

Cost-Effective Determination of Inorganic Anions and Cations in Municipal Drinking Water Using Capillary Ion Chromatography

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

Introduction

The determination of common inorganic anions and cations in drinking water is one of the most important applications of ion chromatography (IC) worldwide. IC has been approved for compliance monitoring of inorganic anions in United States (U.S.) drinking water since the mid-1980s, as described in U.S. Environmental Protection Agency (EPA) Method 300.0.¹ Many other industrialized countries have similar health and environmental standards and a considerable number of regulatory IC methods have been published worldwide (e.g., in Germany, France, Italy, and Japan). 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 determination of inorganic anions in drinking water.^{2,3} The concentration of some anions in drinking water are regulated due to their toxic effects. For example, high levels of fluoride can cause skeletal and dental fluorosis, and nitrite and nitrate can cause methemoglobinemia, which can be fatal to infants. Other common anions, such as chloride and sulfate, are considered secondary contaminants and can affect odor, color, and certain aesthetic characteristics in drinking water.

IC methods for dissolved alkali and alkaline earth metals and ammonia in drinking water are also important. Drinking water is frequently monitored for the presence of sodium under the U.S. EPA Safe Drinking Water Act. Ammonium is commonly a required target analyte for wastewater discharge permits, and is monitored in process wastewaters.

This study describes the determination of inorganic anions and cations in drinking water using the Thermo Scientific™ Dionex™ ICS-5000 capillary IC system. Scaling down from standard bore to capillary scale brings many benefits to IC analysts. Capillary Thermo Scientific Dionex Reagent-Free™ IC systems deliver fast turnaround from sample submission to results by reducing eluent preparation, system startup, and equilibration times. Perhaps most importantly, the system can be left on and ready for analysis at any time because of its low consumption of eluent (15 mL of source water a day). Having the system always on and ready for analysis significantly streamlines the workflow in IC. An always-on system maintains stability and requires less frequent calibrations. The amount of waste generated is significantly decreased and the eluent generation cartridge producing the eluent lasts 18 months under continuous operation mode, which translates into reduced overall cost of ownership.

Figure 1 shows the determination of inorganic anions in drinking water using capillary IC. The inorganic anions were separated on the Thermo Scientific Dionex IonPac™ AS19 capillary column and detected by suppressed conductivity detection. All anions were separated and eluted within 13 min. The relative standard deviation of peak area for each analyte was 0.6% when 60 injections were evaluated within 24 h.

Figure 2 shows the determination of inorganic cations in drinking water using capillary IC. The inorganic cations were separated on the Dionex IonPac CS12A capillary column and detected by suppressed conductivity detection. All cations of interests were determined within 12 min.

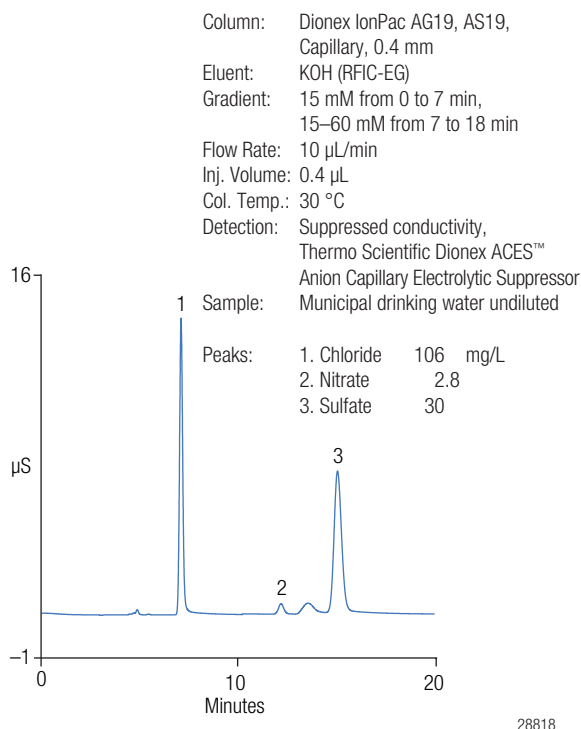


Figure 1. Anions in municipal drinking water on the Dionex IonPac AS19 capillary column.

