preservatives, citric acid and caffeine. This analysis determines eight major ingredients in a single run. The Acclaim OA column is uniquely suited because of its aqueous compatibility for the separation of organic acids, and its hydrophobic selectivity for the separation of preservatives.

Table of Contents

- [Introduction](#)
- [Analytical Technologies](#)
- [Artificial Colorants](#)
 - [Blue Dyes](#)
 - [Caramel](#)
 - [General Dye Method](#)
 - [Global Method](#)
 - [Red Dyes](#)
 - [Sudan Dyes](#)
 - [Yellow Dyes](#)
- [Natural Colorants](#)
 - [Turmeric](#)
- [Flavorants](#)
 - [Amadori Compounds](#)
 - [Capsaicin](#)
 - [Quinine](#)
 - [Vanilla](#)
- [Miscellaneous](#)
 - [Caffeine](#)
 - [Choline](#)
 - [Glucosamine](#)
 - [Iodide](#)
 - [Polyphosphates](#)
- [Preservatives](#)
 - [Benzoate](#)
 - [Nitrate and Nitrite](#)
 - [Sulfites](#)
 - [Soft Drinks](#)
- [References](#)

Column:	Acclaim OA 5 µm, 4.0 × 150 mm																					
Flow:	0.8 mL/min																					
Temperature:	30 °C																					
Injection Volume:	10 µL																					
Mobile Phase:	(A) 14.2 g/L Na ₂ SO ₄ + 0.550 mL/L CH ₃ SO ₃ H in water (B) Methanol																					
Gradient:	<table border="0" style="font-size: small;"> <tr> <td>Time</td> <td>0.0</td> <td>1.0</td> <td>2.0</td> <td>6.5</td> <td>7.5</td> <td>15.0</td> </tr> <tr> <td>%A</td> <td>100</td> <td>100</td> <td>69</td> <td>69</td> <td>60</td> <td>60</td> </tr> <tr> <td>%B</td> <td>0</td> <td>0</td> <td>31</td> <td>31</td> <td>40</td> <td>40</td> </tr> </table>	Time	0.0	1.0	2.0	6.5	7.5	15.0	%A	100	100	69	69	60	60	%B	0	0	31	31	40	40
Time	0.0	1.0	2.0	6.5	7.5	15.0																
%A	100	100	69	69	60	60																
%B	0	0	31	31	40	40																
Detection:	UV at 210 nm; optional 230, 262 nm																					
Samples:	(A) Minute-Maid® Orange Soda (B) Diet Coke® Soda (C) Mountain Dew® Soda (D) Standards																					
Sample Prep.:	Minute-Maid and Diet Coke are registered trademarks of Coca-Cola. Mountain Dew is a registered trademark of PepsiCo. Filter and dilute 5× with mobile phase A																					
Peaks:	<table border="0" style="font-size: small;"> <tr> <td>1. Erythorbic acid</td> <td>50* µg/mL</td> </tr> <tr> <td>2. Citric acid</td> <td>500</td> </tr> <tr> <td>3. Acesulfame-K</td> <td>50</td> </tr> <tr> <td>4. Aspartylphenylalanine</td> <td>50</td> </tr> <tr> <td>5. Caffeine</td> <td>12</td> </tr> <tr> <td>6. Aspartame</td> <td>50</td> </tr> <tr> <td>7. Sorbate potassium</td> <td>50</td> </tr> <tr> <td>8. Benzoic acid</td> <td>50</td> </tr> </table>	1. Erythorbic acid	50* µg/mL	2. Citric acid	500	3. Acesulfame-K	50	4. Aspartylphenylalanine	50	5. Caffeine	12	6. Aspartame	50	7. Sorbate potassium	50	8. Benzoic acid	50					
1. Erythorbic acid	50* µg/mL																					
2. Citric acid	500																					
3. Acesulfame-K	50																					
4. Aspartylphenylalanine	50																					
5. Caffeine	12																					
6. Aspartame	50																					
7. Sorbate potassium	50																					
8. Benzoic acid	50																					



Note: For alternate calibration mixes, use ascorbic acid instead of erythorbic acid, or use saccharin instead of aspartame.

*Listed concentrations are for the standards (chromatogram D).

Figure 4-65. Soft drink ingredients using the Acclaim OA column.

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Additives Including Colorants, Flavorants, and Preservatives



Preservatives: Soft Drinks

Soft drinks can contain many additives including caffeine, preservatives, artificial sweeteners, and other flavor components. Since many of these compounds are weak acids, they are retained on the Acclaim Mixed-Mode WAX-1 column. The pH may be lowered to favor reversed-phase

retention, or raised to promote anion-exchange retention. In this way, the separation can be adjusted to the particular requirements of a specific beverage formulation.

Column: Acclaim Mixed-Mode WAX-1, 5 μ m
Flow: 1 mL/min
Temperature: 30 $^{\circ}$ C
Injection Volume: 2.5 μ L
Dimension: 4.6 \times 150 mm
Mobile Phase: (A) 55/45 v/v Acetonitrile/
0.2 M phosphate buffer, pH6.0
(B) 57/43 v/v Acetonitrile/
0.12 M phosphate buffer, pH2.9
Detection: UV @ 210 nm
Peaks: 1. Caffeine 100 μ g/mL
2. Aspartame 100
3. Acesulfame, potassium 100
4. Saccharin 100
5. Sorbate, potassium 100
6. Benzoic acid 100
7. Citric acid 300

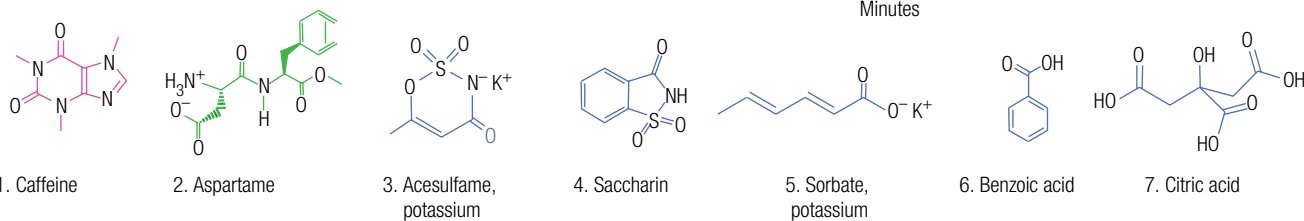
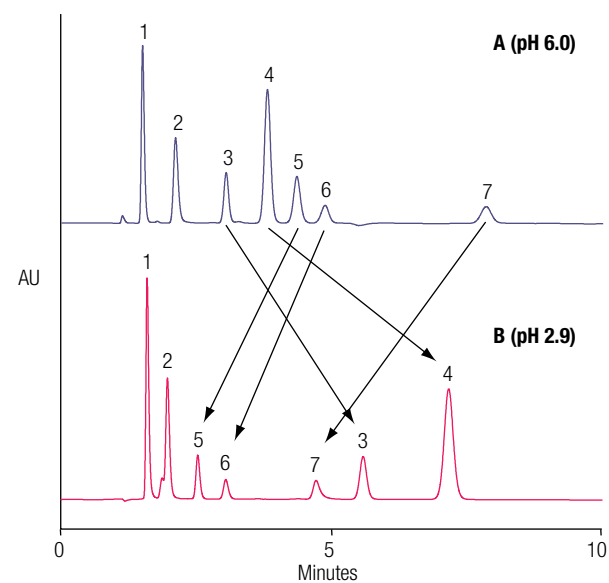


Figure 4-66. Analysis of additives used in soft drinks on Acclaim Mixed-Mode WAX-1 column.



Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Preservatives: Soft Drinks

Rapid analysis is required for quality control testing in highly productive, modern factories. Using the additional leverage provided by its multifunctional chemistry, the Acclaim Mixed-Mode WAX-1 column can be optimized for a 3-minute assay of cola drinks.

Column:	Acclaim Mixed-Mode WAX-1, 5 μ m, 4.6 \times 150 mm	Peaks: 1. Caffeine 2. Aspartame 3. Sorbate 4. Benzoate 5. Citrate 6. Acesulfame
Flow:	2 mL/min	
Temperature:	30 $^{\circ}$ C	
Injection Volume:	2.5 μ L	
Mobile Phase:	57/43 v/v Acetonitrile/ 120 mM phosphate buffer, pH 2.9	
Detection:	UV @ 210 nm	
Sample:	Direct injection of degassed sample	

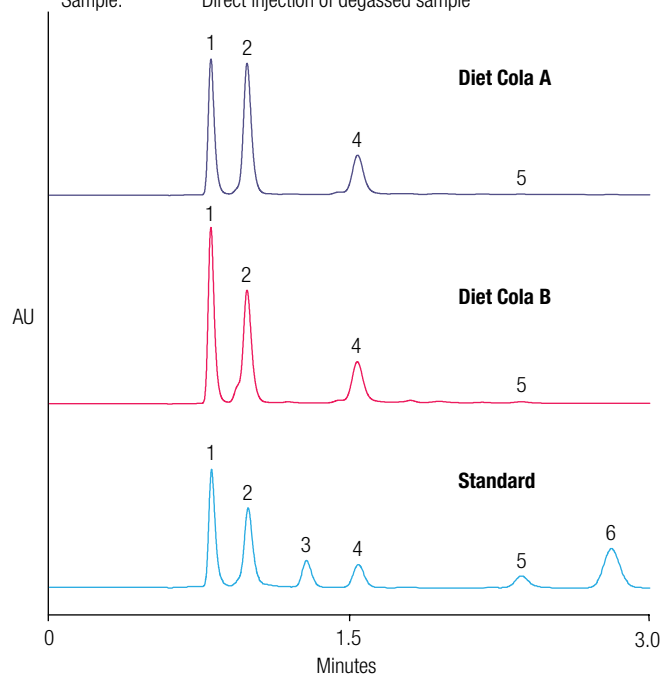


Figure 4-67. Rapid analysis of soft drinks with the Acclaim Mixed-Mode WAX-1 column.



Table of Contents

Introduction

Analytical Technologies

Artificial Colorants

Blue Dyes

Caramel

General Dye Method

Global Method

Red Dyes

Sudan Dyes

Yellow Dyes

Natural Colorants

Turmeric

Flavorants

Amadori Compounds

Capsaicin

Quinine

Vanilla

Miscellaneous

Caffeine

Choline

Glucosamine

Iodide

Polyphosphates

Preservatives

Benzoate

Nitrate and Nitrite

Sulfites

Soft Drinks

References



References



Technical Collateral and Peer Reviewed Journals

Here you'll find a multitude of references using our HPLC, ion chromatography and sample preparation solutions.

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



References



HPLC and UHPLC References

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



References

Peer Reviewed Journals: HPLC and UHPLC Methods

Carbohydrates

Title	Authors	Publication	Publication Date
Carbohydrate and oligosaccharide analysis with a universal HPLC detector.	Asa, D.	<i>American Laboratory</i> 38, 16.	2006
Determination of levoglucosan in atmospheric aerosols using high performance liquid chromatography with aerosol charge detection.	Dixon, R. W.; Baltzell, G.	<i>J. Chromatogr., A.</i> 1109 (2), 214–221	2006 Mar 24
Composition of structural carbohydrates in biomass: Precision of a liquid chromatography method using a neutral detergent extraction and a charged aerosol detector.	Godin, B.; Agneessens, R.; Gerin, P. A.; Delcarte, J.	<i>Talanta</i> 85 (4), 2014–2026	2011 Sep 30
Selectivity issues in targeted metabolomics: Separation of phosphorylated carbohydrate isomers by mixed-mode hydrophilic interaction/weak anion exchange chromatography.	Hinterwirth, H.; Lämmerhofer, M.; Preinerstorfer, B.; Gargano, A.; Reischl, R.; Bicker, W.; Trapp, O.; Brecker, L.; Lindner, W.	<i>J. Sep. Sci.</i> 33 (21), 3273–3282	2010 Nov
Investigation of polar organic solvents compatible with Corona charged aerosol detection and their use for the determination of sugars by hydrophilic interaction liquid chromatography.	Hutchinson, J. P.; Remenyi, T.; Nesterenko, P.; Farrell, W.; Groeber, E.; Szucs, R.; Dicinowski, G.; Haddad, P. R.	<i>Anal. Chim. Acta.</i> 750, 199–206	2012 Oct 31
Characterization of an endoglucanase belonging to a new subfamily of glycoside hydrolase family 45 of the basidiomycete <i>Phanerochaete chrysosporium</i>.	Igarashi, K.; Ishida, T.; Hori, C.; Samejima, M.	<i>Appl. Environ. Microbiol.</i> 74 (18), 5628–5634	2008 Sep
Direct detection method of oligosaccharides by high-performance liquid chromatography with charged aerosol detection.	Inagaki, S.; Min, J. Z.; Toyo'oka, T.	<i>Biomed. Chromatogr.</i> 21 (4), 338–342	2007 Apr
Differential selectivity of the <i>Escherichia coli</i> cell membrane shifts the equilibrium for the enzyme-catalyzed isomerization of galactose to tagatose.	Kim, J. H.; Lim, B. C.; Yeom, S. J.; Kim, Y. S.; Kim, H. J.; Lee, J. K.; Lee, S. H.; Kim, S. W.; Oh, D. K.	<i>Appl. Environ. Microbiol.</i> 74 (8), 2307–2313	2008 Apr
Elution strategies for reversed-phase high-performance liquid chromatography analysis of sucrose alkanolate regioisomers with charged aerosol detection.	Lie, A.; Pedersen, L. H.	<i>J. Chromatogr., A.</i> 1311, 127–133	2013 Oct 11
Design of experiments and multivariate analysis for evaluation of reversed-phase high-performance liquid chromatography with charged aerosol detection of sucrose caprate regioisomers	Lie, A.; Wimmer, R.; Pedersen, L. H.	<i>J. Chromatogr., A.</i> 1281, 67–72	2013 Mar 15
Solvent effects on the retention of oligosaccharides in porous graphitic carbon liquid chromatography	Melmer, M.; Stangler, T.; Premstaller, A.; Lindner, W.	<i>J. Chromatogr., A</i> 1217 (39) 6092–6096	2010 Sep 24
Practical preparation of lacto-N-biose I, a candidate for the bifidus factor in human milk	Nishimoto, M.; Kitaoka, M.	<i>Biosci., Biotechnol., Biochem.</i> 71 (8), 2101–2104	2007 Aug



Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Peer Reviewed Journals: HPLC and UHPLC Methods

Carbohydrates

Title	Authors	Publication	Publication Date
Cellotriose and cellotetraose as inducers of the genes encoding cellobiohydrolases in the basidiomycete <i>Phanerochaete chrysosporium</i>	Suzuki, H.; Igarashi, K.; Samejima, M.	<i>Appl. Environ. Microbiol.</i> 76 (18), 6164–6170	2010 Sep
1,2-alpha-L-Fucosynthase: A glycosynthase derived from an inverting alpha-glycosidase with an unusual reaction mechanism	Wada, J.; Honda, Y.; Nagae, M.; Kato, R.; Wakatsuki, S.; Katayama, T.; Taniguchi, H.; Kumagai, H.; Kitaoka, M.; Yamamoto, K.	<i>FEBS Lett.</i> 582 (27), 3739–3743	2008 Nov 12
Efficient separation of oxidized cello-oligosaccharides generated by cellulose degrading lytic polysaccharide monoxygenases	Westereng, B.; Agger, J. W.; Horn, S. J.; Vaaje-Kolstad, G.; Aachmann, F. L.; Stenström, Y. H.; Eijsink, V. G.	<i>J. Chromatogr., A.</i> 1271 (1), 144–152	2013 Jan 4
Distribution of in vitro fermentation ability of lacto-N-Biose I, a major building block of human milk oligosaccharides, in bifidobacterial strains	Xiao, J. Z.; Takahashi, S.; Nishimoto, M.; Odamaki, T.; Yaeshima, T.; Iwatsuki, K.; Kitaoka, M.	<i>Appl. Environ. Microbiol.</i> 76 (1), 54–59	2010 Jan





Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Peer Reviewed Journals: HPLC and UHPLC Methods

