## thermo scientific



**APPLICATION BRIEF 72480** 

# Detection of common inorganic anions using a compact ion chromatography system coupled with a single quadrupole mass spectrometer

#### Authors

Beibei Huang,<sup>1</sup> Joshua T. Maze,<sup>2</sup> Kyle B. Renfrew,<sup>2</sup> and Jeffrey Rohrer<sup>1</sup>

<sup>1</sup>Thermo Fisher Scientific, Sunnyvale, CA, USA; <sup>2</sup>ThermoFisher Scientific, Austin, TX, USA

#### Keywords

Integrion HPIC, IonPac AS20 column, ISQ EC mass spectrometer, drinking water, IC-MS, inorganic anions

#### Introduction

This application brief describes a method for detecting seven common anions using a Thermo Scientific<sup>™</sup> Dionex<sup>™</sup> Integrion<sup>™</sup> HPIC<sup>™</sup> system coupled to a Thermo Scientific<sup>™</sup> ISQ<sup>™</sup> EC single quadrupole mass spectrometer. The determination of common inorganic anions in drinking water is one of the most important ion chromatography (IC) applications worldwide. In the U.S., the Environmental Protection Agency (EPA) defines the Maximum Contaminant Levels (MCLs) of toxic contaminants, such as fluoride, nitrite, and nitrate in the National Primary Drinking Water Regulations (NPDWR).<sup>1</sup> EPA Methods 300.0<sup>2</sup> and 300.1<sup>3</sup> are IC methods approved by the U.S. EPA for analysis of common anions in the U.S.

As demonstrated in this brief, IC offers the analyst the advantage of selectivity for common low-mass anions. Mass detection adds confidence to inorganic anion identity and sometimes greater sensitivity than provided by conductivity detection. Samples were separated with a Thermo Scientific<sup>™</sup> Dionex<sup>™</sup> IonPac<sup>™</sup> AS20 column set using a Dionex Integrion HPIC system, and an ISQ EC Single Quadrupole Mass Spectrometer. Figure 1 shows chromatograms of a 0.2 mg/L fluoride, 0.3 mg/L chloride, 1 mg/L nitrite, bromide, and nitrate, and 1.5 mg/L phosphate and sulfate standard using suppressed conductivity detection followed by an MS detector in selected ion monitoring (SIM) mode. It is important to note that the SIM ion selected for fluoride is 39.1 *m/z*, which is associated with the fluoride adduct [F(HF)]<sup>-</sup>. The



deprotonated sulfate  $[HSO_4]^-$  and phosphate  $[H_2PO_4]^$ ions which have the same mass-to-charge ratio (*m/z*) at 97.1 are shown as two peaks in SIM channel 97.1, and can also be accurately quantified because they are chromatographically separated. The differentiation of sulfate from phosphate can be confirmed by the isotope peak 99.1 ( $[H^{34}SO_4]^-$ ) at ~5% relative peak intensity of peak 97.1. A smaller peak in SIM channel 97.1, which may be a single water complex of bromide,  $[Br(H_2O)]^-$ , appears at the same retention time as bromide.

Representative municipal drinking water and bottled water samples were analyzed by IC-MS (Figures 2 and 3). Using IC, the anionic components in water samples can be identified by comparing their retention times to those of the standards. To confirm these identifications, a mass selective detector can record SIM channels at the appropriate mass-to-charge ratios. In Figure 2, six anions (fluoride, chloride, nitrite, nitrate, sulfate, and phosphate) were identified in municipal drinking water on the basis of peak retention time and mass confirmation. Bromide was not detected by either detector. Although two peaks showed in the SIM channel for bromide, they did not match the retention time of bromide. Figure 3 shows the conductivity and SIM chromatogram of a bottled water sample. Neither nitrite nor bromide was detected in this sample. With the use of very specific MS SIM detection, ppb-level analyte identification can be achieved with greater confidence on a wide range of samples, such as bottled drinking water.

