

Determination of Inorganic Anions in Municipal Water Using High-Pressure Modular Capillary Ion Chromatography

Carl Fisher, Fei Pang, Terri Christison, Linda Lopez
Thermo Fisher Scientific, Sunnyvale, CA, USA

Key words

Environmental water analysis, HPIC, Wastewater, ICS-5000+, 4 μm particle column, IonPac AS18-4 μm

Goal

Demonstrate fast separation of inorganic anions in municipal water samples by increasing flow rates using a high-pressure capillary IC system and 4 μm resin particle column.

Introduction

Ion chromatography (IC) has been widely adopted to determine inorganic anions in environmental waters, including surface, ground, drinking and wastewaters. In the U.S., the contamination level of wastewater is regulated by the Clean Water Act to minimize discharge of pollutants into the environment. The integrity and safety of drinking water is ensured through the Safe Drinking Water Act. Many other countries (e.g., Germany, France, Italy, China and Japan) have similar health and environmental standards and a considerable number of regulatory IC methods have been published. In addition, many standards organizations (including the International Organization for Standardization (ISO), American Society for Testing and Materials (ASTM), and American Water Works Association (AWWA) have validated IC methods for the analysis of inorganic anions in drinking water.^{1,2} For some anions, their concentration in drinking water is regulated due to their toxic effects. For example, high levels of fluoride cause skeletal and dental fluorosis³, while nitrite and nitrate can cause methemoglobinemia⁴, which can be fatal to infants. Other inorganic anions, such as chloride and sulfate, are considered secondary contaminants and can affect odor, color, and other aesthetic characteristics.



This Technical Note describes the determination of inorganic anions in municipal drinking water samples using the Thermo Scientific™ Dionex™ ICS-5000+ HPIC™ capillary system in combination with the Thermo Scientific™ Dionex™ IonPac™ AS18-4 μm column, which is the column of choice for compliance monitoring of water samples in accordance with U.S. EPA Methods 300.0 (A)⁵ and 300.1.⁶ By increasing the flow rate, multiple samples can be quickly analyzed, resulting in high throughput. Scaling down from standard bore (4 mm i.d.) to capillary format (0.4 mm i.d.) brings many benefits to IC users. The system can be always on and ready for analysis because of its low eluent consumption (15 mL of water a day at 0.01 mL/min flow rate). Additionally, the amount of waste generated is significantly reduced and the eluent generation cartridge (EGC) can last 18 months under standard continuous operation, which translates into reduced overall cost of ownership.

Equipment

- Dionex High-Pressure ICS-5000+ HPIC Reagent-Free™ (RFIC™) capillary IC system
 - Dionex ICS-5000+ DP Dual Pump module with high-pressure capillary pumps
 - Dionex ICS-5000+ EG Eluent Generator module
 - Dionex ICS-5000+ DC Detector/Chromatography module with a Thermo Scientific™ Dionex™ IC Cube™ and high-pressure degas cartridge
 - Thermo Scientific Dionex AS-AP Autosampler
- Thermo Scientific™ Dionex™ Chromeleon™ Chromatography Data System (CDS) software

Reagents and Standards

- 18.2 MΩ-cm degassed deionized water
- Anion Standard solution made from 1000 mg/L stocks

Samples

- Municipal drinking and wastewater diluted as indicated with deionized water

Conditions

Column:	Dionex IonPac AS18-4µm, 0.4 × 150 mm
Eluent Source:	Thermo Scientific Dionex EGC-KOH Cartridge (Capillary) with Thermo Scientific Dionex CR-ATC Continuously Regenerated Anion Trap Column
Eluent:	23 mM KOH
Flow Rate:	0.025 mL/min
Column Temp.:	30 °C
Inj. Volume:	0.4 µL
Detection:	Suppressed conductivity, Thermo Scientific™ Dionex™ ACES™ 300 Anion Capillary Electrolytic Suppressor, recycle mode
Background Conductance:	0.3–0.5 µS
Noise:	Less than 3 nS
System Backpressure:	~3500 psi

*Part numbers of the consumables used in this document are listed in Table 1.

Table 1. Consumables list.