Food, Nutrition, Natural Products, and Supplements

Title	Authors	Publication	Publication Date
Characterization of phenolic compounds in strawberry (<i>Fragaria x ananassa</i>) fruits by different HPLC detectors and contribution of individual compounds to total antioxidant capacity	Aaby, K.; Ekeberg, D.; Skrede, G.	<i>J. Agric. Food Chem.</i> 55 (11), 4395–4406	2007 May 30
Analysis of flavonoids and other phenolic compounds using high-performance liquid chromatography with coulometric array detection: relationship to antioxidant activity	Aaby, K.; Hvattum, E.; Skrede, G.	<i>J. Agric. Food Chem.</i> 52 (15), 4595–4603	2004 Jul 28
Aqueous extract of Astragali Radix induces human natriuresis through enhancement of renal response to atrial natriuretic peptide	Ai, P.; Yong, G.; Dingkun, G.; Qiuyu, Z.; Kaiyuan, Z.; Shanyan, L.	<i>J. Ethnopharmacol.</i> 116 (13), 413–421	2008 Mar 28
Antioxidant, α-amylase inhibitory and oxidative DNA damage protective property of <i>Boerhaavia diffusa</i> (Linn.) root	Akhter, F.; Hashim, A.; Khan, M. S.; Ahmad, S.; Iqbal, D.; Srivastava, A. K.; Siddiqui, M. H.	<i>S. Afr. J. Bot.</i> 88, 265–272	2013 Sep
Antioxidant activity and metabolite profile of quercetin in vitamin-E-depleted rats.	Ameho, C. K.; Chen, C. Y. O.; Smith, D.; Sánchez-Moreno, C.; Milbury, P. E.; Blumberg, J. B.	<i>J. Nutr. Biochem.</i> 19 (7), p.467–474	2008 Jul
Evaluation of tolerable levels of dietary quercetin for exerting its antioxidative effect in high cholesterol-fed rats	Azuma, K.; Ippoushi, K.; Terao, J.	<i>Food Chem. Toxicol.</i> 48 (4), 1117–1122	2010 Apr
Recent methodology in ginseng analysis	Baek, S.; Bae, O.; Park, J.	<i>J. Ginseng Res.</i> 36 (2), 119–134	2012 Apr
Sensitive determination of saponins in radix et rhizoma notoginseng by charged aerosol detector coupled with HPLC	Bai, C.; Han, S.; Chai, X.; Jiang, Y.; Li, P.; Tu, P.	<i>J. Liq. Chromatogr. Relat. Technol.</i> 32 (2), 242–260	2010 Aug 27
Comprehensive analysis of polyphenols in 55 extra virgin olive oils by HPLC-ECD and their correlation with antioxidant activities	Bayram, B.; Esatbeyoglu, T.; Schulze, N.; Ozcelik, B.; Frank, J.; Rimbach, G.	<i>Plant Foods Hum. Nutr. (N. Y., NY, U.S.)</i> 67 (4), 326–336	2012 Dec
Hydrogen sulfide mediates the vasoactivity of garlic	Benavides, G. A.; Squadrito, G. L.; Mills, R. W.; Patel, H. D.; Isbell, T. S.; Patel, R. P.; Darley-Usmar, V. M.; Doeller, J. E.; Kraus, D. W.	<i>Proc. Natl. Acad. Sci. U.S.A.</i> 104 (46), 17977–17982	2007 Nov
Analysis of selected stilbenes in <i>Polygonum cuspidatum</i> by HPLC coupled with CoulArray detection	Benová, B.; Adam, M.; Onderková, K.; Královský, J.; Krájček, M.	<i>J. Sep. Sci.</i> 31 (13), 2404–2409	2008 Jul
Rapid and complete extraction of phenols from olive oil and determination by means of a coulometric electrode array system	Brenes, M.; García, A.; García, P.; Garrido, A.	<i>J. Agric. Food Chem.</i> 48 (11), 5178–5183	2000 Nov
The real nature of the indole alkaloids in <i>Cortinarius infractus</i>: Evaluation of artifact formation through solvent extraction method development	Brondz, I.; Ekeberg, D.; Høiland, K.; Bell, D.; Annino, A.	<i>J. Chromatogr., A</i> 1148 (1), 1–7	2007 Apr 27



Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Peer Reviewed Journals: HPLC and UHPLC Methods

Food, Nutrition, Natural Products, and Supplements

Title	Authors	Publication	Publication Date
Chemotaxonomic differentiation between <i>Cortinarius infractus</i> and <i>Cortinarius subtortus</i> by supercritical fluid chromatography connected to a multi-detection system	Brondz, I.; Høiland, K.	<i>Trends Chromatogr.</i> 4, 79–87	2008
Carotenoid bioavailability is higher from salads ingested with full-fat than with fat-reduced salad dressings as measured with electrochemical detection	Brown, M. J.; Ferruzzi, M. G.; Nguyen, M. L.; Cooper, D. A.; Eldridge, A. L.; Schwartz, S. J.; White, W. S.	<i>Am. J. Clin. Nutr.</i> 80 (2), 396–403	2004 Aug
Naringenin from cooked tomato paste is bioavailable in men	Bugianesi, R.; Catasta, G.; Spigno, P.; D'Uva, A.; Maiani, G.	<i>J. Nutr.</i> 132 (11), 3349–3352	2002 Nov
"Dilute-and-shoot" triple parallel mass spectrometry method for analysis of vitamin D and triacylglycerols in dietary supplements	Byrdwell, W. C.	<i>Anal. Bioanal. Chem.</i> 401 (10), 3317–3334	2011 Dec
Human skeletal muscle ascorbate is highly responsive to changes in vitamin C intake and plasma concentrations	Carr, A. C.; Bozonet, S. M.; Pullar, J. M.; Simcock, J. W.; Vissers, M. C.	<i>Am. J. Clin. Nutr.</i> 97 (4), 800–807	2013 Apr
Utilization of RP-HPLC fingerprinting analysis for the identification of diterpene glycosides from <i>Stevia rebaudiana</i>	Chaturvedula, V.; Prakash, I.	<i>Int. J. Res. Phytochem. Pharmacol.</i> 1 (2), 88–92	2011 Jun 9
Acid and alkaline hydrolysis studies of stevioside and rebaudioside A	Chaturvedula, V.; Prakash, I.	<i>J. Appl. Pharm. Sci.</i> 1 (8), 104–108	2011 Oct
Spectral analysis and chemical studies of the sweet constituent, rebaudioside A	Chaturvedula, V.; Prakash, I.	<i>Eur. J. Med. Plants</i> 2 (1), 57–65	2012 Feb
Flavonoids from almond skins are bioavailable and act synergistically with vitamins C and E to enhance hamster and human LDL resistance to oxidation	Chen, C.; Milbury, P. E.; Lapsley, K.; Blumberg, J. B.	<i>J. Nutr.</i> 135 (6), 1366–1373	2005 Jun 1
Photostability of rebaudioside A and stevioside in beverages	Clos, J. F.; Dubois, G. E.; Prakash, I.	<i>J. Agric. Food Chem.</i> 56 (18), 8507–8513	2008 Sep 24
CoulArray electrochemical evaluation of tocopherol and tocotrienol isomers in barley, oat and spelt grains	Colombo, M. L.; Marangon, K.; Bugatti, C.	<i>Nat. Prod. Commun.</i> 4 (2), 251–254	2009 Feb
Composition and stability of phytochemicals in five varieties of black soybeans (<i>Glycine max</i>)	Correa, C. R.; Li, L.; Aldini, G.; Carini, M.; Oliver Chen, C. Y.; Chun, H.; Cho, S.; Park, K.; Russell, R. M.; Blumberg, J. B.; Yeum, K.	<i>Food Chem.</i> 123 (4), 1176–1184	2010 Dec 15
Effect of UV-B light and different cutting styles on antioxidant enhancement of commercial fresh-cut carrot products	Du, W.; Avena-Bustillos, R. J.; Breksa, A. P., III.; McHugh, T. H.	<i>Food Chem.</i> 134 (4), 1862–1869	2012 Oct 15



Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Peer Reviewed Journals: HPLC and UHPLC Methods

Food, Nutrition, Natural Products, and Supplements

Title	Authors	Publication	Publication Date
Phenols, lignans and antioxidant properties of legume and sweet chestnut flours	Durazzo, A.; Turfani, V.; Azzini, E.; Maiani, G.; Carcea, M.	<i>Food Chem.</i> 140 (4), 666–671	2013 Oct 15
alpha-Lipoic acid in dietary supplements: development and comparison of HPLC-CEAD and HPLC-ESI-MS methods	Durrani, A. I.; Schwartz, H.; Schmid, W.; Sontag, G.	<i>J. Pharm. Biomed. Anal.</i> 45 (4), 694–699	2007 Nov 30
Comparison between evaporative light scattering detection and charged aerosol detection for the analysis of saikosaponins	Eom, H. Y.; Park, S. Y.; Kim, M. K.; Suh, J. H.; Yeom, H.; Min, J. W.; Kim, U.; Lee, J.; Youm, J. R.; Han, S. B.	<i>J. Chromatogr., A.</i> 1217 (26), 4347–4354	2010 Jun 25
Assessment of microcystin purity using charged aerosol detection	Edwards, C.; Lawton, L. A.	<i>J. Chromatogr., A.</i> 1217 (32), 5233–5238	2010 Aug 6
Analysis of lycopene geometrical isomers in biological microsamples by liquid chromatography with coulometric array detection	Ferruzzi, M. G.; Nguyen, M. L.; Sander, L. C.; Rock, C. L.; Schwartz, S. J.	<i>J. Chromatogr., B: Biomed. Sci. Appl.</i> 760 (2), 289–299	2001 Sep 5
Charged aerosol detection to characterize components of dispersed-phase formulations	Fox, C. B.; Sivananthan, S. J.; Mikasa, T. J.; Lin, S.; Parker, S. C.	<i>Adv. Colloid Interface Sci.</i> 199–200, 59–65	2013 Nov
HPLC with charged aerosol detection for the measurement of natural products	Fukushima, K.; Kanedai, Y.; Hirose, K.; Matsumoto, T.; Hashiguchi, K.; Senda, M.; et al.	<i>Chromatography 27 (Suppl. 1)</i> , 83–86	2006
Determination of heterocyclic aromatic amines in beef extract, cooked meat and rat urine by liquid chromatography with coulometric electrode array detection	Gerbl, U.; Cichna, M.; Zsivkovits, M.; Knasmüller, S.; Sontag, G.	<i>J. Chromatogr., B: Anal. Technol. Biomed. Life Sci.</i> 802 (1), 107–113	2004 Mar 25
Determination of macrolide antibiotics in porcine and bovine urine by high-performance liquid chromatography coupled to coulometric detection	González de la Huebra, M. J.; Vincent, U.; Bordin, G.; Rodríguez, A. R.	<i>Anal. Bioanal. Chem.</i> 382 (2), 433–439	2005 May
Development and validation of HPLC-DAD-CAD-MS3 method for qualitative and quantitative standardization of polyphenols in <i>Agrimoniae eupatoriæ herba</i> (Ph. Eur)	Granica, S.; Krupa, K.; Klebowska, A.; Kiss, A. K.	<i>J. Pharm. Biomed. Anal.</i> 86, 112–122	2013 Dec
Total reducing capacity of fresh sweet peppers and five different Italian pepper recipes	Greco, L.; Riccio, R.; Bergero, S.; Del Re, A. A. M.; Trevisan, M.	<i>Food Chem.</i> 103 (4), 1127–1133	2007 Jan
Urinary 3-(3,5-dihydroxyphenyl)-1-propanoic acid, an alkylresorcinol metabolite, is a potential biomarker of whole-grain intake in a U.S. population	Guymon, L. A.; Adlercreutz, H.; Koskela, A.; Li, L.; Beresford, S. A.; Lampe, J. W.	<i>J. Nutr.</i> 138 (10), 1957–1962	2008 Oct
Multidimensional LC x LC analysis of phenolic and flavone natural antioxidants with UV-electrochemical coulometric and MS detection	Hájek, T.; Skeríková, V.; Cesla, P.; Vynuchalová, K.; Jandera, P.	<i>J. Sep. Sci.</i> 31 (19), 3309–3328	2008 Oct
Determination of the urinary aglycone metabolites of vitamin K by HPLC with redox-mode electrochemical detection	Harrington, D. J.; Soper, R.; Edwards, C.; Savidge, G. F.; Hodges, S. J.; Shearer, M. J.	<i>J. Lipid Res.</i> 46 (5), 1053–1060	2005 May



Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Peer Reviewed Journals: HPLC and UHPLC Methods

Food, Nutrition, Natural Products, and Supplements

Title	Authors	Publication	Publication Date
Bioavailability and antioxidant effect of epigallocatechin gallate administered in purified form versus as green tea extract in healthy individuals	Henning, S. M.; Niu, Y.; Liu, Y.; Lee, N. H.; Hara, Y.; Thames, G. D.; Minutti, R. R.; Carpenter, C. L.; Wang, H.; Heber, D.	<i>J. Nutr. Biochem.</i> 16 (10), 610–616	2005 Oct
Procyanidin dimer B₂ [epicatechin-(4β-8)-epicatechin] in human plasma after the consumption of a flavanol-rich cocoa	Holt, R. R.; Lazarus, S. A.; Sullards, M. C.; Zhu, Q. Y.; Schramm, D. D.; Hammerstone, J. F.; Fraga, C. G.; Schmitz, H. H.; Keen, C. L.	<i>Am. J. Clin. Nutr.</i> 76 (4), 798–804	2002 Oct
Effects of natural (RRR α-tocopherol acetate) or synthetic (all-rac α-tocopherol acetate) vitamin E supplementation on reproductive efficiency in beef cows	Horn, M.; Gunn, P.; Van Emon, M.; Lemenager, R.; Burgess, J.; Pyatt, N. A.; Lake, S. L.	<i>J. Anim. Sci. (Savoy, IL, U.S.)</i> 88 (9), 3121–3127	2010 Sep
RP-HPLC analysis of phenolic compounds and flavonoids in beverages and plant extracts using a CoulArray detector	Jandera, P.; Skeifíková, V.; Rehová, L.; Hájek, T.; Baldríanová, L.; Skopová, G.; Kellner, V.; Horna, A.	<i>J. Sep. Sci.</i> 28 (9–10), 1005–1022	2005 Jun
A new application of charged aerosol detection in liquid chromatography for the simultaneous determination of polar and less polar ginsenosides in ginseng products	Jia, S.; Li, J.; Yunusova, N.; Park, J. H.; Kwon, S. W.; Lee, J.	<i>Phytochem. Anal.</i> 24 (4), 374–380	2013 Jul–Aug
A combination of aspirin and γ-tocopherol is superior to that of aspirin and α-tocopherol in anti-inflammatory action and attenuation of aspirin-induced adverse effects	Jiang, Q.; Moreland, M.; Ames, B. N.; Yin, X.	<i>J. Nutr. Biochem.</i> 20 (11), 894–900	2009 Nov
HPLC analysis of rosmarinic acid in feed enriched with aerial parts of <i>Prunella vulgaris</i> and its metabolites in pig plasma using dual-channel coulometric detection	Jirovský, D.; Kosina, P.; Myslíňová, M.; Stýskála, J.; Ulrichová, J.; Simánek V.	<i>J. Agric. Food Chem.</i> 55 (19), 7631–7637	2007 Sep 19
Molar absorptivities and reducing capacity of pyranoanthocyanins and other anthocyanins	Jordheim, M.; Aaby, K.; Fossen, T.; Skrede, G.; Andersen, Ø. M.	<i>J. Agric. Food Chem.</i> 55 (26), 10591–10598	2007 Dec 26
Sensitive electrochemical detection method for alpha-acids, beta-acids and xanthohumol in hops (<i>Humulus lupulus</i> L.)	Kac, J.; Vovk, T.	<i>J. Chromatogr., B: Anal. Technol. Biomed. Life Sci.</i> 850 (1–2), 531–537	2007 May 1
Determination of phenolic compounds and hydroxymethylfurfural in meads using high performance liquid chromatography with coulometric-array and UV detection	Kahoun, D.; Rezková, S.; Veskrnová, K.; Králůvský, J.; Holcapek, M.	<i>J. Chromatogr., A</i> 1202 (1), 19–33	2008 Aug 15
Analysis of terpene lactones in a Ginkgo leaf extract by high-performance liquid chromatography using charged aerosol detection	Kakigi, Y.; Mochizuki, N.; Icho, T.; Hakamatsuka, T.; Goda, Y.	<i>Biosci., Biotechnol., Biochem.</i> 74 (3), 590–594	2010
Linear aglycones are the substrates for glycosyltransferase DesVII in methymycin biosynthesis: analysis and implications	Kao, C.; Borisova, S.; Kim, H.; Liu, H.	<i>J. Am. Chem. Soc.</i> 128 (17), 5606–5607	2006 May 3

