

**Environmental**

# Determination of cations and ammonium in environmental waters using a compact RFIC ion chromatography system

## Authors

Yukiko Kawahara, Jeffrey Rohrer  
Thermo Fisher Scientific,  
Sunnyvale, CA, USA

## Keywords

Dionex Inuvion IC, Dionex IonPac CS16-4 $\mu$ m column, drinking water, wastewater

## Introduction

The amount of alkali and alkaline earth cations and ammonium in drinking and environmental waters can vary significantly depending on the sample. A high-capacity cation exchange column installed on a reagent-free ion chromatography (RFIC™) system with suppressed conductivity detection provides the accuracy and sensitivity as well as a wide calibration range for determining these analytes.

In this application proof note, the performance of the easy-to-use, compact Thermo Scientific™ Dionex™ Inuvion™ ion chromatography system equipped with a Thermo Scientific™ Dionex™ IonPac™ CS16-4 $\mu$ m (4 × 250 mm) cation-exchange column was demonstrated and compared to previously published Thermo Scientific Application Update 204<sup>1</sup> to show the new instrument is suitable for this application.

## Method

### Reagents and standards

- Deionized (DI) water, Type I reagent grade, 18 M $\Omega$ ·cm resistance or better
- Calcium chloride dihydrate, Fisher Chemical (P/N C79)
- Magnesium chloride hexahydrate, Sigma-Aldrich (P/N M9272)
- Potassium chloride, Fisher Chemical (P/N P330)
- Ammonium Standard, 1,000  $\mu$ g/mL, Agilent Technologies (P/N ICC-101)
- Sodium chloride, Sigma-Aldrich (P/N S9888)
- Lithium chloride, Fisher Chemical (P/N L121)
- Thermo Scientific™ Dionex™ Combined Six Cation Standard-II (P/N 046070)

### Instrument method parameters

Instrument	Dionex Inuvion IC system (P/N 22185-60108), including column heater, pump degas module, and eluent generation
Autosampler	Thermo Scientific™ Dionex™ AS-DV autosampler (P/N 068907) with 5 mL Thermo Scientific™ Dionex™ PolyVials™ and filter caps (P/N 038141)
Columns	Dionex IonPac CS16-4 $\mu$ m, Analytical, 4 $\times$ 250 mm (P/N 088584) Dionex IonPac CG16-4 $\mu$ m, Guard, 4 $\times$ 50 mm (P/N 088585)
Eluent	30 mM MSA
Eluent source	Thermo Scientific™ Dionex™ EGC 500 MSA cartridge (P/N 075779), Thermo Scientific™ Dionex™ CR-CTC 600 (P/N 088663), Thermo Scientific™ Dionex™ RFIC™ eluent degasser module
Flow rate	0.90 mL/min
Column temp.	40 °C
Injection volume	10 $\mu$ L
Detection	Suppressed conductivity, Thermo Scientific™ Dionex™ CDRS 600 (4 mm) suppressor (P/N 088668CMD or 088668), 79 mA, recycle mode
System backpressure	~4,100 psi (100 psi = 689.5 kPa)
Background conductance	<0.5 $\mu$ S/cm
Noise	~0.2 nS/cm
Run time	22 min
Software	Thermo Scientific™ Chromeleon™ Chromatography Data System (CDS) software version 7.3.2

## Results and discussion

The analysis method in Application Update 204 was followed to demonstrate the performance of a Dionex Inuvion ion chromatography system.

Table 1 shows the calibration range and MDLs for each cation.

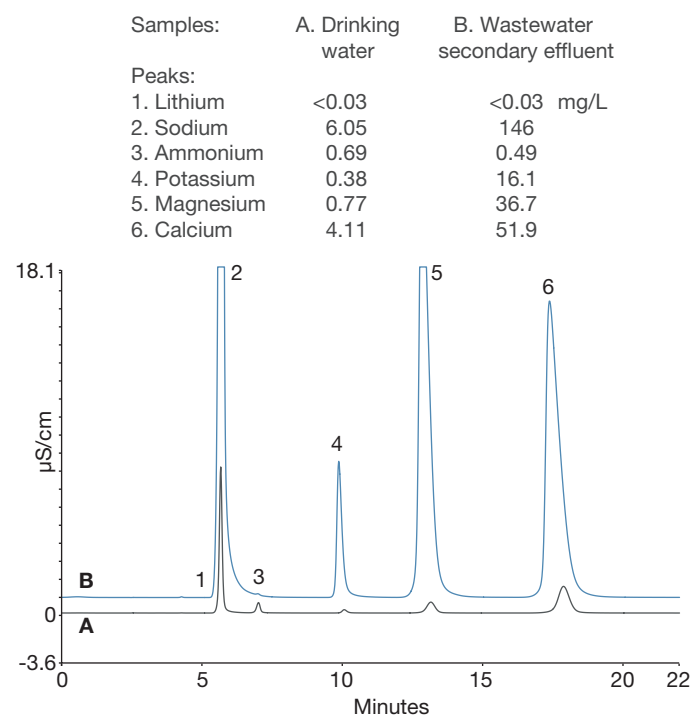
**Table 1. Working range and MDLs for cations and ammonium**

Cation	Range (mg/L)	Coefficient of determination ( $r^2$ )	MDL standard ( $\mu$ g/L)	Calculated MDL <sup>a</sup> ( $\mu$ g/L)
Li <sup>+</sup>	0.03–80	1.0000	1	0.14
Na <sup>+</sup>	0.1–1000	0.9999	4	0.59
NH <sub>4</sub> <sup>+</sup> <sup>b</sup>	0.02–40	0.9965	5	1.76
K <sup>+</sup>	0.03–80	0.9996	10	1.31
Mg <sup>2+</sup>	0.03–80	0.9999	5	0.11
Ca <sup>2+</sup>	0.03–80	0.9997	10	11.88

<sup>a</sup>MDL = (t)  $\times$  (S), where t = Student's t value for a 99% confidence level and a standard deviation estimate with n-1 degrees of freedom (t = 3.14 for seven replicates of the MDL standard), and S = standard deviation of the replicate analysis.

<sup>b</sup>Quadratic fit

Two samples, drinking water and wastewater secondary effluent, were tested for cations and ammonium. The wastewater sample was filtered and injected without dilution (Figure 1).



**Figure 1. Determination of inorganic cations and ammonium in environmental waters**

## Conclusions

The study successfully demonstrates the performance of a Dionex Inuvion IC system using a Dionex IonPac CS16-4 $\mu$ m column for determining cations and ammonium in environmental waters.

## Reference

1. Thermo Scientific Application Update 204: Analysis of Environmental Waters for Cations and Ammonium Using a Compact Ion Chromatography System. 2016 (accessed June 2023)

 Learn more at [thermofisher.com/inuvion](https://thermofisher.com/inuvion)