

Now sold under the
Thermo Scientific brand

Thermo
SCIENTIFIC

Determination of Nitrate and Sulfate Collected on Air Filters

PREFACE

Since the initial development of this Application Note in May 1978, there have been several updates in suggested IC conditions. These updates have improved the speed and resolution of IC. The latest recommended conditions from Dionex are listed below. See Figure 3.

Column:	IonPac® AS14 Analytical and AG14 Guard
Eluent:	3.5 mM Na ₂ CO ₃ /1.0 mM NaHCO ₃
Flow Rate:	1.2 mL/min
Suppressor:	Anion Self-Regenerating Suppressor (ASRS), recycle mode
Detection:	Conductivity
Injection Volume:	50 µL

The remainder of this Note is similar to when it was originally published.

INTRODUCTION

The classical method of analysis for nitrate and sulfate is wet chemistry. This report describes a comparison between IC and wet chemistry methods for the determination of nitrate and sulfate collected on filters used in air pollution monitoring studies. The filters were provided by J. Mulik of the United States Environmental Protection Agency (U.S. EPA) in Research Triangle Park, North Carolina.

RECOMMENDED EQUIPMENT

Dionex DX-120 or
DX 500 Ion Chromatography System consisting of:
GP40 Gradient Pump
CD20 Conductivity Detector or
ED40 Electrochemical Detector

Autosampler
PeakNet Chromatography Workstation

PROCEDURE

Four 20- x 1.5-cm strips of labeled filter paper were submitted for analysis. Each contained a known weight of nitrate and sulfate. Similar strips were analyzed by wet chemistry techniques for nitrate and sulfate.

Filter strips 6054 and 6052 contained 465 mg of sulfate and 1381 mg of nitrate determined by known addition. The following procedure was used to extract and analyze the strips for nitrate and sulfate: Immerse each strip in 50.0 mL of a buffer (or eluent) at 50 ± 3 °C for 7 minutes, agitating the solution during this period. (*Note:* More recently the eluent used has been 0.003 M NaHCO₃/0.0024 M Na₂CO₃ and the extractions were performed by letting the filter strips stand for 10 minutes or sonicating them for 10 minutes at room temperature without heat. Decant the buffer and inject directly into a Dionex Ion Chromatograph.

The strips were allowed to stay in contact with the buffer for an additional 4 hours at room temperature (covered to prevent evaporation) and then analyzed to ensure that all of the nitrate and sulfate had been extracted.