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



References

Peer Reviewed Journals: HPLC and UHPLC Methods

Food, Nutrition, Natural Products, and Supplements

Title	Authors	Publication	Publication Date
Antioxidant-rich food intakes and their association with blood total antioxidant status and vitamin C and E levels in community-dwelling seniors from the Quebec longitudinal study NuAge	Khalil, A.; Gaudreau, P.; Cherki, M.; Wagner, R.; Tessier, D. M.; Fulop, T.; Shatenstein, B.	<i>Exp. Gerontol.</i> 46 (6), 475–481	2011 Jun
Certification of a pure reference material for the ginsenoside Rg1	Kim, D.; Chang, J.; Sohn, H.; Cho, B.; Ko, S.; Nho, K.; Jang, D.; Lee, S.	<i>Accredit. Qual. Assur.</i> 15 (2), 81–87	2009 Sep
Optimization of pressurized liquid extraction for spicatoside A in <i>Liriope platyphylla</i>	Kim, S. H.; Kim, H. K.; Yang, E. S.; Lee, K. Y.; Kim, S. D.; Kim, Y. C.; Sung, S. H.	<i>Sep. Purif. Technol.</i> 71 (2), 168–172	2010
Production of surfactin and iturin by <i>Bacillus licheniformis</i> N1 responsible for plant disease control activity	Kong, H. G.; Kim, J. C.; Choi, G. J.; Lee, K. Y.; Kim, H. J.; Hwang, E. C.; Moon, B. J.; Lee, S. W.	<i>Plant Pathol. J.</i> 26 (2), 170–177	2010
Transepithelial transport of microbial metabolites of quercetin in intestinal Caco-2 cell monolayers	Konishi, Y.	<i>J. Agric. Food Chem.</i> 53 (3), 601–607	2005 Feb 9
Absorption and bioavailability of artemisin C in rats after oral administration	Konishi, Y.; Hitomi, Y.; Yoshida, M.; Yoshioka, E.	<i>J. Agric. Food Chem.</i> 53 (26), 9928–9933	2005 Dec 28
Pharmacokinetic study of caffeic and rosmarinic acids in rats after oral administration	Konishi, Y.; Hitomi, Y.; Yoshida, M.; Yoshioka, E.	<i>J. Agric. Food Chem.</i> 53 (12), 4740–4746	2005 Jun 15
Intestinal absorption of <i>p</i>-coumaric and gallic acids in rats after oral administration	Konishi, Y.; Hitomi, Y.; Yoshioka, E.	<i>J. Agric. Food Chem.</i> 52 (9), 2527–2532	2004 May 5
Microbial metabolites of ingested caffeic acid are absorbed by the monocarboxylic acid transporter (MCT) in intestinal Caco-2 cell monolayers	Konishi, Y.; Kobayashi, S.	<i>J. Agric. Food Chem.</i> 52 (21), 6418–6424	2004 Oct 20
Transepithelial transport of rosmarinic acid in intestinal Caco-2 cell monolayers	Konishi, Y.; Kobayashi, S.	<i>Biosci., Biotechnol., Biochem.</i> 69 (3), 583–591	2005 Mar
Effects of various doses of selenite on stinging nettle (<i>Urtica dioica</i> L.)	Krystofova, O.; Adam, V.; Babula, P.; Zehnalek, J.; Beklova, M.; Havel, L.; Kizek, R.	<i>Int. J. Environ. Res. Public Health</i> 7 (10), 3804–3815	2010 Oct
Biofortified cassava increases β-carotene and vitamin A concentrations in the TAG-rich plasma layer of American women	La Frano, M. R.; Woodhouse, L. R.; Burnett, D. J.; Burri, B. J.	<i>Br. J. Nutr.</i> 110 (2), 310–320	2013 Jul 28
Chlorogenic acid is absorbed in its intact form in the stomach of rats	Lafay, S.; Gil-Izquierdo, A.; Manach, C.; Morand, C.; Besson, C.; Scalbert, A.	<i>J. Nutr.</i> 136 (5), 1192–1197	2006 May
Determination of 4-ethylcatechol in wine by high-performance liquid chromatography-coulometric electrochemical array detection	Larcher, R.; Nicolini, G.; Bertoldi, D.; Nardin, T.	<i>Anal. Chim. Acta</i> 609 (2), 235–240	2008 Feb 25
Determination of volatile phenols in wine using high-performance liquid chromatography with a coulometric array detector	Larcher, R.; Nicolini, G.; Puecher, C.; Bertoldi, D.; Moser, S.; Favaro, G.	<i>Anal. Chim. Acta</i> 582 (1), 55–60	2007 Jan 16



Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Peer Reviewed Journals: HPLC and UHPLC Methods

Food, Nutrition, Natural Products, and Supplements

Title	Authors	Publication	Publication Date
Acute, quercetin-induced reductions in blood pressure in hypertensive individuals are not secondary to lower plasma angiotensin-converting enzyme activity or endothelin-1: nitric oxide	Larson, A.; Witman, M. A. H.; Guo, Y.; Ives, S.; Richardson, R. S.; Bruno, R. S.; Jalili, T.; Symons, J. D.	<i>Nutr. Res. (N. Y., NY, U.S.)</i> 32 (8), 557–564	2012 Aug
High-performance liquid chromatography method for the determination of folic acid in fortified food products	Lebiedzinska, A.; Dałbrowska, M.; Szefer, P.; Marszałł M.	<i>Toxicol. Mech. Methods</i> 18 (6), 463–467	2008 Jul
Reversed-phase high-performance liquid chromatography method with coulometric electrochemical and ultraviolet detection for the quantification of vitamins B(1) (thiamine), B(6) (pyridoxamine, pyridoxal and pyridoxine) and B(12) in animal and plant foods	Lebiedzinska, A.; Marszałł, M. L.; Kuta, J.; Szefer, P.	<i>J. Chromatogr., A</i> 1173 (1–2), 71–80	2007 Nov 30
An improved method for the determination of green and black tea polyphenols in biomatrices by high-performance liquid chromatography with coulometric array detection	Lee, M. J.; Prabhu, S.; Meng, X.; Li, C.; Yang, C. S.	<i>Anal. Biochem.</i> 279 (2), 164–169	2000 Mar 15
Characterisation, extraction efficiency, stability and antioxidant activity of phytonutrients in <i>Angelica keiskei</i>	Li, L.; Aldini, G.; Carini, M.; Chen, C. Y. O.; Chun, H.; Cho, S.; Park, K.; Correa, C. R.; Russell, R. M.; Blumberg, J. B.; Yeum, K.	<i>Food Chem.</i> 115 (1), 227–232	2009 Jul
Vitamin A equivalence of the β-carotene in β-carotene-biofortified maize porridge consumed by women	Li, S.; Nugroho, A.; Rocheford, T.; White, W. S.	<i>Am. J. Clin. Nutr.</i> 92 (5), 1105–1112	2010 Nov
Phase IIa chemoprevention trial of green tea polyphenols in high-risk individuals of liver cancer: modulation of urinary excretion of green tea polyphenols and 8-hydroxydeoxyguanosine	Luo, H.; Tang, L.; Tang, M.; Billam, M.; Huang, T.; Yu, J.; Wei, Z.; Liang, Y.; Wang, K.; Zhang, Z. Q.; Zhang, L.; Wang, J. S.	<i>Carcinogenesis</i> 27 (2), 262–268	2006 Feb
Determination of four water-soluble compounds in <i>Salvia miltiorrhiza</i> Bunge by high-performance liquid chromatography with a coulometric electrode array system	Ma, L.; Zhang, X.; Guo, H.; Gan, Y.	<i>J. Chromatogr., B: Anal. Technol. Biomed. Life Sci.</i> 833 (2), 260–263	2006 Apr 3
Effect of green tea powder (<i>Camellia sinensis</i> L. cv. Benifuuki) particle size on O-methylated EGCG absorption in rats. The Kakegawa Study	Maeda-Yamamoto, M.; Ema, K.; Tokuda, Y.; Monobe, M.; Tachibana, H.; Sameshima, Y.; Kuriyama, S.	<i>Cytotechnology</i> 63 (2), 171–179	2011 Mar
Supplementation of a γ-tocopherol-rich mixture of tocopherols in healthy men protects against vascular endothelial dysfunction induced by postprandial hyperglycemia	Mah, E.; Noh, S. K.; Ballard, K. D.; Park, H. J.; Volek, J. S.; Bruno, R. S.	<i>J. Nutr. Biochem.</i> 24 (1), 196–203	2013 Jan

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



References

Peer Reviewed Journals: HPLC and UHPLC Methods

Food, Nutrition, Natural Products, and Supplements

Title	Authors	Publication	Publication Date
Mediterranean diet reduces endothelial damage and improves the regenerative capacity of endothelium	Marin, C.; Ramirez, R.; Delgado-Lista, J.; Yubero-Serrano, E. M.; Perez-Martinez, P.; Carracedo, J.; Garcia-Rios, A.; Rodriguez, F.; Gutierrez-Mariscal, F. M.; Gomez, P.; Perez-Jimenez, F.; Lopez-Miranda, J.	<i>Am. J. Clin. Nutr.</i> 93 (2), 267–274	2011 Feb
Photodiode array (PDA) and other detection methods in HPLC of plant metabolites	Markowski, W.; Waksmundzka-Hajnos, M.	Chapter 13 in <i>High Performance Liquid Chromatography in Phytochemical Analysis</i> , Chromatographic Science Series, Markowski, W., Sherma, J., Eds.; Taylor & Francis Group, LLC: Boca Raton, FL; 331–350	2010 Nov
Determination of water-soluble vitamins in infant milk and dietary supplement using a liquid chromatography on-line coupled to a corona-charged aerosol detector	Márquez-Sillero, I.; Cárdenas, S.; Valcárcel, M.	<i>J. Chromatogr., A.</i> 1313C, 253–258	2013 Oct 25
Sensitive high-performance liquid chromatographic method using coulometric electrode array detection for measurement of phytoestrogens in dried blood spots	Melby, M. K.; Watanabe, S.; Whitten, P. L.; Worthman, C. M.	<i>J. Chromatogr., B: Anal. Technol. Biomed. Life Sci.</i> 826 (1–2), 81–90	2005 Nov 5
Phenolic acids from beer are absorbed and extensively metabolized in humans	Nardini, M.; Natella, F.; Scaccini, C.; Ghiselli, A.	<i>J. Nutr. Biochem.</i> 17 (1), 14–22	2006 Jan
High-performance liquid chromatography analysis of plant saponins: An update 2005-2010	Negi, J. S.; Singh, P.; Pant, G. J.; Rawat, M. S.	<i>Pharmacogn. Rev.</i> 5 (10), 155–158	2011 Jul
Physicochemical effect of pH and antioxidants on mono- and triglutamate forms of 5-methyltetrahydrofolate, and evaluation of vitamin stability in human gastric juice: Implications for folate bioavailability	Ng, X.; Lucock, M.; Veysey, M.	<i>Food Chem.</i> 106 (1), 200–210	2008 Jan
Practical preparation of lacto-N-biose I, a candidate for the bifidus factor in human milk	Nishimoto, M.; Kitaoka, M.	<i>Biosci., Biotechnol., Biochem.</i> 71 (8), 2101-2104	2007 Aug
Hydrophilic interaction liquid chromatography—charged aerosol detection as a straightforward solution for simultaneous analysis of ascorbic acid and dehydroascorbic acid	Nováková, L.; Solichová, D.; Solich, P.	<i>J. Chromatogr., A.</i> 1216 (21), 4574–4581	2009 May 22



Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Peer Reviewed Journals: HPLC and UHPLC Methods

Food, Nutrition, Natural Products, and Supplements

Title	Authors	Publication	Publication Date
No effect on adenoma formation in Min mice after moderate amount of flaxseed	Oikarinen, S.; Heinonen, S. M.; Nurmi, T.; Adlercreutz, H.; Mutanen, M.	<i>Eur. J. Nutr.</i> 44 (5), 273–280	2005 Aug
Measurement of isoflavones using liquid chromatography with multi-channel coulometric electrochemical detection	Ouchi, K.; Gamache, P.; Acworth, I.; Watanabe, S.	<i>BioFactors.</i> 22 (1–4), 353–356	2004
Quantitation of clovamide-type phenylpropenoic acid amides in cells and plasma using high-performance liquid chromatography with a coulometric electrochemical detector	Park, J. B.	<i>J. Agric. Food Chem.</i> 53 (21), 8135–8140	2005 Oct 19
Synthesis, HPLC measurement and bioavailability of the phenolic amide amkamide	Park, J. B.	<i>J. Chromatogr. Sci.</i> [Epub ahead of print]	2013 May 27
Synthesis of safflomide and its HPLC measurement in mouse plasma after oral administration	Park, J. B.; Chen, P.	<i>J. Chromatogr., B: Anal. Technol. Biomed. Life Sci.</i> 852 (1–2), 398–402	2007 Jun 1
Determination of lignans in human plasma by liquid chromatography with coulometric electrode array detection	Peñalvo, J. L.; Nurmi, T.; Haajanen, K.; Al-Maharik, N.; Botting, N.; Adlercreutz, H.	<i>Anal. Biochem.</i> 332 (2), 384–393	2004 Sep 15
Supercritical antisolvent fractionation of lignans from the ethanol extract of flaxseed	Perretti, G.; Virgili, C.; Troilo, A.; Marconi, O.; Regnicoli, G. F.; Fantozzi, P.	<i>J. Supercrit. Fluids</i> 75, 94–100	2013 Mar
Analysis of flavonoids in honey by HPLC coupled with coulometric electrode array detection and electrospray ionization mass spectrometry	Petrus, K.; Schwartz, H.; Sontag, G.	<i>Anal. Bioanal. Chem.</i> 400 (8), 2555–2563	2011 Jun
High-dose supplementation with natural α-tocopherol does neither alter the pharmacodynamics of atorvastatin nor its phase I metabolism in guinea pigs	Podszun, M. C.; Grebenstein, N.; Hofmann, U.; Frank, J.	<i>Toxicol. Appl. Pharmacol.</i> 266 (3), 452–458	2013 Feb 1
Application of high-performance liquid chromatography with charged aerosol detection for universal quantitation of undeclared phosphodiesterase-5 inhibitors in herbal dietary supplements	Poplawska, M.; Blazewicz, A.; Bukowska, K.; Fijalek, Z.	<i>J. Pharm. Biomed. Anal.</i> 84, 232–243	2013 Oct
Isolation and analysis of ginseng: advances and challenges	Qi, L.; Wang, C.; Yuan, C.	<i>Nat. Prod. Rep.</i> 28 (3), 467–495	2011 Mar
Folate analysis in complex food matrices: Use of a recombinant Arabidopsis γ-glutamyl hydrolase for folate deglutamylation	Ramos-Parra, P. A.; Urrea-López, R.; Díaz de la Garza, R. I.	<i>Food Res. Int.</i> 54 (1), 177–185	2013 Nov
Optimisation of gradient HPLC analysis of phenolic compounds and flavonoids in beer using a coularray detector	Rehová, L.; Skeríková, V.; Jandera, P.	<i>J. Sep. Sci.</i> 27 (15–16), 1345–1359	2004 Nov
Chiral separation of (+)/(-)-catechin from sulfated and glucuronidated metabolites in human plasma after cocoa consumption	Ritter, C.; Zimmermann, B. F.; Galensa, R.	<i>Anal. Bioanal. Chem.</i> 397 (2), 723–730	2010 May



Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Peer Reviewed Journals: HPLC and UHPLC Methods