Conditions

The Dionex ICS-5000 capillary system, Thermo Scientific Dionex AS-AP Autosampler, and Thermo Scientific Dionex Chromeleon™ Chromatography Data System software are used in this experiment. All experimental parameters are listed in Figures 1 and 2.

Sample Preparation

Analyze municipal drinking water by capillary IC without sample pretreatment.

Conclusion

The introduction of the capillary Reagent-Free IC systems redefine the IC workflow for determination of inorganic anions and cations, providing enhanced mass sensitivity and ease of use.⁴ These systems are a great solution for routine characterization of water samples with the always on, always ready capability simplifying the overall IC workflow.

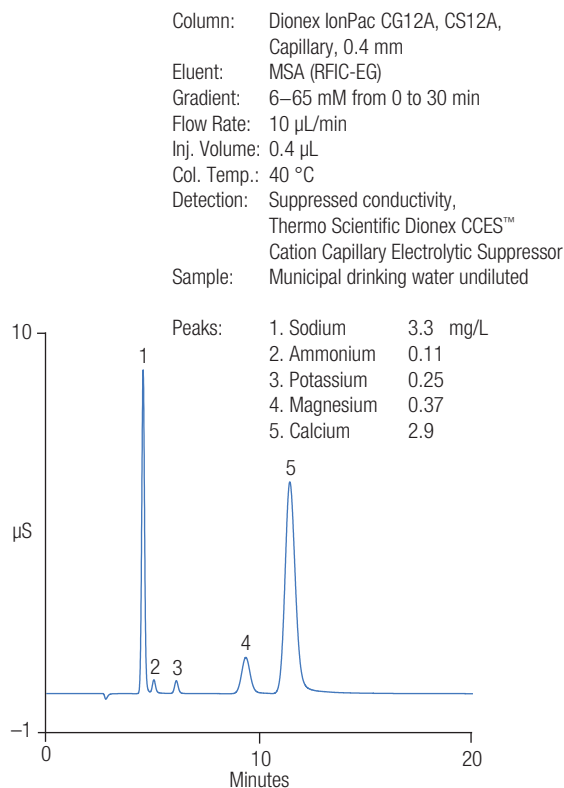


Figure 2. Cations in municipal drinking water on the Dionex IonPac CS12A capillary column.

References

1. *The Determination of Inorganic Anions in Water by Ion Chromatography; Method 300.0*, Revision 2.1; U.S. Environmental Protection Agency: Cincinnati, OH, 1993.
2. Greenberg, A. E.; Clesceri, L. S.; Eaton, A. D., Eds.; *Standard Methods for the Examination of Water and Wastewater*, 18th ed.; American Public Health Association: Washington, DC, 1992.
3. *Standard Test Methods for Anions in Water by Chemically Suppressed Ion Chromatography*; D4327-97, Vol. 11.01; American Society for Testing and Materials; West Conshohocken, PA; 1999; pp 420–427.
4. Dionex Corporation (now part of Thermo Scientific), *Mass Sensitivity of Capillary IC Systems Explained*. Technical Note 90, LPN 2649, Sunnyvale, CA, 2011.

www.thermofisher.com/dionex

©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 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.

Australia +61 3 9757 4486
Austria +43 1 333 50 34 0
Belgium +32 53 73 42 41
Brazil +55 11 3731 5140
China +852 2428 3282

Denmark +45 70 23 62 60
France +33 1 60 92 48 00
Germany +49 6126 991 0
India +91 22 2764 2735
Italy +39 02 51 62 1267

Japan +81 6 6885 1213
Korea +82 2 3420 8600
Netherlands +31 76 579 55 55
Singapore +65 6289 1190
Sweden +46 8 473 3380

Switzerland +41 62 205 9966
Taiwan +886 2 8751 6655
UK/Ireland +44 1442 233555
USA and Canada +847 295 7500

Thermo
 SCIENTIFIC

Part of Thermo Fisher Scientific