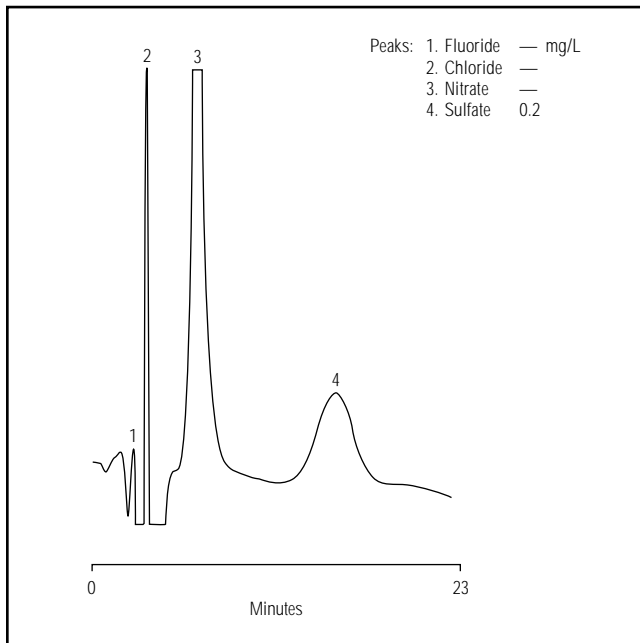


Figure 1 EPA Standard

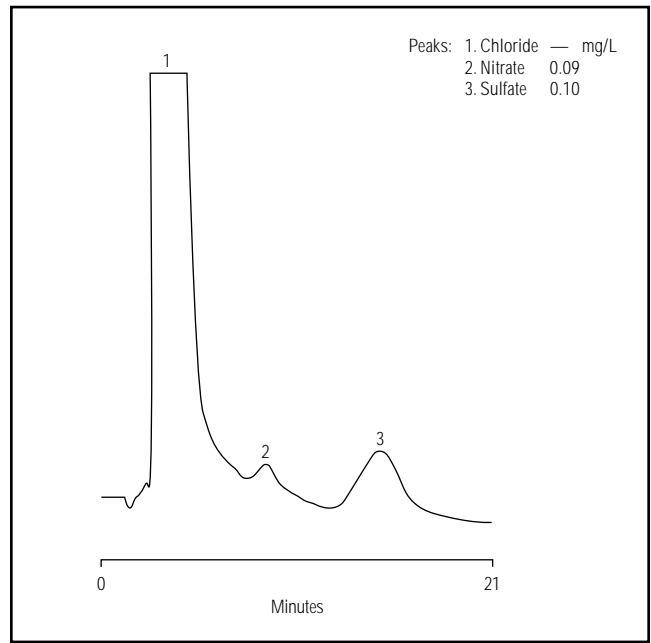


Figure 2 Sample 5163

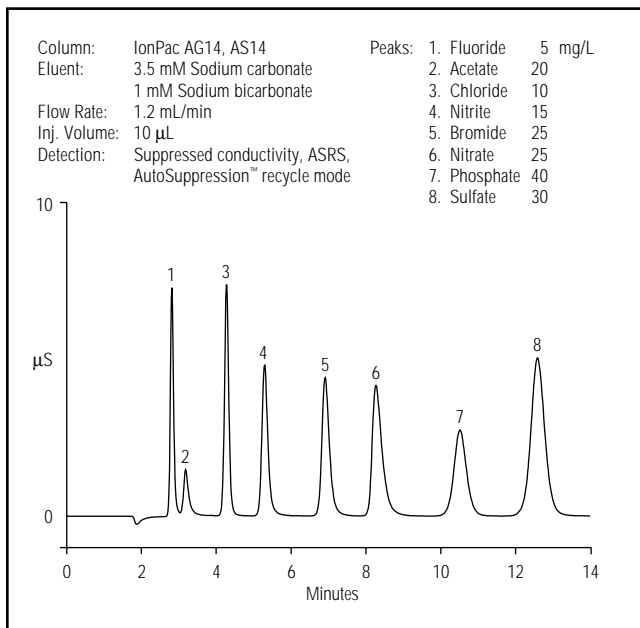


Figure 3 Anion standards chromatographed on the IonPac AS14 anion-exchange column

CONDITIONS

Eluent:	1.8 mM Na ₂ CO ₃ /1.7 mM NaHCO ₃
Flow Rate:	2.0 mL/min
Column:	IonPac AS1 (3 x 250 mm)
Detection:	Conductivity
Suppressor:	AMMS

Table 1 gives the results of the analysis for nitrate and sulfate. Notice that it would be straightforward to determine fluoride and chloride in addition to nitrate and sulfate.

The data clearly show that determination of sulfate by IC is more accurate than by wet chemistry methods. It is known that chloride ion interferes to give a positive variance in the barium chloranilate analysis¹. Sample 6054 contains over 138 mg/L chloride, suggesting one source of error. The sulfate results for samples 5162 and 5163 indicate an even larger error for the wet chemistry method, as the concentration of sulfate is reduced. The nitrate determination by IC compares favorably with wet chemistry, particularly at the higher concentration. No increase in nitrate or sulfate levels was observed even after soaking the filters in buffer for four additional hours. This, along with the good overall experimental reproducibility, implies that the extraction is complete.

CONCLUSION

IC offers a convenient and more accurate method of sulfate analysis than wet chemical methods. Also, nitrate analysis by IC is comparable in accuracy and faster than wet chemistry methods (total analysis time was 20 minutes). Since fluoride, chloride, nitrate, sulfate, and other ions such as nitrite, phosphate, and acetate may be simultaneously analyzed, IC represents a powerful analytical tool for the rapid and accurate analysis of anions for air pollution.

REFERENCE

1. Personal communication, Jim Mulik, U.S. EPA.

Table 1 Analysis of Filter Extracts

Samples	Nitrate Concentration (mg/L)			Sulfate Concentration (mg/L)		
	IC ¹	Wet Chemistry ²	Known	IC ¹	Wet Chemistry ²	Known
6052, 6054	25.4 ± 0.6 ³	27.5	27.6	9.25 ± 0.5 ³	13.6	9.3
5162, 5163	0.09 ± 0.005	0.45	—	0.1 ± 0.01	11.6	—

¹ Average of the average of each sample. Each individual analysis was repeated at least three times giving a relative standard deviation of 3% for 6054 and 6052. The RSD for 5162 and 5163 was not calculated but was higher due to poorer signal to noise. Only one analysis per sample was made in this case.

² Error analysis is unknown.

³ No increase in concentration was observed after 4 hours of soaking the filter in buffer.

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
Salt Lake City, UT
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 (856) 596-06009

Dionex International Subsidiaries
Austria (01) 616 51 25 *Belgium* (32) 3-353 42 94 *Canada* (905) 844-9650 *China* (852) 2428 3282 *Denmark* (45) 36 36 90 90
France 01 39 30 01 10 *Germany* 06126-991-0 *Italy* (06) 66 51 50 52 *Japan* (06) 6885-1213 *The Netherlands* (0161) 43 43 03
Switzerland (062) 205 99 66 *United Kingdom* (01276) 691722
* Designed, developed, and manufactured under an NSAI registered ISO 9001 Quality System.