Food, Nutrition, Natural Products, and Supplements

Title	Authors	Publication	Publication Date
Analysis of alkylresorcinols in cereal grains and products using ultrahigh-pressure liquid chromatography with fluorescence, ultraviolet, and CoulArray electrochemical detection	Ross, A. B.	<i>J. Agric. Food Chem.</i> 60 (36), 8954–8962	2012 Sep 12
Rapid and sensitive analysis of alkylresorcinols from cereal grains and products using HPLC-CoulArray-based electrochemical detection	Ross, A. B.; Kochhar, S.	<i>J. Agric. Food Chem.</i> 57 (12), 5187–5193	2009 Jun 24
Analysis of soy isoflavone plasma levels using HPLC with coulometric detection in postmenopausal women	Saracino, M. A.; Raggi, M. A.	<i>J. Pharm. Biomed. Anal.</i> 53 (3), 682–687	2010 Nov 2
A biosynthetic pathway for BE-7585A, a 2-thiosugar-containing angucycline-type natural product	Sasaki, E.; Ogasawara, Y.; Liu, H. W.	<i>J. Am. Chem. Soc.</i> 132 (21), 7405–7417	2010 Jun 2
The senescence-accelerated mouse-prone 8 is not a suitable model for the investigation of cardiac inflammation and oxidative stress and their modulation by dietary phytochemicals	Schiborr, C.; Schwamm, D.; Kocher, A.; Rimbach, G.; Eckert, G. P.; Frank, J.	<i>Pharmacol. Res.</i> 74, 113–120	2013 Aug
Comprehensive impurity profiling of nutritional infusion solutions by multidimensional off-line reversed-phase liquid chromatography × hydrophilic interaction chromatography-ion trap mass-spectrometry and charged aerosol detection with universal calibration	Schiesel, S.; Lämmerhofer, M.; Lindner, W.	<i>J. Chromatogr., A.</i> 1259, 100–10	2012 Oct 12
The effect of α-tocopherol supplementation on training-induced elevation of S100B protein in sera of basketball players	Schulpis, K. H.; Moukas, M.; Parthimos, T.; Tsakiris, T.; Parthimos, N.; Tsakiris, S.	<i>Clin. Biochem.</i> 40 (12), 900–906	2007 Aug
Determination of secoisolariciresinol, lariciresinol and isolariciresinol in plant foods by high performance liquid chromatography coupled with coulometric electrode array detection	Schwartz, H.; Sontag, G.	<i>J. Chromatogr., B: Anal. Technol. Biomed. Life Sci.</i> 838 (2), 78–85	2006 Jul 11
Assessment of probiotic strains ability to reduce the bioaccessibility of aflatoxin M₁ in artificially contaminated milk using an in vitro digestive model	Serrano-Niño, J. C.; Cavazos-Garduño, A.; Hernandez-Mendoza, A.; Applegate, B.; Ferruzzi, M. G.; San Martin-González, M. F.; García, H. S.	<i>Food Control</i> 31 (1), 202–207	2013 May
Intestinal uptake of quercetin-3-glucoside in rats involves hydrolysis by lactase phlorizin hydrolase	Sesink, A. L.; Arts, I. C.; Faassen-Peters, M.; Hollman, P. C.	<i>J. Nutr.</i> 133 (3), 773–776	2003 Mar
Quercetin glucuronides but not glucosides are present in human plasma after consumption of quercetin-3-glucoside or quercetin-4'-glucoside	Sesink, A. L.; O'Leary, K. A.; Hollman, P. C.	<i>J. Nutr.</i> 131 (7), 1938–1941	2001 Jul
Co-administration of quercetin and catechin in rats alters their absorption but not their metabolism	Silberberg, M.; Morand, C.; Manach, C.; Scalbert, A.; Remesy, C.	<i>Life Sci.</i> 77 (25), 3156–3167	2005 Nov 4
Nutritional status is altered in the self-neglecting elderly	Smith, S. M.; Mathews Oliver, S. A.; Zwart, S. R.; Kala, G.; Kelly, P. A.; Goodwin, J. S.; Dyer, C. B.	<i>J. Nutr.</i> 136 (10), 2534–2541	2006 Oct



Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Peer Reviewed Journals: HPLC and UHPLC Methods

Food, Nutrition, Natural Products, and Supplements

Title	Authors	Publication	Publication Date
Binding of heterocyclic aromatic amines by lactic acid bacteria: results of a comprehensive screening trial	Stidl, R.; Sontag, G.; Koller, V.; Knasmüller, S.	<i>Mol. Nutr. Food Res.</i> 52 (3), 322–329	2008 Mar
Direct separation and detection of biogenic amines by ion-pair liquid chromatography with chemiluminescent nitrogen detector	Sun, J.; Guo, H. X.; Semin, D.; Cheetham, J.	<i>J. Chromatogr., A.</i> 1218 (29), 4689–4697	2011 Jul 22
Rapid purification method for fumonisin B1 using centrifugal partition chromatography	Szekeres, A.; Lorántfy, L.; Bencsik, O.; Kecskeméti, A.; Szécsi, Á.; Mesterházy, Á.; Vágvölgyi, C.	<i>Food Addit. Contam.</i> 30 (1), 147–155	2013
Determination of coenzyme Q10 in over-the-counter dietary supplements by high-performance liquid chromatography with coulometric detection	Tang, P. H.	<i>J. AOAC Int.</i> 89 (1), 35–39	2006 Jan–Feb
α-Tocopherol supplementation restores the reduction of erythrocyte glucose-6-phosphate dehydrogenase activity induced by forced training	Tsakiris, S.; Reclus, G. J.; Parthimos, T.; Tsakiris, T.; Parthimos, N.; Schulpis, K. H.	<i>Pharmacol. Res.</i> 54 (5), 373–379	2006 Nov
Tissue distribution of isoflavones in ewes after consumption of red clover silage	Urpi-Sarda, M.; Morand, C.; Besson, C.; Kraft, G.; Viala, D.; Scalbert, A.; Besle, J. M.; Manach, C.	<i>Arch. Biochem. Biophys.</i> 476 (2), 205–210	2008 Aug 15
Performance evaluation of charged aerosol and evaporative light scattering detection for the determination of ginsenosides by LC	Wang, L.; He, W. S.; Yan, H. X.; Jiang, Y.; Bi, K. S.; Tu, P. F.	<i>Chromatographia</i> 70 (3–4), 603–608	2009 Aug
Catechins are bioavailable in men and women drinking black tea throughout the day	Warden, B. A.; Smith, L. S.; Beecher, G. R.; Balentine, D. A.; Clevidence, B. A.	<i>J. Nutr.</i> 131 (6), 1731–1737	2001 Jun
Identification and quantification of polyphenol phytoestrogens in foods and human biological fluids	Wilkinson, A. P.; Wähälä, K.; Williamson, G.	<i>J. Chromatogr., B: Anal. Technol. Biomed. Life Sci.</i> 777 (1–2), 93–109	2002 Sep 25
Bioavailability and pharmacokinetics of caffeoylquinic acids and flavonoids after oral administration of Artichoke leaf extracts in humans	Wittemer, S. M.; Ploch, M.; Windeck, T.; Müller, S. C.; Drewelow, B.; Derendorf, H.; Veit, M.	<i>Phytomedicine</i> 12 (1–2), 28–38	2005 Jan
Validated method for the determination of six metabolites derived from artichoke leaf extract in human plasma by high-performance liquid chromatography-coulometric-array detection	Wittemer, S. M.; Veit, M.	<i>J. Chromatogr., B: Anal. Technol. Biomed. Life Sci.</i> 793 (2), 367–375	2003 Aug 15
HPLC in natural product analysis: The detection issue	Wolfender, J. L.	<i>Planta Med.</i> 75 (07), 719–734	2009 Jun
Simultaneous determination of isoflavones and bisphenol A in rat serum by high-performance liquid chromatography coupled with coulometric array detection	Yasuda, S.; Wu, P. S.; Hattori, E.; Tachibana, H.; Yamada, K.	<i>Biosci., Biotechnol., Biochem.</i> 68 (1), 51–58	2004 Jan
Impurities from polypropylene microcentrifuge tubes as a potential source of interference in simultaneous analysis of multiple lipid-soluble antioxidants by HPLC with electrochemical detection	Yen, H. C.; Hsu, Y. T.	<i>Clin. Chem. Lab. Med.</i> 42 (4), 390–395	2004 Apr



Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Peer Reviewed Journals: HPLC and UHPLC Methods

Food, Nutrition, Natural Products, and Supplements

Title	Authors	Publication	Publication Date
Simultaneous determination of triterpenoid saponins from <i>pulsatilla koreana</i> using high performance liquid chromatography coupled with a charged aerosol detector (HPLC-CAD)	Yeom, H.; Suh, J. H.; Youm, J. R.; Han, S. B.	<i>Bull. Korean Chem. Soc.</i> 31 (5), 1159–1164	2010
DPPH radical scavenging activities of 31 flavonoids and phenolic acids and 10 extracts of Chinese materia medica	Yuan, Y.; Chen, C.; Yang, B.; Kusu, F.; Kotani, A.	<i>Zhongguo Zhongyao Zazhi</i> 34 (13), 1695–1700	2009 Jul
Determination of residual clenbuterol in pork meat and liver by HPLC with electrochemical detection	Zhang, X. Z.; Gan, Y. R.; Zhao, F. N.	<i>Yaoxue Xuebao</i> 39 (4), 276–280	2004 Apr
Identification of equol producers in a Japanese population by high-performance liquid chromatography with coulometric array for determining serum isoflavones	Zhao, J. H.; Sun, S. J.; Arai, Y.; Oguma, E.; Yamada, K.; Horiguchi, H.; Kayama, F.	<i>Phytomedicine</i> 13 (5), 304–309	2006 May
Simultaneous sampling of volatile and non-volatile analytes in beer for fast fingerprinting by extractive electrospray ionization mass spectrometry	Zhu, L.; Hu, Z.; Gamez, G.; Law, W. S.; Chen, H.; Yang, S.; Chingin, K.; Balabin, R. M.; Wang, R.; Zhang, T.; Zenobi, R.	<i>Anal. Bioanal. Chem.</i> 398 (1), 405–413	2010 Sep
Comparison of various easy-to-use procedures for extraction of phenols from apricot fruits	Zitka, O.; Sochor, J.; Rop, O.; Skalickova, S.; Sobrova, P.; Zehnalek, J.; Beklova, M.; Krska, B.; Adam, V.; Kizek, R.	<i>Molecules</i> 16 (4), 2914–2936	2011 Apr 4



Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



References

Peer Reviewed Journals: HPLC and UHPLC Methods

Lipids

Title	Authors	Publication	Publication Date
Development of analytical procedures to study changes in the composition of meat phospholipids caused by induced oxidation	Cascone, A.; Eerola, S.; Ritieni, A.; Rizzo, A.	<i>J. Chromatogr., A</i> 1120 (1–2), 211–220	2006 Jul 7
Evaporative light scattering and charged aerosol detector.	Chaminade, P.	Chapter 5. In <i>Hyphenated and Alternative Methods of Detection in Chromatography</i> , Chromatographic Science Series; Shalliker, A., Ed.; Taylor & Francis Group, LLC: Boca Raton, FL.; 145–160	2012
Simple and efficient profiling of phospholipids in phospholipase D-modified soy lecithin by HPLC with charged aerosol detection	Damjanovic, J.; Nakano, H.; Iwasaki, Y.	<i>J. Am. Oil Chem. Soc.</i> 90 (7), 951–957	2013 Jul
Discriminating olive and non-olive oils using HPLC-CAD and chemometrics	de la Mata-Espinosa, P.; Bosque-Sendra, J. M.; Bro, R.; Cuadros-Rodríguez, L.	<i>Anal. Bioanal. Chem.</i> 399 (6), 2083–2092	2011 Feb
Olive oil quantification of edible vegetable oil blends using triacylglycerols chromatographic fingerprints and chemometric tools	de la Mata-Espinosa, P.; Bosque-Sendra, J. M.; Bro, R.; Cuadros-Rodríguez, L.	<i>Talanta</i> 85 (1), 177–182	2011 Jul 15
Quantification of triacylglycerols in olive oils using HPLC-CAD	de la Mata-Espinosa, P.; Bosque-Sendra, J.; Cuadros-Rodríguez, L.	<i>Food Analytical Methods</i> 4 (4), 574–581	2011 Dec
Quantification of pegylated phospholipids decorating polymeric microcapsules of perfluorooctyl bromide by reverse phase HPLC with a charged aerosol detector	Díaz-López, R.; Libong, D.; Tsapis, N.; Fattal, E.; Chaminade, P.	<i>J. Pharm. Biomed. Anal.</i> 48 (3), 702–707	2008 Nov 4
Squalene emulsions for parenteral vaccine and drug delivery	Fox, C. B.	<i>Molecules</i> 14 (9), 3286–3312	2009 Sep 1
Interactions between parenteral lipid emulsions and container surfaces	Gonyon, T.; Tomaso, A.; Kotha, P.; Owen, H.; Patel, D.; Carter, P.; Cronin, J.; Green, J.	<i>PDA J. Pharm. Sci. and Tech.</i> 67 (3), 247–254	2013 May–Jun
Composition analysis of positional isomers of phosphatidylinositol by high-performance liquid chromatography	Iwasaki, Y.; Masayama, A.; Mori, A.; Ikeda, C.; Nakano, H.	<i>J. Chromatogr., A</i> 1216 (32), 6077–6080	2009 Aug 7
Determination of phospholipid and its degradation products in liposomes for injection by HPLC-charged aerosol detection (CAD)	Jiang, Q.; Yang, R.; Mei, X.	<i>Chinese Pharmaceutical Journal (Zhongguo Yaoxue Zazhi, Beijing, China)</i> 42 (23), 1794–1796	2007



Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Peer Reviewed Journals: HPLC and UHPLC Methods

Lipids

Title	Authors	Publication	Publication Date
Rapid quantification of yeast lipid using microwave-assisted total lipid extraction and HPLC-CAD	Khoomrung, S.; Chumnanpuen, P.; Jansa-Ard, S.; Ståhlman, M.; Nookaew, I.; Borén, J.; Nielsen, J.	<i>Anal. Chem.</i> 85 (10), 4912–4919	2013 May 21
A new liquid chromatography method with charge aerosol detector (CAD) for the determination of phospholipid classes. Application to milk phospholipids	Kiełbowicz, G.; Micek, P.; Wawrzenczyk, C.	<i>Talanta</i> 105, 28–33	2013 Feb 15
An LC method for the analysis of phosphatidylcholine hydrolysis products and its application to the monitoring of the acyl migration process	Kiełbowicz, G.; Smuga, D.; Gładkowski, W.; Chojnacka, A.; Wawrzenczyk, C.	<i>Talanta</i> 94, 22–29	2012 May 30
Separation of acylglycerols, FAME and FFA in biodiesel by size exclusion chromatography	Kittirattanapiboon, K.; Krisnangkura, K.	<i>Eur. J. Lipid Sci. Technol.</i> 110 (5), 422–427	2008 Mar 17
Quantitation of triacylglycerols from plant oils using charged aerosol detection with gradient compensation	Lísa, M.; Lynen, F.; Holčápek, M.; Sandra, P.	<i>J. Chromatogr., A.</i> 1176 (1–2), 135–142	2007 Dec 28
Quantitative study of the stratum corneum lipid classes by normal phase liquid chromatography: comparison between two universal detectors	Merle, C.; Laugel, C.; Chaminade, P.; Baillet-Guffroy, A.	<i>J. Liq. Chromatogr. Relat. Technol.</i> 33, 629–644	2010 Mar
The analysis of lipids via HPLC with a charged aerosol detector	Moreau, R. A.	<i>Lipids</i> 41 (7), 727–34	2006 Jul
Lipid analysis via HPLC with a charged aerosol detector	Moreau, R. A.	<i>Lipid Technol.</i> 21 (8–9), 191–194	2009 Oct 23
Extraction and analysis of food lipids	Moreau, R. A.; Winkler-Moser, J. K.	Chapter 6 in <i>Methods of Analysis of Food Components and Additives</i> , Second Edition; Ötles, S., Ed.; Taylor & Francis Group, LLC: Boca Raton, FL.; 115–134	2011 Nov
Aerosol based detectors for the investigation of phospholipid hydrolysis in a pharmaceutical suspension formulation	Nair, L.; Werling, J.	<i>J. Pharm. Biomed. Anal.</i> 49 (1), 95–99	2009 Jan 15
Structure/function relationships of adipose phospholipase A2 containing a cys-his-his catalytic triad	Pang, X. Y.; Cao, J.; Addington, L.; Lovell, S.; Battaile, K. P.; Zhang, Rao, J. L.; Dennis, E. A.; Moise, A. R.	<i>J. Biol. Chem.</i> 287 (42), 35260–35274	2012 Oct 12
Simultaneous assessment of lipid classes and bile acids in human intestinal fluid by solid-phase extraction and HPLC methods	Persson, E.; Löfgren, L.; Hansson, G.; Abrahamsson, B.; Lennernäs, H.; Nilsson, R.	<i>J. Lipid Res.</i> 48 (1), 242–251	2007 Jan



Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Peer Reviewed Journals: HPLC and UHPLC Methods

Lipids

Title	Authors	Publication	Publication Date
The use of charged aerosol detection with HPLC for the measurement of lipids	Plante, M.; Bailey, B.; Acworth, I.	<i>Methods Mol. Biol.</i> (Totowa, NJ, U.S.) 579, 469–482	2009
Comparison between charged aerosol detection and light scattering detection for the analysis of Leishmania membrane phospholipids	Ramos, R. G.; Libong, D.; Rakotomanga, M.; Gaudin, K.; Loiseau, P. M.; Chaminade, P.	<i>J. Chromatogr., A.</i> 1209 (1–2), 88–94	2008 Oct 31
Authentication of geographical origin of palm oil by chromatographic fingerprinting of triacylglycerols and partial least square-discriminant analysis	Ruiz-Samblás, C.; Arrebola-Pascual, C.; Tres, A.; van Ruth, S.; Cuadros-Rodríguez, L.	<i>Talanta.</i> 116, 788–793	2013 Nov 15
Simple and precise detection of lipid compounds present within liposomal formulations using a charged aerosol detector	Schönherr, C.; Touchene, S.; Wilser, G.; Peschka-Süss, R.; Francese, G.	<i>J. Chromatogr., A.</i> 1216 (5), 781–786	2009 Jan 30
Determination of intralumenal individual bile acids by HPLC with charged aerosol detection	Vertzoni, M.; Archontaki, H.; Reppas, C.	<i>J. Lipid Res.</i> 49 (12), 2690–2695	2008 Dec
Neurolipids and the use of a charged aerosol detector	Waraska, J.; Acworth, I.	<i>Am. Biotechnol. Lab.</i> 26 (1), 12–13	2008





Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Technical Collateral: HPLC and UHPLC Methods

Product Number	Technique	Title
AB 119	UV	Rapid Separation of Paclitaxel and Related Compounds in Paclitaxel Injection
AB 134	MS	LC-MS Analysis of Anthocyanins in Bilberry Extract
AB 139	UV	Separation of Schizandrin, Schizandrin A, and Schizandrin B in a Tablet Sample
AB 153	UV	Save the Flavor – Robust Iso- α -Acids Assaying in Beer within Ten Minutes
AB 155	UV	Monitor the Brewing Process with LC-Transformation of Hop alpha-Acids into Beer Iso-alpha-Acids
AN 109	FLD	Determination of Glyphosate by Cation-Exchange Chromatography with Postcolumn Derivatization
AN 156	UV	The Everlasting Paradigm-Keep Beer Tradition or Prevent Beer from a Skunky Off-Flavor?
AN 196	FLD	Determination of Polycyclic Aromatic Hydrocarbons (PAHs) in Edible Oils by Donor-Acceptor Complex Chromatography (DACC)-HPLC with Fluorescent Detection
AN 207	UV	Chromatographic Fingerprinting of <i>Flos Chrysanthema indicis</i> Using HPLC
AN 213	UV/FLD	Determination of Polycyclic Aromatic Hydrocarbons (PAHs) in Tap Water Using on-Line Solid-Phase Extraction Followed by HPLC with UV and Fluorescence Detections
AN 216	UV	Determination of Water- and Fat-Soluble Vitamins in Functional Waters by HPLC with UV-PDA Detection
AN 224	UV	Determination of Melamine in Milk Powder by Reversed-Phase HPLC with UV Detection
AN 232	UV	Determination of Anthraquinones and Stilbenes in Giant Knotweed Rhizome by HPLC with UV Detection
AN 236	UV	Determination of Iodide and Iodate in Seawater and Iodized Table Salt by HPLC-UV Detection
AN 245	UV	Fast Analysis of Dyes in Foods and Beverages
AN 251	UV	Determination of Water- and Fat-Soluble Vitamins in Nutritional Supplements by HPLC with UV Detection
AN 252	UV	HPLC Assay of Water-Soluble Vitamins, Fat-Soluble Vitamins, and a Preservative in Dry Syrup Multivitamin Formulation
AN 261	UV	Sensitive Determination of Microcystins in Drinking and Environmental Waters
AN 264	UV	Fast Determination of Anthocyanins in Pomegranate Juice
AN 266	FLD	Determination of Sialic Acids Using UHPLC with Fluorescence Detection
AN 272	FLD	Faster Yet Sensitive Determination of N-Methylcarbamates in Rice, Potato, and Corn by HPLC
AN 275	UV	Sensitive Determination of Catechins in Tea by HPLC
AN 287	UV	Two-Dimensional HPLC Combined with On-Line SPE for Determination of Sudan Dyes I-IV in Chili Oil
AN 292	UV	Determination of Aniline and Nitroanilines in Environmental and Drinking Waters by On-Line SPE
AN 293	CAD and UV	Steviol Glycoside Determination by HPLC with Charged Aerosol and UV Detections Using the Acclaim Trinity P1 Column
AN 299	UV	HPLC Analysis of Six Active Components of <i>Caulis Ionicerae</i> Using a Phenyl-1 Column
AN 1008	UV	Determination of Nitidine Chloride, Toddalolactone, and Chelerythrine Chloride by HPLC

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



References

Technical Collateral: HPLC and UHPLC Methods

Product Number	Technique	Title
AN 1020	EC, UV	Chalcinoids and Bitter Acids in Beer by HPLC with UV and ECD
AN 1023	UV	Determination of Sudan Dyes I-IV in Curry Paste
AN 1026	CAD	Fatty Acid Esters at Low Nanogram Levels
AN 1027	CAD	Ginseng
AN 1028	CAD	Ginkgo biloba
AN 1029	CAD	Black Cohosh
AN 1030	CAD	Soy Saponins
AN 1032	CAD	Unsaturated Fatty Acid: Arachidonic, Linoleic, Linolenic and Oleic Acids
AN 1033	CAD	Corn Syrup
AN 1034	CAD	Honey Sugars
AN 1035	CAD	Phenolic Acids
AN 1036	CAD	Water-Soluble Antioxidants: Ascorbic Acid, Glutathione and Uric Acid
AN 1037	CAD	Artificial Sweeteners-Global Method
AN 1039	CAD	Simultaneous Measurement of Glycerides (Mono-, Di- and Triglycerides) and Free Fatty Acids in Palm Oil
AN 1040	CAD	Analysis of Commercially Available Products Containing Stevia
AN 1041	CAD	Phytosterols
AN 1042	UV	Rapid Separation of Anthocyanins in Cranberry and Bilberry Extracts Using a Core-Shell Particle Column
AN 1045	UV	Determination of Phthalates in Drinking Water by UHPLC with UV Detection
AN 1046	UV	Determination of Phenylurea Compounds in Tap Water and Bottled Green Tea
AN 1055	CAD	Determination of Virginiamycin, Erythromycin, and Penicillin in Dried Distillers Grains with Solubles
AN 1063	ECD	Targeted Analyses of Secondary Metabolites in Herbs, Spices, and Beverages Using a Novel Spectro-Electro Array Platform
AN 1064	ECD	Product Authentication and Adulteration Determination Using a Novel Spectro-Electro Array Platform
AN 1067	UV	Determination of Carbendazim in Orange Juice
AN 1069	UV	Two-Dimensional HPLC Determination of Water-Soluble Vitamins in a Nutritional Drink
AN 1070	UV	Determination of Inositol Phosphates in Dried Distillers Grains and Solubles
AN 20583	UV	Determination of Catechins and Phenolic Acids in Red Wine by Solid Phase Extraction and HPLC
AN 20610	UV	Fast Analysis of Coffee Bean Extracts Using a Solid Core HPLC Column
AN 20663	CAD	Comparative Analysis of Cooking Oils Using a Solid Core HPLC Column
AN 20847	CAD	Analysis of a Sports Beverage for Electrolytes and Sugars Using Multi-Mode Chromatography with Charged Aerosol Detection



Technical Collateral: HPLC and UHPLC Methods

Product Number	Technique	Title
AN 70158	CAD	Novel Universal Approach for the Measurement of Natural Products in a Variety of Botanicals and Supplements
AN 70277	CAD	Simultaneous Analysis of Glycerides and Fatty Acids in Palm Oil
AU 144	UV	Determination of Hexavalent Chromium in Drinking Water Using Ion Chromatography
AU 170	UV	Fast Determination of Vanillin and its Synthesis Precursor by HPLC
AU 182	CAD	Measuring Lactose in Milk: A Validated Method
AU 184	CAD, UV	Mogroside V Determination by HPLC with Charged Aerosol and UV Detections
CAN 106	UV	Determination of the Punicalagins Found in Pomegranate by High Performance Liquid Chromatography
CAN 111	CAD	Determination of Triterpenes in <i>Centella asiatica</i> (Gotu Kola) by HPLC-CAD
CAN 112	CAD	Determination of Ginsenosides in Panax ginseng by HPLC-CAD
CAN 115	FLD	Clean-Up and Analysis of Aflatoxins and Ochratoxin A in Herbs and Spices
LPN 2062	MS	Profiling Analysis of 15 Prominent Naturally Occurring Phenolic Acids by LC-MS
LPN 2069	FLD	Fast and Effective Determination of Aflatoxins in Grains or Food Using Accelerated Solvent Extraction followed by HPLC
LPN 2421	UV	Achieving Maximum Productivity by Combining UHPLC with Advanced Chromatographic Techniques
LPN 2818	CAD	Analysis of Fat-Soluble Vitamins and Antioxidants in Supplements by RP-HPLC
LPN 2870	FLD	Benefits of High-Speed Wavelength Switching in UHPLC Methods Using Fluorescence Detection
LPN 2930	CAD	Determination of the Composition of Natural Products by HPLC with Charged Aerosol Detection
LPN 2923	CAD	Simple and Direct Analysis of Falcarinol and Other Polyacetylenic Oxylipins in Carrots by Reversed-Phase HPLC and Charged Aerosol Detection
LPN 2931	CAD	Quantification of Underivatized Omega-3 and Omega-6 Fatty Acids in Foods by HPLC CAD
LPN 2932	ECD	A Versatile Detector for the Sensitive and Selective Measurement of Numerous Fat-Soluble Vitamins and Antioxidants in Human Plasma and Plant Extracts
LPN 2934	CAD	Sensitive Analysis of Commonly Used Artificial and Natural Sweeteners Including Stevia and Their Impurities and Degradation Products
LPN 2991	CAD	Evaluation of Methods for the Characterization and Quantification of Polysorbates and Impurities Along with Other Surfactants and Emulsifiers Used in the Food and Pharmaceutical Industries
PN 70026	CAD	Carbohydrate Analysis Using PAD, FLD, CAD and MS Detectors
PN 70037	CAD	Sensitive HPLC Method for Triterpenoid Analysis Using Charged Aerosol Detection with Improved Resolution
PN 70055	CAD	Direct Analysis of Surfactants using HPLC with Charged Aerosol Detection
PN 70138	UV	Rapid Determination of Polyphenol Antioxidants in Green Tea and Cranberry Extract Using Core Shell Columns
PN 70538	CAD	Analysis of Silicone Oils by HPLC-CAD
PN 70540	CAD, ECD	Profiling <i>Hoodia</i> Extracts by HPLC with CAD, ECD, Principal Component Analysis

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Table of Contents

Introduction

Analytical Technologies

Artificial Colorants

Blue Dyes

Caramel

General Dye Method

Global Method

Red Dyes

Sudan Dyes

Yellow Dyes

Natural Colorants

Turmeric

Flavorants

Amadori Compounds

Capsaicin

Quinine

Vanilla

Miscellaneous

Caffeine

Choline

Glucosamine

Iodide

Polyphosphates

Preservatives

Benzoate

Nitrate and Nitrite

Sulfites

Soft Drinks

References



References



Ion Chromatography References



Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Technical Collateral: Ion Chromatography Methods

Product Number	Technique	Title
AB 127	IC-PAD	Determination of Carbohydrates in Fruit Juice Using Capillary High-Performance Anion-Exchange Chromatography
AB 135	IC-SC	Determination of Anions and Organic Acids in Brewed Coffee Samples Using Capillary IC
AB 137	IC-SC	Determination of Inorganic and Organic Acids in Apple and Orange Juice Samples Using Capillary IC
AN 25	IC-SC	Determination of Inorganic Ions and Organic Acids in Non-Alcoholic Carbonated Beverages
AN 37	IC-PAD	Determination of Iodide and Iodate in Soy- and Mil-Based Infant Formulas
AN 46	IC-PAD	Ion Chromatography: A Versatile Technique for the Analysis of Beer
AN 54	IC-PAD	Determination of Total and Free Sulfite in Foods and Beverages
AN 67	IC-PAD	Determination of Plant-Derived Neutral Oligo- and Polysaccharides
AN 81	IC-SC	Ion Chromatographic Determination of Oxyhalides and Bromide at Trace Level Concentrations in Drinking Water Using direct Injection
AN 82	IC-PAD	Analysis of Fruit Juice Adulterated with Medium Invert Sugar from Beets
AN 87	IC-PAD	Determination of Sugar Alcohols in Confections and Fruit Juices by High-Performance Anion-Exchange Chromatography with Pulsed Amperometric Detection
AN 101	IC-SC	Trace Level Determination of Bromate in Ozonated Drinking Water Using Ion Chromatography
AN 112	IC-UV	Determination of Nitrate and Nitrite in Meat Using High-Performance Anion-Exchange Chromatography
AN 121	IC-SC	Analysis of Low Concentrations of Perchlorate in Drinking Water and Ground Water by Ion Chromatography
AN 123	IC-SC	Determination of Inorganic Anions and Organic Acids in Fermentation Broths
AN 133	IC-SC	Determination of Inorganic Anions in Drinking Water by Ion Chromatography
AN 136	IC-SC and IC-UV	Determination of Inorganic Oxyhalide Disinfection Byproduct Anions and Bromide in Drinking Water Using Ion Chromatography with the Addition of a Postcolumn Reagent for Trace Bromate Analysis
AN 140	IC-SC	Fast Analysis of Anions in Drinking Water by Ion Chromatography
AN 143	IC-SC	Determination of Organic Acids in Fruit Juices
AN 149	IC-SC	Determination of Chlorite, Bromate, Bromide, and Chlorate in Drinking Water by Ion Chromatography with an On-Line-Generated Postcolumn Reagent for Sub- $\mu\text{g/L}$ Bromate Analysis
AN 150	IC-PAD	Determination of Amino Acids in Cell Cultures and Fermentation Broths
AN 154	IC-SC	Determination of Inorganic Anions in Environmental Waters Using a Hydroxide-Selective Column
AN 155	IC-PAD	Determination of Trans-Galactooligosaccharides in Foods by AOAC Method 2001.02