Product Name	Description	Dionex Part Number
Dionex EGC-KOH (Capillary)	Eluent generator cartridge	072076
Dionex CR-ATC (Capillary)	Electrolytic trap column	072078
Dionex IonPac AS18-4µm	Separation column	082314
Dionex IonPac AG18-4µm	Guard column	076033
Thermo Scientific Dionex CRD Bypass cartridge	Bypass (needed for flow path)	072056
Dionex ACES 300	Suppressor	072052
Dionex HP fittings (blue)	Bolts / Ferrules	074449/ 074373
Thermo Scientific™ Dionex™ CRD 180 Carbonate Removal Cartridge	For 4 µm particle capillary columns	079960
EG Degas cartridge (Capillary)	High-pressure degas cartridge, up to 5000 psi	AAA-074459
Dionex AS-AP Autosampler vials	Package of 100, 10 mL, polystyrene vials, caps, blue septa	074228

Standard and Sample Preparation

The anion standards were prepared by diluting ~1000 mg/L stock solutions with 18 MΩ-cm degassed deionized water. The municipal wastewater samples were diluted 1000-fold and drinking water 5-fold with 18 MΩ-cm degassed deionized water and filtered with a 0.2 µm syringe filter prior to injection.

It is important to use 18 MΩ-cm resistivity, deionized water for standard, eluent, and autosampler flush solutions to avoid system contamination resulting in decreased sensitivity and poor calibration. Degassing the deionized water by vacuum filtration prior to use is a good practice. It prevents the formation of gas bubbles and increases the reproducibility of results.

Instrument Setup and Installation

To achieve the best chromatography with capillary IC, it is important to minimize void volumes in each connection by using precision cut tubing, high-pressure connectors and fittings (blue color), and seating the ferrule > 2 mm above the end of the tubing. Extra care should be used to prevent air in all consumables or tubing by observing steady liquid flow before installing the next device in line. A thorough discussion can be found in Technical Note 113, “Practical Guidance for Capillary IC”.⁷

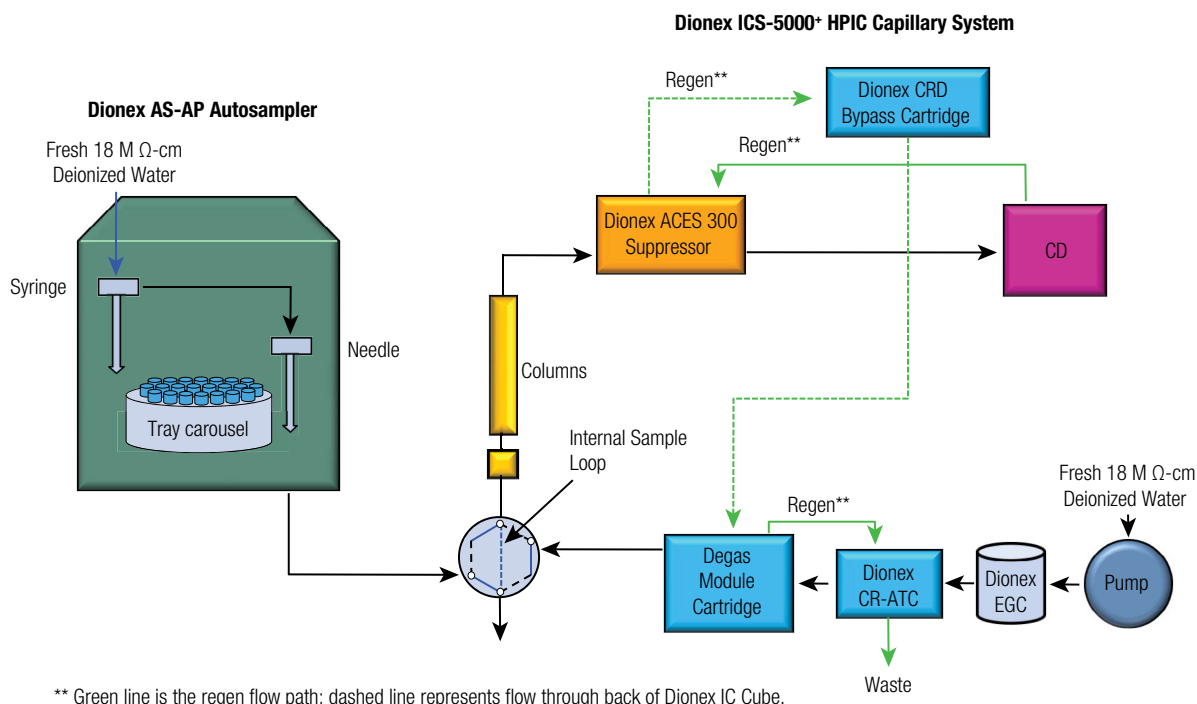


Figure 1. Flow diagram for the Dionex ICS-5000⁺ HPIC capillary system.