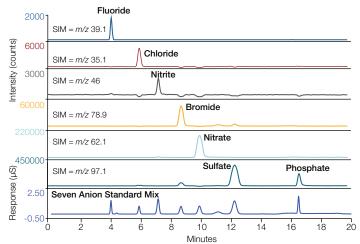
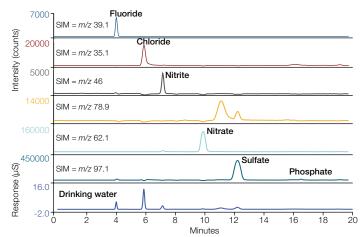
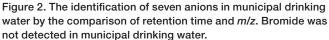


Figure 1. Conductivity and SIM chromatograms of a 0.2 mg/L fluoride, 0.3 mg/L chloride, 1 mg/L nitrite, bromide, nitrate and 1.5 mg/L phosphate and sulfate standard.





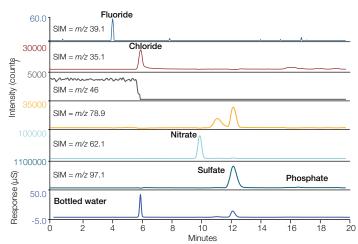


Figure 3. The identification of seven anions in bottled water by the comparison of retention time and m/z. Nitrite and bromide were not detected in this bottled water sample.

#### Method

Method						
Conditions						
IC System:	Dionex Integrion HPIC system					
MS Detector:	ISQ EC single quadrupole mass spectrometer					
Columns:	Dionex IonPac AG20 Guard, 2 × 50 mm Dionex IonPac AS20 Analytical, 2 × 250 mm Thereas Scientifica™ Diversu™ ECO 500 I/OLL Elwart Consultant Contriduc with Thereas Scientifica™					
Eluent Source:	Thermo Scientific <sup>™</sup> Dionex <sup>™</sup> EGC 500 KOH Eluent Generator Cartridge with Thermo Scientific <sup>™</sup> Dionex <sup>™</sup> CR-ATC 600 Continuously Rengenerated Anion Trap Column					
Gradient:	15 mM KOH (0–11 min), 45 mM KOH (11.10–15 min), 15 mM KOH (15–20 min)					
Flow Rate:	0.25 mL/min					
Injection Volume:	10 µL					
Temperature:	30 °C (column compartment), 20 °C (detector compartment)					
System Backpressure:	~2350 psi					
Detection:	Suppressed Conductivity, Dionex AERS 500e Anion Electrolytically Regenerated Suppressor (2 mm), AutoSuppression, 28 mA, external water mode via Thermo Scientific <sup>™</sup> AXP <sup>™</sup> Auxilliary Pump, external water flow rate (0.5 mL/min)					
Background Conductance:	~ 0.3 µS					
Run Time:	20 min					
Mass Spectrometric De	tection					
Ionization Interface:	Electrospray ionization (ESI), negative mode					
Gas Control:	Sheath gas pressure: 50 psi Aux gas pressure: 5 psi Sweep gas pressure: 0.0 psi					
Source Voltage:	-2500 V					
Vaporizer Temperature:	450 °C					
Ion Transfer Tube Temperature:	200 °C					
Advanced Scan Mode:	Scan Name Fluoride Chloride Nitrite Bromide Nitrate Sulfate, Phosphate	Mass list (amu) 39.1 35.1 46 78.9 62.1 97.1	Dwell or Scan Times (sec): 0.3 0.3 0.3 0.3 0.3 0.3 0.3	SIM Widths (amu): 0.1 0.1 0.1 0.1 0.1 0.1	lon Polarity: Negative Negative Negative Negative Negative Negative	Source CID Voltage: 20 20 20 20 20 20 20
Groups:	Chrom. Filter Peak Width (sec): 30					

#### References

- National Primary Drinking Water Regulations. Code of Federal Regulations, 40 CFR, Part 141; U.S. Environmental Protection Agency: Cincinnati, OH, 1998.
- 2.U.S. EPA Method 300.0., U.S. Environmental Protection Agency; Cincinnati, Ohio, 1993.
- 3.U.S. EPA Method 300.1, U.S. Environmental Protection Agency, Cincinnati, Ohio, 1997.

### 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. **AB72480-EN 0817S** 