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



References

Technical Collateral: Ion Chromatography Methods

Product Number	Technique	Title
AN 165	IC-SC	Determination of Benzoate in Liquid Food Products by Reagent-Free Ion Chromatography
AN 167	IC-SC	Determination of Trace Concentrations of Oxyhalides and Bromide in Municipal and Bottled Waters Using a Hydroxide-Selective Column with a Reagent-Free Ion Chromatography System
AN 168	IC-UV	Determination of Trace Concentrations of Disinfection By-Product Anions and Bromide in Drinking Water Using Reagent-Free Ion Chromatography Followed by Postcolumn Addition of Iol-Dianisidine for Trace Bromate Analysis
AN 169	IC-SC	Rapid Determination of Phosphate and Citrate in Carbonated Soft Drinks Using a Reagent-Free Ion Chromatography System
AN 172	IC-SC	Determination of Azide in Aqueous Samples by Ion Chromatography with Suppressed Conductivity Detection
AN 173	IC-PAD	Direct Determination of Cyanide in Drinking Water by Ion Chromatography with Pulsed Amperometric Detection (PAD)
AN 178	IC-SC	Improved Determination of Trace Concentrations of Perchlorate in Drinking Water Using Preconcentration with Two-Dimensional Ion Chromatography and Suppressed Conductivity Detection
AN 182	IC-SC and IC-PAD	Determination of Biogenic Amines in Alcoholic Beverages by Ion Chromatography with Suppressed Conductivity and Integrated Pulsed Amperometric Detections
AN 183	IC-SC and IC-PAD	Determination of Biogenic Amines in Fermented and Non-Fermented Foods Using Ion Chromatography with Suppressed Conductivity and Integrated Pulsed Amperometric Detections
AN 187	IC-SC	Determination of sub-µg/L Bromate in Municipal Waters Using Preconcentration with Two-Dimensional Ion Chromatography and Suppressed Conductivity Detection
AN1 88	IC-PAD	Determination of Glycols and Alcohols in Fermentation Broths Using Ion-Exclusion Chromatography and Pulsed Amperometric Detection
AN 197	IC-PAD	Determination of Glucosamine in Dietary Supplements Using HPAE-PAD
AN 227	ICE-PAD	Determination of Total Cyanide in Municipal Wastewater and Drinking Water Using Ion-Exclusion Chromatography with Pulsed Amperometric Detection (ICE-PAD)
AN 248	IC-PAD	Determination of Lactose in Lactose-Free Milk Products by High-Performance Anion-Exchange Chromatography with Pulsed Amperometric Detection
AN 253	IC-PAD	HPAE-PAD Determination of Infant Formula Sialic Acids
AN 270	IC-PAD	Determination of Hydroxymethylfurfural in Honey and Biomass
AN 273	IC-SC	Determination of Organic Acids in Fruit Juices and Wines by High-Pressure IC
AN 279	IC-SC	Time Savings and Improved Reproducibility of Nitrate and Nitrite Ion Chromatography Determination in Milk Samples
AN 280	IC-PAD	Carbohydrates in Coffee: AOAC Method 995.13 vs a New Fast Ion Chromatography Method
AN 295	IC-SC	Determination of Phytic Acid in Soybeans and Black Sesame Seeds
AN 1007	IC-SC	Determination of Mono-, Di-, and Triphosphates and Citrate in Shrimp by Ion Chromatography



Technical Collateral: Ion Chromatography Methods

Product Number	Technique	Title
AN 1044	IC-SC	Determination of Anions in Dried Distillers Grains with Solubles
AN 1068	IC-SC	Determination of Organic Acids in Fruit Juices and Wines by High-Pressure IC
AU 132	IC-UV	Determination of Nitrite and Nitrate in drinking Water by Ion Chromatography with Direct UV Detection
AU 144	IC-UV	Determination of Hexavalent Chromium in Drinking Water Using Ion Chromatography
AU 148	IC-SC	Determination of Perchlorate in Drinking Water Using Reagent-Free Ion Chromatography
AU 150	IC-PAD	Determination of Plant-Derived Neutral Oligo- and Polysaccharides Using the CarboPac PA200
AU 151	IC-PAD	Determination of Sucralose in Reduced- Carbohydrate Colas using High-Performance Anion-Exchange Chromatography with Pulsed Amperometric Detection
AU 189	IC-SC	Determination of Choline in Infant Formula and Other Food Samples by IC
LPN 2982	IC-SC	Determination of Inorganic Anions and Organic Acids in Beverages Using a Capillary IC on a Monolith Anion-Exchange Column
PN 70743	IC-SC	Determination of Perchlorate Levels in Food and Soil Samples Using Accelerated Solvent Extraction and Ion Chromatography
TN 20	IC-PAD	Analysis of Carbohydrates by High-Performance Anion-Exchange Chromatography with Pulsed Amperometric Detection (HPAE-PAD)
TN 126	IC-SC	Determination of Organic Acids in Beer Samples Using a High-Pressure Ion Chromatography System
TN 135	IC-PAD	Determinations of Monosaccharides and Disaccharides in Beverages by Capillary HPAE-PAD

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Table of Contents

Introduction

Analytical Technologies

Artificial Colorants

Blue Dyes

Caramel

General Dye Method

Global Method

Red Dyes

Sudan Dyes

Yellow Dyes

Natural Colorants

Turmeric

Flavorants

Amadori Compounds

Capsaicin

Quinine

Vanilla

Miscellaneous

Caffeine

Choline

Glucosamine

Iodide

Polyphosphates

Preservatives

Benzoate

Nitrate and Nitrite

Sulfites

Soft Drinks

References



References



Sample Preparation References

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



References

Peer Reviewed Journals: Sample Preparation Methods

Title	Authors	Publication	Publication Date
Accelerated, microwave-assisted, and conventional solvent extraction methods affect anthocyanin composition from colored grains	Abdel-Aal el-SM; Akhtar, H.; Rabalski, I.; Bryan, M.	<i>J. Food Sci.</i> 79 (2), C138–46	2014 Feb
Multiresidue method for the analysis of pesticide residues in fruits and vegetables by accelerated solvent extraction and capillary gas chromatography	Adou, K.; Bontoyan, W. R.; Sweeney, P. J.	<i>J. Agric. Food Chem.</i> 49 (9), 4153–4160	2001 Sep
The development of an optimized sample preparation for trace level detection of 17α-ethinylestradiol and estrone in whole fish tissue	Al-Ansari, A. M.; Saleem, A.; Kimpe, L. E.; Trudeau, V. L.; Blais, J. M.	<i>J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.</i> 879 (30), 3649–52	2011 Nov
Determination of polyphenolic profiles of basque cider apple varieties using accelerated solvent extraction	Alonso-Salces, R. M.; Korta, E.; Barranco, A.; Berrueta, L.A.; Gallo, B.; Vicent, F.	<i>J. Agric. Food Chem.</i> 49 (8), 3761–376	2001
Pressurized liquid extraction for the determination of polyphenols in apple	Alonso-Salces, R. M.; Korta, E.; Barranco, A.; Berrueta, L. A.; Gallo, B.; Vicente, F.;	<i>J. Chromatogr., A.</i> 933 (1–2), 37–43	2001 Nov
Methods for extraction and determination of phenolic acids in medicinal plants: a review	Arceusz, A.; Wesolowski, M.; Konieczynski, P.	<i>Nat. Prod. Commun.</i> 8 (12), 1821–9	2013 Dec
Study of an accelerated solvent extraction procedure for the determination of acaricide residues in honey by high-performance liquid chromatography-diode array detector	Bakkali, A.; Korta, E.; Berrueta, L. A.	<i>J. Food Protection</i> 65 (1), 161–166	2002
Pressurized liquid extraction of medicinal plants	Benthin, B.; Danz, H.; Hamburger, M.	<i>J. Chromatogr., A.</i> 837 (1-2), 211–9	1999 Apr
Comparison of the chemical composition of extracts from <i>Scutellaria lateriflora</i> using accelerated solvent extraction and supercritical fluid extraction versus standard hot water or 70% ethanol extraction	Bergeron, C.; Gafner, S.; Clausen, E.; Carrier, D. J.	<i>J. Agric. Food Chem.</i> 53 (8), 3076–80	2005 Apr
Polybrominated diphenyl ethers (PBDEs) in Mediterranean mussels (<i>Mytilus gallo-provincialis</i>) from selected Apulia coastal sites evaluated by GC-HRMS	Bianco, G.; Novario, G.; Anzilotta, G.; Palma, A.; Mangone, A.; Cataldi, T. R.	<i>J. Mass Spectrom.</i> 45 (9), 1046–55	2010 Sep
Free and bound phenolic compounds in barley (<i>Hordeum vulgare</i> L.) flours. evaluation of the extraction capability of different solvent mixtures and pressurized liquid methods by micellar electrokinetic chromatography and spectrophotometry	Bonoli, M.; Marconi, E.; Caboni, M. F.	<i>J. Chromatogr., A.</i> 19; 1057 (1-2), 1–12	2004 Nov
Pressurized liquid extraction of lipids for the determination of oxysterols in egg-containing food	Boselli, E.; Velazco, V.; Caboni, M. F.; Lercker, G.	<i>J. Chromatogr., A.</i> 11; 917 (1-2), 239–44	2001 May
Optimisation of accelerated solvent extraction of cocaine and benzoylecgonine from coca leaves	Brachet, A.; Rudaz, S.; Mateus, L.; Christen, P.; Veuthey, J-L.	<i>J. Sep. Sci.</i> 24 (10-11), 865–873	2001 Nov

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



References

Peer Reviewed Journals: Sample Preparation Methods

Title	Authors	Publication	Publication Date
Multi-residue determination of 130 multiclass pesticides in fruits and vegetables by gas chromatography coupled to triple quadrupole tandem mass spectrometry	Cervera, M.I.; Medina, C.; Portolés, T.; Pitarch, E; Beltrán, J.; Serrahima, E.; Pineda, L.; Muñoz, G.; Centrich, F.; Hernández, F.	<i>Anal. Bioanal. Chem.</i> 397 (7), 2873–91	2010 Aug
Influence of extraction methodologies on the analysis of five major volatile aromatic compounds of citronella grass (<i>Cymbopogon nardus</i>) and lemongrass (<i>Cymbopogon citratus</i>) grown in Thailand	Chanthai, S.; Prachakoll, S.; Ruangviriyachai, C.; Luthria, D. L.	<i>J. AOAC Int.</i> 95 (3), 763–72	2012 May-Jun
Accelerated solvent extraction of vitamin K₁ in medical foods in conjunction with matrix solid-phase dispersion	Chase, G. W.; Thompson, B.	<i>J. AOAC Int.</i> 83 (2), 407–10	2000
Development of a liquid chromatography-tandem mass spectrometry with pressurized liquid extraction method for the determination of benzimidazole residues in edible tissues	Chen, D.; Tao, Y.; Zhang, H.; Pan, Y.; Liu, Z.; Huang, L.; Wang, Y.; Peng, D.; Wang, X.; Dai, M.; Yuan, Z.	<i>J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.</i> 879 (19), 1659–67	2011 Jun
Determination of 88 pesticide residues in tea using gas chromatography-tandem mass spectrometry	Chen, H.; Liu, X.; Wang, Q.; Jiang, Y.	<i>Se Pu.</i> 29 (5), 409–16	2011 May
Optimization of accelerated solvent extraction for the determination of chlorinated pesticides from animal feed	Chen, S.; Gfrerer, M.; Lankmayr, E.; Quan, X.; Yang, F.	<i>Chromatographia</i> 58, 631–636	2003
Uptake of oxytetracycline, sulfamethoxazole and ketoconazole from fertilised soils by plants	Chitescu, C. L.; Nicolau, A. I.; Stolker, A. A.	<i>Food Addit. Contam. Part A Chem. Anal. Control Expo. Risk Assess.</i> 30 (6), 1138–46	2013
Ultrasonic or accelerated solvent extraction followed by U-HPLC-high mass accuracy MS for screening of pharmaceuticals and fungicides in soil and plant samples	Chitescu, C. L.; Oosterink, E.; de Jong, J.; Stolker, A. A.	<i>Talanta</i> 2012; 88, 653–62	2011 Jan
Evaluation of analytical methods for determining pesticides in baby foods and adult duplicate-diet samples	Chuang, J. C.; Hart, K.; Chang, J. S.; Boman, L. E.; Van Emon, J. M.; Reed, A. W.	<i>Anal. Chim. Acta.</i> 444 (1), 87–95	2001 Oct
Comparison of extraction techniques and modeling of accelerated solvent extraction for the authentication of natural vanilla flavors	Cicchetti, E.; Chaintreau, A.	<i>J. Sep. Sci.</i> 32 (11), 1957–64	2009 Jun
Development of a fast and convenient method for the isolation of triterpene saponins from <i>Actaea racemosa</i> by high-speed countercurrent chromatography coupled with evaporative light scattering detection	Cicek, S. S.; Schwaiger, S.; Ellmerer, E. P.; Stuppner, H.	<i>Planta. Med.</i> 76 (5), 467–73	2010 Mar
Extraction of bitter acids from hops and hop products using pressurized solvent extraction (PSE)	Culik, J.; Jurková, M.; Horák, T.; Cejka, P.; Kellner, V.; Dvorák, J.; Karásek, P.; Roth, M.	<i>J. Inst. Brew.</i> 115 (3), 220–225	2009
Comparison of methods for extraction of flavanones and xanthenes from the root bark of the osage orange tree using liquid chromatography	da Costa, C. T.; Margolis, S. A.; Benner, Jr. B.A.; Horton, D.	<i>J. Chromatogr., A.</i> 831 (2), 167–178	1999 Jan
Pressurized liquid extraction prior to liquid chromatography with electrochemical detection for the analysis of vitamin E isomers in seeds and nuts	Delgado-Zamarreño, M. M.; Bustamante-Rangel, M.; Sánchez-Pérez, A.; Carabias-Martínez, R.	<i>J. Chromatogr., A.</i> 12; 1056 (1-2), 249–52	2004 Nov

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



References

Peer Reviewed Journals: Sample Preparation Methods

Title	Authors	Publication	Publication Date
Pressurized fluid extraction of carotenoids from <i>Haematococcus pluvialis</i> and <i>Dunaliella salina</i> and kavalactones from <i>Piper methysticum</i>	Denery, J. R.; Dragull, K.; Tang, C. S.; Li, Q. X.	<i>Anal. Chim. Acta.</i> 501 (2), 175–181	2004 Jan
Development and comparison of two multiresidue methods for the analysis of 17 mycotoxins in cereals by liquid chromatography electrospray ionization tandem mass spectrometry	Desmarchelier, A.; Oberson, J. M.; Tella, P.; Gremaud, E.; Seefelder, W.; Mottier, P.	<i>J. Agric. Food Chem.</i> 58 (13), 7510–9	2010 Jul
Identification, extraction and quantification of the synthetic cannabinoid JWH-018 from commercially available herbal marijuana alternatives	Dunham, S. J.; Hooker, P. D.; Hyde, R. M.	<i>Forensic Sci. Int.</i> 223 (1-3), 241–4	2012 Nov
Evaluation of polyphenol contents in differently processed apricots using accelerated solvent extraction followed by high-performance liquid chromatography-diode array detector	Erdogan, S.; Erdemoglu, S.	<i>Int. J. Food Sci. Nutr.</i> 62 (7), 729–39	2011 Nov
Determination of 2,4,6-trichloroanisole and guaiacol in cork stoppers by pressurised fluid extraction and gas chromatography–mass spectrometry	Ezquerro, Ó.; Garrido-López, Á.; Tena, M. T.	<i>J. Chromatogr., A.</i> 1102 (12), 18–24	2006 Jan
Multiwalled carbon nanotubes as matrix solid-phase dispersion extraction absorbents to determine 31 pesticides in agriculture samples by gas chromatography-mass spectrometry	Fang, G.; Min, G.; He, J.; Zhang, C.; Qian, K.; Wang, S.	<i>J. Agric. Food Chem.</i> 57 (8), 3040–5	2009 Apr
High-anthocyanin strawberries through cultivar selection	Fredericks, C. H.; Fanning, K. J.; Gidley, M. J.; Netzel, G.; Zabar, D.; Herrington, M.; Netzel, M.	<i>J. Sci. Food Agric.</i> 93 (4), 846–52	2013 Mar
Optimal extraction and fingerprint analysis of <i>Cnidii fructus</i> by accelerated solvent extraction and high performance liquid chromatographic analysis with photodiode array and mass spectrometry detections	Gao, F.; Hu, Y.; Ye, X.; Li, J.; Chen, Z.; Fan, G.	<i>Food Chem.</i> 141 (3), 1962–71	2013 Dec
Simultaneous analysis of seven alkaloids in <i>Coptis-evodia</i> herb couple and Zuojin pill by UPLC with accelerated solvent extraction	Gao, X.; Yang, X. W.; Marriott, P. J.	<i>J. Sep. Sci.</i> 33 (17-18), 2714–22	2010 Sep
Determination of chromones in <i>Dysophylla stellata</i> by HPLC: method development, validation and comparison of different extraction methods	Gautam, R.; Srivastava, A.; Jachak, S. M.	<i>Nat. Prod. Commun.</i> 5 (4), 555–8	2010 Apr
Comparison of different extraction techniques for the determination of chlorinated pesticides in animal feed	Gfrerer, M.; Chen, S.; Lankmayr, E.; Xie, Q.; Yang, F.	<i>Anal. Bioanal. Chem.</i> 378 (7), 1861–1867	2004
Speciation analysis of selenium compounds in yeasts using pressurised liquid extraction and liquid chromatography–microwave-assisted digestion–hydride generation–atomic fluorescence spectrometry	Gómez-Ariza, J. L.; Caro de la Torre, M. A.; Giráldez, I.; Morales, E.	<i>Anal. Chim. Acta.</i> 524, (1–2), 305–314	2004 Oct