To set up this application, plumb the consumables and modules of the Dionex ICS-5000⁺ HPIC capillary system according to Figure 1. Install and hydrate the Dionex EGC-KOH cartridge, Dionex CR-ATC column, and suppressor following the instructions in the product manuals. The Dionex CRD Bypass Cartridge was installed for all experiments in this technical note. To achieve optimal column efficiencies and greatest removal of carbonate introduced by samples, install a Dionex CRD 180 Carbonate Removal Device. This device is optimized for 4 μm particle separation columns. Use high-pressure connectors and ferrules (blue) for all of the fittings from the pump to prior to the suppressor. Standard pressure fittings can be used for the suppressor and detector connections. Detailed instructions are described in Technical Note 131, the product manuals, and the instrument installation and operator's manuals.⁸⁻¹²

Results and Discussion

As illustrated in Figure 2, the time to separate nine inorganic anions was reduced from seven to less than three minutes by increasing the flow rate from 0.010 to 0.025 mL/min. This shortened run time enables increased throughput when analyzing large numbers of samples. At this higher flow rate, the backpressure of the system was about 3800 psi, which is above the 3000 psi limit of many IC systems, but well below the limit of the Dionex ICS-5000⁺ HPIC system (5000 psi). The use of higher flow rates, while maintaining resolution, is made possible by the 4 μm particle size of the Dionex IonPac AS18-4μm resin, which is about half the diameter of the previously available resin (7.5 μm). These 4 μm columns possess both higher capacity and efficiency. Despite the use of a faster flow rate, there is still only a modest amount of water consumed and waste generated (36 mL/day) keeping labor and waste disposal costs to a minimum.

Column:	Dionex IonPac AG18, Dionex IonPac AS18-4μm (0.4 × 150 mm)	Peaks:	1. Fluoride 0.1 mg/L
Eluent Source:	Dionex EGC-KOH Cartridge (Capillary)	2. Chlorite 1.0	3. Chloride 0.6
Eluent:	23 mM KOH	4. Nitrite 1.0	5. Carbonate –
Inj. Volume:	0.4 μL	6. Bromide 2.0	7. Sulfate 2.0
Col. Temp.:	30 °C	8. Nitrate 2.0	9. Chlorate 2.0
Flow Rate:	A. 0.010 mL/min B. 0.015 mL/min C. 0.025 mL/min		
Detection:	Suppressed conductivity, Dionex ACES 300, recycle mode		

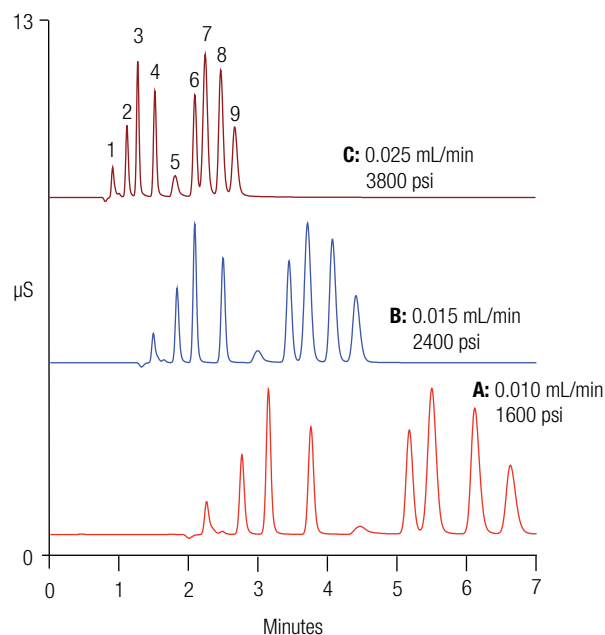


Figure 2. Rapid determination of inorganic anions using high-pressure capillary IC.

Figure 3 shows that inorganic anions in municipal drinking water samples were separated in less than three minutes when a flow rate of 0.025 mL/min was used. The concentration of fluoride that was measured ranged from 0.05 to over 1 mg/L, a 20-fold difference. Nitrate had the largest variance among the samples analyzed ranging from 0.16 for the lowest concentration detected to 17 mg/L for the highest, a difference of > 100-fold.

