thermoscientific



Fast determination of inorganic cations and low mass amines in a spoiled cranberry juice sample using IC-MS

Authors

Terri Christison, Alexander N. Semyonov, Jeffrey S. Rohrer Thermo Scientific, Sunnyvale, CA

Keywords

HPIC, Integrion, Dionex IonPac CS12A-5µm column, ISQ EC mass spectrometer, single quadrupole mass spectrometer, Dionex CERS 500e Suppressor, suppressed conductivity detection

Introduction

Reliable analytical methods for food and beverage samples are required in order to report ingredients for food labeling requirements, to maintain product quality, and to ensure the absence of food contamination and spoilage. Food and beverage samples are almost invariably one of the most complex sample matrices and, therefore, are among the most challenging samples to analyze.

Ion chromatography (IC), using eluent generation and suppressed conductivity detection, provides chromatographic selectivity, low chemical noise, and high compatibility with a mass spectrometer (MS).^{1,2} Additionally, the analytes leave the IC system as ions, further increasing MS compatibility.

Mass spectrometry coupled to IC provides higher selectivity and often, better detection limits. In addition, MS is one of the most universal detectors, providing powerful screening, structural, and confirmatory information.² The heated electrospray ionization (HESI), with applied high temperature aiding voltage, transforms the aqueous IC stream into a fine spray that can enter the MS detector.¹



In this application brief, separations of four inorganic cations and diethylamine in a spoiled cranberry juice sample were achieved within six minutes. The analytes were detected serially, by suppressed conductivity and single quadrupole mass spectrometry in full scan (FS) and SIM (select ion monitoring) modes without the aid of a desolvation agent during separation or ionization.

This application brief is an update to Application Note 269.³ Here, the application is performed using a Thermo Scientific[™] Dionex[™] Integrion[™] HPIC[™] system coupled to a Thermo Scientific[™] ISQ[™] EC single quadrupole mass spectrometer.

Experimental

Ion Chromatography

- Dionex Integrion HPIC system, RFIC[™] model with a second six-port high-pressure divert valve and Conductivity Detector
- Thermo Scientific[™] Dionex[™] AS-AP autosampler
- Thermo Scientific[™] Dionex[™] AXP-MS auxiliary pump to supply water for the suppressor

Mass Spectrometry

- ISQ EC single quadrupole mass spectrometer
- Thermo Scientific syringe pump for method optimization
- HESI II probe

Software

Thermo Scientific[™] Chromeleon[™] Chromatography Data System (CDS) software, 7.2 SR 6

Methods

Columns:	Thermo Scientific™ Dionex™ IonPac™ CG12A-5µm, CS12A-5µm, 3 mm i.d.		
Eluent:	33 mM N	lethanesulfoni	c acid (MSA)
Eluent Source:	MSA cart Dionex™ (ridge, Thermo	trap column, high
Flow Rate:	0.5 mL/min		
Injection Volume:	100 µL		
Detection 1:	CERS 50	0.7 mL/min by	r, external water
Typical Conductance Background:	< 1 µS-m	in	
MS Detection:*	+ESI, +3000 V, Full Scan, 18-250 <i>m/z</i> and SIM, HESI II		
Temperatures:	Vaporizer: 250 °C; Ion Transfer: 300 °C		
Flow (N ₂):	Sheath: 60 psi, Aux: 26 psi; Sweep: 0.5 psi		
Desolvation agent:	None		
SIM mode:		lon <i>m/z</i>	CID (V)
Sodium as Na•2H ₂ C)	59	10
Ammonium as $NH_4 \bullet H_2O$		36	2
Potassium		39	45
Magnesium as Mg ₂ •H ₂ O		66	5
Calcium		40	45
Dimethylamine		46	10
Ethylamine		46	45
Monoethanolamine		62	15
Diethylamine		74	15
Triethanolamine		150	25

*Note: The optimum HESI ion source settings and responses may vary between instruments.

Results

Figure 1 shows the IC-MS results of inorganic cations and amines in a diluted spoiled cranberry juice sample with no visible mold. The ions were detected in SIM mode from m/z 39 to 150 as native ions or water-adducts. The IC and SIM peaks are symmetrical and have strong responses, E2 to E5 ion counts (in SIM mode). Bare, unsolvated calcium m/z 40 had the lowest response at 3e2 (300) counts. Notice that Peak 4, diethylamine at m/z 74, was detected only by the MS detector, thereby demonstrating the sensitivity and selective detection possible by MS. This application brief demonstrates a fast and accurate method using ion chromatography separations and mass spectrometry detection to identify spoilage in juice samples. More information can be found in Thermo Scientific[™] AppLabs Library of Analytical Applications.⁴

References

- Wang, J. and Schnute, W.C. Optimizing mass spectrometric detection for ion chromatographic analysis. I. Common anions and selected organic acids. *Rapid Commun. Mass Spectrom.*, 2009, 23 (21), 3439–3447. DOI:10.1002/rcm.4263
- Thermo Scientific White Paper 70424: Specific and Selective Detection for Food and Beverage Analysis by Ion Chromatography-Mass Spectrometric Detection. 2016. <u>http://tools.thermofisher.com/content/sfs/brochures/</u> <u>WP70424-specific-and-selective-detection-for-food-and-beverage-analysis-by-ionchromatography-mass-spectrometric-detection.pdf (accessed Jun. 9, 2017).</u>
- Thermo Scientific Application Note 269: Identification and Quantification at ppb Levels of Common Cations and Amines by IC-MS. <u>https://appslab.thermofisher.com/</u> <u>Search?SearchText=an269</u>
- 4. Thermo Scientific AppsLab Library of Analytical Applications. <u>https://appslab.</u> thermofisher.com



urce:	Dionex IonPac CG12A, CS12A-5 µm, 3 mm 33 mM Methanesulfonic Acid (MSA) Dionex EGC 500 MSA cartridge, Dionex CR-CTC 600 trap column 0.5 mL/min
ime:	100 μL
perature:	30 °C
1:	Suppressed conductivity, Dionex CERS 500e suppressor, 49 mA, 20 °C, external water mode, 0.7 mL/min by Dionex AXP-MS pump
2:	ISQ EC, +ESI, +3000 V source, HESI II
de:	Full scan: 18-250 m/z, SIM
solvent:	none
mp.:	Vaporizer 250 °C , Ion Transfer 300 °C
ow (psi):	Sheath 60, Aux 26, Sweep 0.5
rep.:	1000-fold dilution with DI

Figure 1. Determination of co-eluting amines in a two-year-old cranberry juice sample.

Find out more at www.thermofisher.com/ISQEC

©2017 Thermo Fisher Scientific Inc. All rights reserved. All 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 manners 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 representatives for details. **AB72404EN 0617S**