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



References

Peer Reviewed Journals: Sample Preparation Methods

Title	Authors	Publication	Publication Date
Multianalysis of 35 mycotoxins in traditional Chinese medicines by ultra-high-performance liquid chromatography-tandem mass spectrometry coupled with accelerated solvent extraction	Han, Z.; Ren Y.; Zhu, J.; Cai, Z.; Chen, Y.; Luan, L.; Wu, Y.	<i>J. Agric. Food Chem.</i> 60 (33), 8233-47.	2012 Aug
Pressurized liquid extraction-capillary electrophoresis-mass spectrometry for the analysis of polar antioxidants in rosemary extracts	Herrero, M.; Arráez-Román, D.; Segura A.; Kennler, E.; Gius, B.; Raggid, M. A.; Ibáñez, E.; Cifuentes, A.	<i>J. Chromatogr., A.</i> 1084 (1-2), 54-62.	2005 Aug
Accelerated solvent extraction of alkylresorcinols in food products containing uncooked and cooked wheat	Holt, M D.; Moreau, R A.; DerMarderosian, A.; McKeown, N.; Jacques, P. F.	<i>J. Agric. Food Chem.</i> 60 (19), 4799-802	2012 May
Application of response surface methodology to optimize pressurized liquid extraction of antioxidant compounds from sage (<i>Salvia officinalis</i> L.), basil (<i>Ocimum basilicum</i> L.) and thyme (<i>Thymus vulgaris</i> L.)	Hossain, M. B.; Brunton, N. P.; Martin-Diana, A. B.; Barry-Ryan, C.	<i>Food Funct.</i> 1(3), 269-77	2010 Dec
A review of modern sample-preparation techniques for the extraction and analysis of medicinal plants	Huie, C. W.	<i>Anal. Bioanal. Chem.</i> 373 (1-2), 23-30.	2002 May
Polychlorinated dioxins, furans, and biphenyls, and polybrominated diphenyl ethers in a U.S. meat market basket and estimates of dietary intake	Huwe, J. K.; Larsen, G. L.	<i>Environ. Sci. Technol.</i> 39 (15), 5606-5611	2005
Study of the effect of sample preparation and cooking on the selenium speciation of selenized potatoes by HPLC with ICP-MS and electrospray ionization MS/MS	Infante, H. G.; Borrego, A. A.; Peachey, E.; Hearn, R.; O'Connor, G.; Barrera, T G.; Ariza, J. L.	<i>J. Agric. Food Chem.</i> 57(1), 38-45.	2009 Jan
Pentacyclic triterpene distribution in various plants – rich sources for a new group of multi-potent plant extracts	Jäger, S.; Trojan, H.; Kopp, T.; Laszczyk, M. N.; Scheffler, A.	<i>Molecules.</i> 14 (6), 2016-31.	2009 Jun
Comprehensive multiresidue method for the simultaneous determination of 74 pesticides and metabolites in traditional Chinese herbal medicines by accelerated solvent extraction with high-performance liquid chromatography/tandem mass spectrometry	Jia, Z.; Mao, X.; Chen, K.; Wang, K.; Ji S.	<i>J. AOAC Int.</i> ; 93(5), 1570-88.	2010 Sep-Oct
Gas chromatography-mass spectrometry (GC-MS) method for the determination of 16 European priority polycyclic aromatic hydrocarbons in smoked meat products and edible oils	Jira, W.; Ziegenhals, K.; Speer, K.	<i>Food Addit. Contam. Part A Chem. Anal. Control Expo. Risk Assess.</i> 25 (6), 704-13.	2008 Jun
Assessing pressurized liquid extraction for the high-throughput extraction of marine-sponge-derived natural products	Johnson, T. A.; Morgan, M. V.; Aratow, N. A.; Estee, S. A.; Sashidhara, K. V.; Loveridge, S. T.; Segraves, N L.; Crews, P.	<i>J. Nat. Prod.</i> 73 (3), 359-64.	2010 Mar
Lipophilic stinging nettle extracts possess potent anti-inflammatory activity, are not cytotoxic and may be superior to traditional tinctures for treating inflammatory disorders	Johnson, T. A.; Sohn, J.; Inman, W. D.; Bjeldanes, L. F.; Rayburn, K.	<i>Phytomedicine</i> 20(2), 143-7.	2013 Jan
Effects of solvent and temperature on pressurized liquid extraction of anthocyanins and total phenolics from dried red grape skin	Ju Z. Y.; Howard, L. R.	<i>J. Agric. Food Chem.</i> 51 (18), 5207-13.	2003 Aug

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



References

Peer Reviewed Journals: Sample Preparation Methods

Title	Authors	Publication	Publication Date
Accelerated solvent extraction of ochratoxin A from rice samples	Juan, C.; González, L.; Soriano, J. M.; Moltó, J. C.; Mañes, J.	<i>J. Agric. Food Chem.</i> 53 (24), 9348–9351	2005
Accelerated solvent extraction of paclitaxel and related compounds from the bark of <i>Taxus cuspidate</i>	Kawamura, F.; Kikuchi, Y.; Ohira, T.; Yatagai, M.	<i>J. Nat. Prod.</i> 62 (2), 244–7.	1999 Feb
Determination of polybromodiphenyl ethers (PBDEs) in milk cream by gas chromatography-mass spectrometry	Kinani, S.; Bouchonnet, S.; Abjean, J.; Campargue, C.	<i>Food Addit. Contam. Part A Chem. Anal. Control Expo. Risk Assess.</i> 25 (8), 1007–14	2008 Aug
Determination of isoflavones in soy bits by fast column high-performance liquid chromatography coupled with UV-visible diode-array detection	Klejduš, B.; Mikelová, R.; Petřlová, J.; Potešil, D.; Adam, V.; Stiborová, J.; Hodek, P.; Vacek, J.; Kizek, R.; Kubán, V.	<i>J. Chromatogr., A.</i> 1084 (1–2), 19, 71–79	2005 Aug
Accelerated solvent extraction of lignin from <i>Aleurites moluccana</i> (candlenut) nutshells	Klein, A. P.; Beach, E. S.; Emerson, J. W.; Zimmerman, J. B.	<i>J. Agric. Food Chem.</i> 58 (18), 10045–8	2010 Sep
Application of TLC method with video scanning in estimation of daily dietary intake of specific flavonoids – preliminary studies	Koch, W.; Kukuła-Koch, W.; Marzec, Z.; Marc, D.	<i>Acta Pol. Pharm.</i> 70 (4), 611–20	2013 Jul-Aug
Evaluation of a fibrous cellulose drying agent in supercritical fluid extraction and pressurized liquid extraction of diverse pesticides	Lehotay, S. J.; Lee, C. H.	<i>J. Chromatogr., A.</i> 785 (1–2), 313–27	1997 Oct
Application of accelerated solvent extraction to the investigation of saikosaponins from the roots of <i>Bupleurum falcatum</i>	Li, W.; Liu, Z.; Wang, Z.; Chen, L.; Sun, Y.; Hou, J.; Zheng, Y.	<i>J. Sep. Sci.</i> 33 (12), 1870–6	2010 Jun
Applicability of accelerated solvent extraction for synthetic colorants analysis in meat products with ultrahigh performance liquid chromatography-photodiode array detection	Liao, Q. G.; Li, W. H.; Luo, L. G.	<i>Anal. Chim. Acta.</i> 716, 128–32	2012 Feb
Extraction, isolation, and purification of analytes from samples of marine origin – a multivariate task	Liguori, L.; Bjørsvik, H. R.	<i>J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.</i> 910, 46–53	2012 Dec
Investigation on levels of polybrominated diphenyl ethers in retail fish and egg products in Shenzhen	Liu, B.; Zhang, L. S.; Zhang, J. Q.; Jiang, Y. S.; Zhou, J.; Huang, H. Y.	<i>Zhonghua Yu Fang Yi Xue Za Zhi.</i> 45 (12), 1068–72	2011 Dec
Characterization of secondary volatile profiles in <i>Nigella sativa</i> seeds from two different origins using accelerated solvent extraction and gas chromatography-mass spectrometry	Liu, X.; Abd El-Aty, A. M.; Cho, S. K.; Yang, A.; Park, J. H.; Shim, J. H.	<i>Biomed. Chromatogr.</i> 26 (10), 1157–62	2012 Oct
Accelerated solvent extraction of monacolin K from red yeast rice and purification by high-speed counter-current chromatography	Liu, Y.; Guo, X.; Duan, W.; Wang, X.; Du, J.	<i>J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.</i> 878 (28), 2881–5	2010 Oct
Multiresidue determination of organophosphorus pesticides in ginkgo leaves by accelerated solvent extraction and gas chromatography with flame photometric detection	Lu, Y.; Yi, X.	<i>J. AOAC Int.</i> 88 (3), 729–735	2005

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



References

Peer Reviewed Journals: Sample Preparation Methods

Title	Authors	Publication	Publication Date
Influence of sample preparation on assay of phenolic acids from eggplant	Luthria, DL.; Mukhopadhyay, S.	<i>J. Agric. Food Chem.</i> 54 (1), 41–47	2006
Pressurised solvent extraction for organotin speciation in vegetable matrices	Marcic, C.; Lespe, S G.; Potin-Gautier, M.	<i>Anal. Bioanal. Chem.</i> 382 (7), 1574–83	2005 Aug
Comparison of different methods for the determination of the oil content in oilseeds	Matthäus, B.; Brühl, L.	<i>J. AOCS</i> 78 95–102.	2001 Jan
A comparison of automated and traditional methods for the extraction of arsenicals from fish	McKiernan, J. W.; Creed, J. T.; Brockhoff, C. A.; Caruso, J. A.; Lorenzana, R. M.	<i>J. Anal. At. Spectrom.</i> 14, 607–613	1999
Subcritical solvent extraction of anthocyanins from dried red grape pomace	Monrad, J. K.; Howard, L. R.; King, J.; Srinivas, K.; Mauromoustakos, A.	<i>J. Agric. Food Chem.</i> 58 (5), 2862–8	2010 Mar
Subcritical solvent extraction of procyanidins from dried red grape pomace	Monrad, J. K.; Howard, L. R.; King, J. W.; Srinivas, K.; Mauromoustakos, A.	<i>J. Agric. Food Chem.</i> 58 (7), 4014–21	2010 Apr
Pressurized liquid extraction of polar and nonpolar lipids in corn and oats with hexane, methylene chloride, isopropanol, and ethanol	Moreau, R. A.; Powell, M. J.; Singh, V.	<i>J. Oil Fat Industr.</i> 80 (11), 1063–1067	2003 Jan
Accelerated solvent extraction for natural products isolation	Mottaleb, M. A.; Sarker, S. D.	<i>Methods Mol. Biol.</i> 864, 75–87	2012
Optimization of extraction process for phenolic acids from black cohosh (<i>Cimicifuga racemosa</i>) by pressurized liquid extraction	Mukhopadhyay, S.; Luthria, D. L.; Robbins, R. J.	<i>J. Sci. Food Agric.</i> 86 (1), 156–162, 15	2006 Jan
Anxiolytic activity of a supercritical carbon dioxide extract of <i>Souroubea sympetala</i> (Marcgraviaceae)	Mullally, M.; Kramp, K.; Cayer, C.; Saleem, A.; Ahmed, F; McRae, C.; Baker, J.; Goulah, A.; Otorola, M.; Sanchez, P.; Garcia, M.; Poveda, L.; Merali, Z.; Durst, T.; Trudeau, V. L.; Arnason, J. T.	<i>Phytother. Res.</i> 25 (2), 264–70	2011 Feb
On-line clean-up of pressurized liquid extracts for the determination of polychlorinated biphenyls in feedingstuffs and food matrices using gas chromatography–mass spectrometry	Müller, A.; Björklund, E.; von Holst, C.	<i>J. Chromatogr., A.</i> 925 (1–2), 197–205	2001 Aug
Analysis of multiple herbicides in soybeans using pressurized liquid extraction and capillary electrophoresis	Nemoto, S.; Lehotay, S. J.	<i>J. Agric. Food Chem.</i> ; 46 (6), 2190–2199	1998
Comparison of sample preparation methods, validation of an UPLC-MS/MS procedure for the quantification of tetrodotoxin present in marine gastropods and analysis of pufferfish	Nzoughet, J. K.; Campbell, K.; Barnes, P.; Cooper, K. M.; Chevallier, O. P; Elliott, C. T.	<i>Food Chem.</i> 15; 136 (3-4), 1584–9	2013 Feb
Multiresidue analysis of pesticides in vegetables and fruits using two-layered column with graphitized carbon and water absorbent polymer	Obana, H.; Akutsu, K.; Okihashi, M.; Hori, S.	<i>The Analyst</i> 123, 711–714	1998
Analysis of 2-alkylcyclobutanones with accelerated solvent extraction to detect irradiated meat and fish	Obana, H.; Furuta, M.; Tanaka, Y.	<i>J. Agric. Food Chem.</i> 53 (17), 6603–8	2005 Aug

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



References

Peer Reviewed Journals: Sample Preparation Methods

Title	Authors	Publication	Publication Date
Determination of organophosphorus pesticides in foods using an accelerated solvent extraction system	Obana, H.; Kikuchi, K.; Okihashi, M.; Hori, S.	<i>Analyst</i> 122 (3), 217–20	1997 Mar
Pressurized hot water extraction of berberine, baicalin and glycyrrhizin in medicinal plants	Ong, E. S.; Shea Mei, L.	<i>Anal. Chim. Acta.</i> 482 (1), 81–89	2003 Apr
Pressurized liquid extraction of berberine and aristolochic acids in medicinal plants	Ong E. S.; Woo S. O.; Yong, Y. K.	<i>J. Chromatogr., A.</i> 904 (1), 57–6422	2000 Dec
Rapid determination of pesticide multiresidues in vegetables and fruits by accelerated solvent extraction coupled with online gel permeation chromatography-gas chromatography-mass spectrometry	Ouyang, Y.; Tang, H.; Wu, Y.; Li, G.	<i>Se Pu.</i> 30(7), 654–9	2012 Jul
Determination of zearalenone from wheat and corn by pressurized liquid extraction and liquid chromatography-electrospray mass spectrometry	Pallaroni, L.; von Holst, C.	<i>J. Chromatogr., A.</i> 993, 39–45	2003
Development of an extraction method for the determination of zearalenone in corn using less organic solvents	Pallaroni, L.; von Holst, C.	<i>J. Chromatogr., A.</i> 5 1055 (1-2), 247–9	2004 Nov
Stability of phenolic compounds during extraction with superheated solvents	Palma, M.; Piñeiro, Z.; Barroso, C. G.	<i>J. Chromatogr., A.</i> 6 921 (2), 169–74	2001 Jul
Extraction and analysis of trace amounts of cyclonite (RDX) and its nitroso-metabolites in animal liver tissue using gas chromatography with electron capture detection (GC-ECD)	Pan, X.; Zhang, B.; Cobb, G. P.	<i>Talanta</i> 67 (4), 816–23	2005 Oct
Simultaneous determination of 405 pesticide residues in grain by accelerated solvent extraction then gas chromatography-mass spectrometry or liquid chromatography-tandem mass spectrometry	Pang, G.; Liu, Y.; Fan, C.; Zhang, J.; Cao, Y.; Li, X.; Li, Z.; Wu, Y.; Guo, T.	<i>Anal. Bioanal. Chem.</i> 384, 1366–1408	2006 Mar
Automated sample preparation by pressurized liquid extraction-solid-phase extraction for the liquid chromatographic-mass spectrometric investigation of polyphenols in the brewing process	Papagiannopoulos, M.; Mellenthin, A.	<i>J. Chromatogr., A.</i> 8 976 (1-2), 345–8	2002 Nov
Online coupling of pressurized liquid extraction, solid-phase extraction and high-performance liquid chromatography for automated analysis of proanthocyanidins in malt	Papagiannopoulos, M.; Zimmermann, B.; Mellenthin, A.; Krappe, M.; Maio, G.; Galensa, R.	<i>J. Chromatogr., A.</i> 7 958 (1-2), 9–16	2002 Jun
Simultaneous determination of 13 quinolones from feeds using accelerated solvent extraction and liquid chromatography	Pecorelli, I.; Galarini, R.; Bibi, R.; Floridi, A. I.; Casciarri, E.; Floridi, A.	<i>Anal. Chim. Acta.</i> 483 (1-2), 81–89	2003 April
Comparison of soxhlet, ultrasound-assisted and pressurized liquid extraction of terpenes, fatty acids and Vitamin E from <i>Piper gaudichaudianum</i> Kunth	Péres, V. F.; Saffi, J.; Melecchi, M. I.; Abad, F. C.; de Assis Jacques, R.; Martinez, M. M.; Oliveira, E. C.; Caramão, E. B.	<i>J. Chromatogr., A.</i> 1105 (1-2), 115–8	2006 Feb
Pressurised fluid extraction (PFE) as an alternative general method for the determination of pesticide residues in rape seed	Pihlström, T.; Isaac, G.; Waldebäck, M.; Osterdahl, B. G.; Markides, K. E.	<i>Analyst</i> 127 (4), 554–9	2002 Apr
Determination of catechins by means of extraction with pressurized liquids	Piñeiro, Z.; Palma, M.; Barroso C. G.	<i>J. Chromatogr., A.</i> 13 1026 (1-2), 19–23.	2004 Feb



Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)

Peer Reviewed Journals: Sample Preparation Methods

Title	Authors	Publication	Publication Date
An improved clean-up strategy for simultaneous analysis of polychlorinated dibenzo-p-dioxins (PCDD), polychlorinated dibenzofurans (PCDF), and polychlorinated biphenyls (PCB) in fatty food samples	Pirard, C.; Focant, J. F.; De, P. E.	<i>Anal. Bioanal. Chem.</i> 372 (2), 373–81.	2002 Jan
Extraction of polar and hydrophobic pollutants using accelerated solvent extraction (ASE)	Pörschmann, J., Plugge, J.	<i>Fresen. J. Anal. Chem.</i> 364 (7), 643–645	1999
Quantification of the total amount of artemisinin in leaf samples by thin layer chromatography	Quennoz, M.; Bastian, C.; Simonnet, X.; Grogg, A. F.	<i>Chimia (Aarau)</i> 64 (10), 755–7.	2010
Determination of fat in dairy products using pressurized solvent extraction	Richardson, R. K.	<i>J. AOAC Int.</i> 84 (5), 1522–1533	2001
Influence of altitudinal variation on the content of phenolic compounds in wild populations of <i>Calluna vulgaris</i>, <i>Sambucus nigra</i>, and <i>Vaccinium myrtillus</i>	Rieger, G.; Müller, M.; Guttenberger, H.; Bucar, F.	<i>J. Agric. Food Chem.</i> 56 (19), 9080–6.	2008 Oct
Pressurized liquid extraction of isoflavones from soybeans	Rostagno, M. A.; Palma, M.; Barroso, C. G.	<i>Anal. Chim. Acta.</i> 522 (2), 169–177.	2004 Sep
A multi-residue method for the analysis of organophosphorus residues in cooked and polished rice using accelerated solvent extraction and dispersive-solid phase extraction (D-SPE) technique and uncertainty measurement	Sanyal, D.; Rani, A.; Alam, S.	<i>J. Environ. Sci. Health, B</i> 44 (7), 706–16.	2009 Sep
Accelerated solvent extraction of lipids for determining the fatty acid composition of biological material	Schäfer, K.	<i>Anal. Chim. Acta.</i> 358 (1), 69–77	1998 Jan
HPLC analysis of kaempferol and quercetin derivatives isolated by different extraction techniques from plant matrix	Skalicka-Wozniak, K.; Szypowski, J.; Glowniak, K.	<i>J. AOAC Int.</i> 94 (1), 17–21.	Jan-Feb 2011
Statistical evaluation of fatty acid profile and cholesterol content in fish (common carp) lipids obtained by different sample preparation procedures	Spiric, A.; Trbovic, D.; Vranic, D.; Djinic, J.; Petronijevic, R.; Matekalo-Sverak, V.	<i>Anal. Chim. Acta.</i> 672 (1-2), 66–71.	2010 Jul
Application of accelerated solvent extraction in the analysis of organic contaminants, bioactive and nutritional compounds in food and feed	Sun, H.; Ge, X.; Lv, Y.; Wang, A.	<i>J. Chromatogr., A.</i> 1237, 1–23.	2012 May
Development of an accelerated solvent extraction, ultrasonic derivatisation LC-MS/MS method for the determination of the marker residues of nitrofurans in freshwater fish	Tao, Y.; Chen, D.; Wei, H.; Yuanhu, P.; Liu, Z.; Huang, L.; Wang, Y.; Xie, S.; Yuan, Z.	<i>Food Addit. Contam. Part A Chem. Anal. Control Expo. Risk Assess.</i> 29 (5), 736–45.	2012
Simultaneous determination of lincomycin and spectinomycin residues in animal tissues by gas chromatography-nitrogen phosphorus detection and gas chromatography-mass spectrometry with accelerated solvent extraction	Tao, Y.; Chen, D.; Yu, G.; Yu, H.; Pan, Y.; Wang, Y.; Huang, L.; Yuan, Z.	<i>Food Addit. Contam. Part A Chem. Anal. Control Expo. Risk Assess.</i> 28 (2), 145–54.	2011 Feb
Determination of 17 macrolide antibiotics and avermectins residues in meat with accelerated solvent extraction by liquid chromatography-tandem mass spectrometry	Tao, Y.; Yu, G.; Chen, D.; Pan, Y.; Liu, Z.; Wei, H.; Peng, D.; Huang, L.; Wang, Y.; Yuan, Z.	<i>J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.</i> 897, 64–71.	2012 May
Determination of seven toxaphene congeners in ginseng and milkvetch root by gas chromatography tandem mass spectrometry	Tian, S.; Mao, X.; Miao, S.; Jia, Z.; Wang, K.; Ji, S.	<i>Se Pu.</i> 30 (1), 14–20.	2012 Jan

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



References

Peer Reviewed Journals: Sample Preparation Methods

Title	Authors	Publication	Publication Date
A consecutive preparation method based upon accelerated solvent extraction and high-speed counter-current chromatography for isolation of aesculin from <i>Cortex fraxinus</i>	Tong, X.; Zhou, T, Xiao, X.; Li, G.	<i>J. Sep. Sci.</i> 35 (24), 3609–14	2012 Dec
Characterization of anthocyanins and anthocyanidins in purple-fleshed sweetpotatoes by HPLC-DAD/ESI-MS/MS	Truong, V. D.; Deighton, N.; Thompson, R. T.; McFeeters, R. F.; Dean, L. O.; Pecota, K. V.; Yencho, G. C.	<i>J. Agric. Food Chem.</i> 58 (1), 404–10	2010 Jan
Fat extraction from acid- and base-hydrolyzed food samples using accelerated solvent extraction	Ullah, S. M.; Murphy, B.; Dorich, B.; Richter, B.; Srinivasan, K.	<i>J. Agric. Food Chem.</i> 59 (6), 2169–74.	2011 Mar
Analysis of zearalenone in cereal and swine feed samples using an automated flow-through immunosensor	Urraca, J. L.; Benito-Peña, E.; Pérez-Conde, C.; Moreno-Bondi, M. C.; Pestka, J. J.	<i>J. Agric. Food Chem.</i> 53 (9), 3338–3344	2005
Accelerated solvent extraction and gas chromatography/mass spectrometry for determination of polycyclic aromatic hydrocarbons in smoked food samples	Wang, G.; Lee, A. S.; Lewis, M.; Kamath, B.; Archer, R. K.	<i>J. Agric. Food Chem.</i> 47 (3), 1062–6.	1999 Mar
Subcritical water extraction of alkaloids in <i>Sophora flavescens</i> Ait. and determination by capillary electrophoresis with field-amplified sample stacking	Wang, H.; Lu, Y.; Chen, J.; Li, J.; Liu, S.	<i>J. Pharm. Biomed. Anal.</i> 58, 146–51.	2012 Jan
Evaluation of Soxhlet extraction, accelerated solvent extraction and microwave-assisted extraction for the determination of polychlorinated biphenyls and polybrominated diphenyl ethers in soil and fish samples	Wang, P.; Zhang, Q.; Wang, Y.; Wang, T.; Li X.; Ding, L.; Jiang, G.	<i>Anal. Chim. Acta.</i> 663 (1), 43–8.	2010 Mar
Determination of ten pesticides of pyrazoles and pyrroles in tea by accelerated solvent extraction coupled with gas chromatography-tandem mass spectrometry	Xu, D.; Lu, S.; Chen, D.; Lan, J.; Zhang, Z.; Yang, F.; Zhou, Y.	<i>Se Pu.</i> ; 31 (3), 218–22.	2013 Mar
Online cleanup of accelerated solvent extractions for determination of adenosine 5'-triphosphate (ATP), adenosine 5'-diphosphate (ADP), and adenosine 5'-monophosphate (AMP) in royal jelly using high-performance liquid chromatography	Xue, X.; Wang, F.; Zhou, J.; Chen, F.; Li, Y.; Zhao, J.	<i>J. Agric. Food Chem.</i> 57 (11), 4500–5.	2009 Jun
Identification and quantitation of eleven sesquiterpenes in three species of <i>Curcuma</i> rhizomes by pressurized liquid extraction and gas chromatography–mass spectrometry	Yang, F. Q.; Li, S.; Chen, Y.; Lao, S. C.; Wang, Y.T.; Dong, T. T. X.; Tsim, K. W. K.	<i>J. Pharm. Biomed. Anal.</i> 39 (3/4), 552–558	2005 Sep
Dispersive solid-phase extraction cleanup combined with accelerated solvent extraction for the determination of carbamate pesticide residues in <i>Radix glycyrrhizae</i> samples by UPLC-MS-MS	Yang, R. Z.; Wang, J. H.; Wang, M. L.; Zhang, R.; Lu, X. Y.; Liu, W. H.	<i>J. Chromatogr. Sci.</i> 49 (9), 702–8.	2011 Oct
Simultaneous determination of amitraz and its metabolite residue in food animal tissues by gas chromatography-electron capture detector and gas chromatography-mass spectrometry with accelerated solvent extraction	Yu, H.; Tao, Y.; Le, T.; Chen, D.; Ishsan, A.; Liu, Y.; Wang, Y.; Yuan, Z.	<i>J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.</i> 878 (21), 1746–52.	2010 Jul
Simultaneous determination of fluoroquinolones in foods of animal origin by a high performance liquid chromatography and a liquid chromatography tandem mass spectrometry with accelerated solvent extraction	Yu, H.; Tao, Y.; Chen, D.; Pan, Y.; Liu, Z.; Wang, Y.; Huang, L.; Dai, M.; Peng, D.; Wang, X.; Yuan, Z.	<i>J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.</i> 885-886, 150–9.	2012 Feb

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



References

Peer Reviewed Journals: Sample Preparation Methods

Title	Authors	Publication	Publication Date
Determination of pentachlorophenol residue in meat and fish by gas chromatography-electron capture detection and gas chromatography-mass spectrometry with accelerated solvent extraction	Zhao, D.	<i>J. Chromatogr. Sci.</i>	2013 May
Response surface modeling and optimization of accelerated solvent extraction of four lignans from <i>fructus schisandrae</i>	Zhao, L. C.; He, Y. Deng.; X, Yang, G. L.; Li, W.; Liang, J.; Tang, Q. L.	<i>Molecules. 17 (4),</i> 3618–29	2012 Mar
Determination of acetanilide herbicides in cereal crops using accelerated solvent extraction, solid-phase extraction and gas chromatography-electron capture detector	Zhang, Y.; Yang, J.; Shi, R.; Su, Q.; Yao, L.; Li, P.	<i>J. Sep. Sci. 34 (14),</i> 1675–82	2011 Jul
Application of accelerated solvent extraction coupled with high-performance counter-current chromatography to extraction and online isolation of chemical constituents from <i>Hypericum perforatum L</i>	Zhang, Y.; Liu, C.; Yu, M.; Zhang, Z.; Qi, Y.; Wang, J.; Wu, G.; Li, S.; Yu, J.; Hu, Y.	<i>J. Chromatogr., A. 1218 (20),</i> 2827–34	2011 May
Analysis of volatile components in Qingshanlvshui tea using solid-phase microextraction/accelerated solvent extraction-gas chromatography-mass spectrometry	Zhan, J.; Lu, S.; Meng, Z.; Xiang, N.; Cao, Q.; Miao, M.	<i>Se Pu. 26 (3),</i> 301–5.	2008 May





Technical Collateral: Sample Preparation Methods

Product Number	Technique	Title
AN 326	HPLC-UV	Extraction of Drugs from Animal Feeds Using Accelerated Solvent Extraction (ASE)
AN 335	HPLC-UV	Accelerated Solvent Extraction (ASE) of Active Ingredients from Natural Products
AN 356	IC-conductivity	Determination of Perchlorate in Vegetation Samples Using Accelerated Solvent Extraction and Ion Chromatography
AN 357	HPLC	Extraction of Phenolic Acids from Plant Tissue Using Accelerated Solvent Extraction (ASE)
AN 363	HPLC	Extraction of Herbal Marker Compounds Using Accelerated Solvent Extraction Compared to Traditional Pharmacopoeia Protocols

Table of Contents

[Introduction](#)

[Analytical Technologies](#)

[Artificial Colorants](#)

[Blue Dyes](#)

[Caramel](#)

[General Dye Method](#)

[Global Method](#)

[Red Dyes](#)

[Sudan Dyes](#)

[Yellow Dyes](#)

[Natural Colorants](#)

[Turmeric](#)

[Flavorants](#)

[Amadori Compounds](#)

[Capsaicin](#)

[Quinine](#)

[Vanilla](#)

[Miscellaneous](#)

[Caffeine](#)

[Choline](#)

[Glucosamine](#)

[Iodide](#)

[Polyphosphates](#)

[Preservatives](#)

[Benzoate](#)

[Nitrate and Nitrite](#)

[Sulfites](#)

[Soft Drinks](#)

[References](#)



www.thermoscientific.com/liquidchromatography

©2015 Thermo Fisher Scientific Inc. All rights reserved. ISO is a trademark of the International Standards Organization. All other trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific products. It is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.



Thermo Fisher Scientific,
Sunnyvale, CA USA is
ISO 9001:2008 Certified.

Africa +43 1 333 50 34 0
Australia +61 3 9757 4300
Austria +43 810 282 206
Belgium +32 53 73 42 41
Canada +1 800 530 8447
China 800 810 5118 (free call domestic)
400 650 5118

Denmark +45 70 23 62 60
Europe-Other +43 1 333 50 34 0
Finland +358 9 3291 0200
France +33 1 60 92 48 00
Germany +49 6103 408 1014
India +91 22 6742 9494
Italy +39 02 950 591

Japan +81 45 453 9100
Korea +82 2 3420 8600
Latin America +1 561 688 8700
Middle East +43 1 333 50 34 0
Netherlands +31 76 579 55 55
New Zealand +64 9 980 6700
Norway +46 8 556 468 00

Russia/CIS +43 1 333 50 34 0
Singapore +65 6289 1190
Spain +34 914 845 965
Sweden +46 8 556 468 00
Switzerland +41 61 716 77 00
UK +44 1442 233555
USA +1 800 532 4752

Thermo
S C I E N T I F I C

A Thermo Fisher Scientific Brand