Column:	Dionex IonPac AG18-4 μ m, Dionex IonPac AS18-4 μ m, (0.4 \times 150 mm)				
Eluent Source:	Dionex EGC-KOH Cartridge (Capillary)				
Eluent:	23 mM KOH				
Flow Rate:	0.025 mL/min				
Inj. Volume:	0.4 μ L				
Column Temp.:	30 $^{\circ}$ C				
IC Cube Temp.:	15 $^{\circ}$ C				
Detection:	Suppressed conductivity, Dionex ACES 300, Anion Capillary Electrolytic Suppressor recycle mode				
Sample Prep:	A, C: Diluted 5-fold B, D: Undiluted				
Samples:	A: City 1 B: City 2 C: City 3 D: City 4				
Peaks:	A	B	C	D	
1. Fluoride	0.61	0.95	0.05	1.14	mg/L
2. Chloride	8.36	5.76	38.80	13.90	
3. Carbonate	—	—	—	—	
4. Sulfate	50.10	18.00	66.90	20.30	
5. Nitrate	7.09	0.16	17.00	—	

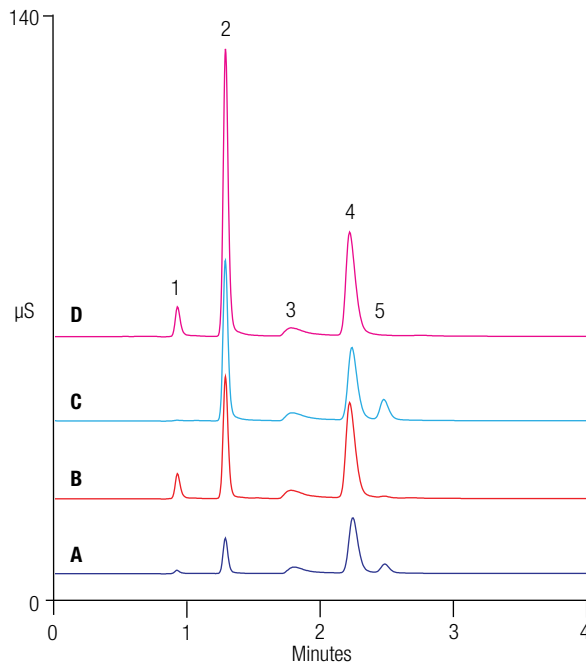


Figure 3. Determination of inorganic anions in municipal drinking water samples using high-pressure capillary IC.

In Figure 4, concentration of the inorganic anions in municipal wastewater were determined. Due to the high concentrations of ions present, these samples were diluted 1000-fold so that the quantities injected were well within the column capacity. Both nitrite and nitrate had a large range of concentrations in samples from different stages of the wastewater treatment process (25- and 18-fold, respectively).

Column:	Dionex IonPac AG18-4 μ m, Dionex IonPac AS18-4 μ m, (0.4 \times 150 mm)				
Eluent Source:	Dionex EGC-KOH Cartridge (Capillary)				
Eluent:	23 mM KOH				
Flow Rate:	0.025 mL/min				
Inj. Volume:	0.4 μ L				
Column Temp.:	30 $^{\circ}$ C				
IC Cube Temp.:	15 $^{\circ}$ C				
Detection:	Suppressed conductivity, Dionex ACES 300, Anion Capillary Electrolytic Suppressor recycle mode				
Sample Prep:	Diluted 1000-fold, filtered, 0.2 μ m				
Samples:	A: Influent B: Primary effluent C: Tricking effluent D: Final effluent				
Peaks:	A	B	C	D	
1. Chloride	76.5	146.0	154.0	130.0	mg/L
2. Nitrite	1.5	2.1	37.4	1.6	
3. Carbonate	—	—	—	—	
4. Sulfate	41.6	88.9	84.8	91.8	
5. Nitrate	28.8	7.2	31.7	128.0	

