

Now sold under the
Thermo Scientific brand

Thermo
SCIENTIFIC

Determination of Nitrite and Nitrate in Drinking Water Using Chemically Suppressed Ion Chromatography

INTRODUCTION

Ion chromatography provides a convenient method for the determination of common inorganic anions in drinking water, including nitrite and nitrate. Using chemically suppressed ion chromatography, trace nitrite and nitrate are accurately and rapidly determined in drinking water. This includes water samples that have been preserved with sulfuric acid. The method will tolerate chloride levels up to 150 ppm and sulfuric acid concentrations as high as 0.5%. For most drinking water samples, the addition of 0.1% by volume sulfuric acid will acidify the sample to $\text{pH} < 2$. This ion chromatographic method, using an IonPac[®] AS9, will tolerate levels of sulfate up to five times that amount.

Figure 1 shows a chromatogram of drinking water spiked to 3.3 ppm nitrite and 22.0 ppm nitrate. The sample is preserved with sulfuric acid to $\text{pH} < 2$.

RECOMMENDED EQUIPMENT

Any Dionex Ion Chromatograph with a conductivity detector.

CONDITIONS

Column: IonPac AS9 analytical with
IonPac AG9 guard
Eluent: 1.8 mM Sodium carbonate
1.7 mM Sodium bicarbonate
Flow Rate: 2.0 mL/min

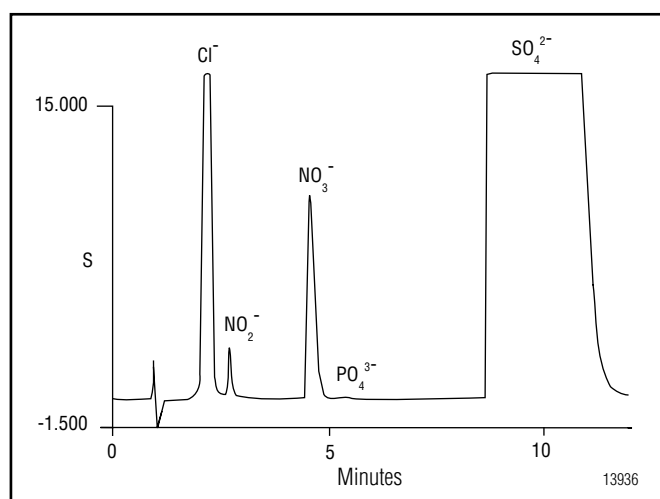


Figure 1 Nitrite and nitrate in acid preserved drinking water.

Suppressor: Anion MicroMembrane[™]
Suppressor (AMMS[®])
Regenerant: 25 mN Sulfuric acid
Regenerant Flow Rate: 6 mL/min using AutoRegen[®]
accessory
Sample Volume: 25 μL
Detection: Suppressed conductivity,
10 μS full scale

REAGENT AND STANDARD PREPARATION

Sodium carbonate/sodium bicarbonate concentrate

(P/N 39513)

Suppressor regenerant concentrate

(4 pack, P/N 37164)

Sodium nitrite, ACS Grade

Sodium nitrate, ACS Grade

ELUENT

To prepare 1.0-L of eluent, dilute 10.0 mL of eluent concentrate to 1000 mL with deionized water.

REGENERANT

To prepare 1.0-L of regenerant, dilute regenerant concentrate to 1000 mL with deionized water as directed on the regenerant concentrate label.

STOCK STANDARDS

1000 ppm Nitrite – Dissolve 1.499 g of sodium nitrite in

1.0 L of deionized water. 1000 ppm Nitrate – Dissolve

1.371 g sodium nitrate in 1.0 L of deionized water.

WORKING STANDARDS

Dilute the stock standards to concentration levels that bracket the concentration level of interest. Prepare working standards fresh from the stock standard prior to analysis.

PERFORMANCE CHARACTERISTICS

Detection limits in drinking water samples using a 25- μ L loop are 10 ppb for nitrite and nitrate. This corresponds to 3.0 ppb nitrogen as nitrite, and 2.3 ppb nitrogen as nitrate. The method is linear for nitrite over the range 10 ppb to 50 ppm. It is linear for nitrate over the range of 20 ppb to 100 ppm.

MicroMembrane is a trademark and AMMS, AutoRegen, and IonPac are registered trademarks of Dionex Corporation.



Printed on recycled and recyclable paper.

Dionex Corporation

1228 Titan Way
P.O. Box 3603
Sunnyvale, CA
94088-3603
(408) 737-0700

Dionex Corporation

Salt Lake City Technical Center
1515 West 2200 South, Suite A
Sunnyvale, CA
84119-1484
(801) 972-9292

Dionex U.S. Regional Offices

Sunnyvale, CA (408) 737-8522
Westmont, IL (630) 789-3660
Houston, TX (281) 847-5652
Atlanta, GA (770) 432-8100
Marlton, NJ (609) 596-0600

Dionex International Subsidiaries

Austria (01) 616 51 25 Belgium (015) 203800 Canada (905) 844-9650 France 01 39 46 08 40 Germany 06126-991-0
Italy (06) 66030052 Japan (06) 885-1213 The Netherlands (0161) 6885-1213 Switzerland (062) 205 99 66 United Kingdom (01276) 691722
* Designed, developed, and manufactured under an NSAI registered ISO 9001 Quality System.

<http://www.dionex.com>



LPN 034504-02 2M 7/98
© 1998 Dionex Corporation