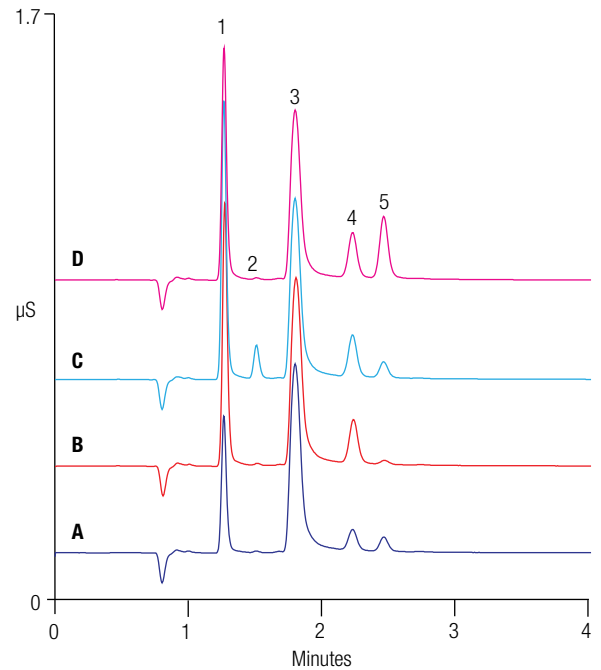


Figure 4. Determination of inorganic anions in municipal wastewater using high-pressure capillary IC.

Conclusion

This Technical Note demonstrates that the Dionex ICS-5000+ capillary HPLC system, when combined with the Dionex IonPac AS18-4 μ m column, provides an ideal solution for the analysis of inorganic anions in municipal water samples. The high-pressure capability of this system permits the use of higher flow rates resulting in short run times, while the 4 μ m resin particle column ensures good peak resolution.

For information on inorganic ion determinations in water using standard and microbore flow rates, see Application Notes 130 and 154.^{13,14}

References

- Greenberg, A. E.; Clesceri, L. S.; Eaton, A. D., Eds.; Standard Methods for the Examination of Water and Wastewater, 18th ed.; Am. Public Health Assoc.; Washington, DC, 1992.
- American Society for Testing and Materials. Standard Test Methods for Anions in Water by Chemically Suppressed Ion Chromatography, D4327-97, Vol. 11.01, West Conshohocken, PA, 1999, 420-427.
- Environmental Protection Agency. Fluoride: Dose-Response Analysis for Non-cancer Effects, EPA Report No. 820-R-10-019, Washington, CD, 2010.
- Environmental Protection Agency. Monitored Natural Attenuation of Inorganic Contaminants in Ground Water, EPA Report No. 600-R-07-140, Cincinnati, OH, 2007.
- U.S. EPA Method 300.0, rev 2.1, Determination of Inorganic Anions in Water by Ion Chromatography, U.S. Environmental protection Agency, Cincinnati, OH, 1993.
- U.S. EPA Method 300.1, rev 1.0, Determination of Inorganic Anions in Drinking Water by Ion Chromatography, U.S. Environmental Protection Agency, Cincinnati, OH, 1997.
- Thermo Fisher Scientific. Technical Note 113, Practical Guidance for Using Capillary Anion Chromatography, Dionex LPN 3043, Sunnyvale, CA, 2012.
- Thermo Fisher Scientific. Technical Note 131, Configuring High-Pressure Capillary IC on the Modular IC System, Document No. TN70352, Sunnyvale, CA, 2012.
- Thermo Fisher Scientific. Dionex ICS-5000+ Ion Chromatography System Installation Instructions. Document No. 065447, Sunnyvale, CA, 2012.
- Thermo Fisher Scientific. Dionex Product Manual for the Continuously Regenerated Trap Column (CR-TC). Doc No. 031910, Sunnyvale, CA, 2010.
- Thermo Fisher Scientific. CES 300 Product Manual. Document No. 065386, Sunnyvale, CA, 2012.
- Thermo Fisher Scientific. AS-AP Operator's Manual. Document No. 065361, Sunnyvale, CA, 2012.
- Thermo Fisher Scientific. Application Note 130, Fast Determinations of Inorganic Ions in Salton Sea Samples Using a High-Pressure IC System, Document No. TN70351, Sunnyvale, CA, 2013.
- Thermo Fisher Scientific. Application Note 154, Determination of Inorganic Anions in Environmental Waters Using a Hydroxide-Selective Column, Dionex LPN 1539, Sunnyvale, CA, 2003.

www.thermofisher.com/chromatography

©2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific Inc. 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 representative for details.

Africa +43 1 333 50 34 0
Australia +61 3 9757 4300
Austria +43 810 282 206
Belgium +32 53 73 42 41
Brazil +55 11 3731 5140
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 6 6885 1213
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
Sweden +46 8 556 468 00
Switzerland +41 61 716 77 00
Taiwan +886 2 8751 6655
UK/Ireland +44 1442 233555
USA +1 800 532 4752

Thermo
 S C I E N T I F I C

A Thermo Fisher Scientific Brand