

Direct Determination of **Existent Chloride** and **Sulfate** in **Denatured Ethanol Samples** Using a **Compact Ion Chromatography System**

Terri Christison and Jeff Rohrer, Thermo Fisher Scientific, Sunnyvale, CA, USA

Key Words

HPIC, Integrion, Dionex IonPac AS22 Column, Dionex AS22 Eluent Concentrate, Dionex AERS 500 Suppressor, Biofuel

Introduction

This application proof note demonstrates the direct determination of existent chloride and sulfate in denatured ethanol samples, as based on the method published in [Application Update 194](#).¹ In this proof note, the method is performed using a Thermo Scientific™ Dionex™ Integrion™ HPIC™ system.

Method

IC System:	Thermo Scientific Dionex Integrion HPIC system with column heater
Columns:	Thermo Scientific™ Dionex™ IonPac™ AS22 Analytical (4 × 250 mm) Thermo Scientific Dionex IonPac AG22 Guard (4 × 50 mm)
Eluent:	4.5 mM Sodium Carbonate/1.4 mM Bicarbonate
Flow Rate:	1.2 mL/min
Injection Volume:	25 µL
Temperature:	30 °C
Detection:	Suppressed conductivity, Thermo Scientific™ Dionex™ AERS™ 500 Carbonate Suppressor, 4 mm, 40 mA, recycle mode

Reference

1. Thermo Scientific Application Note 194: Determination of Existent and Potential Sulfate and Total Inorganic Chloride in Denatured Ethanol by Direct Injection Using an RFIC System, Sunnyvale, CA [Online] <http://www.thermoscientific.com/content/dam/tfs/ATG/CMD/CMD%20Documents/Application%20&%20Technical%20Notes/Chromatography/Ion%20Chromatography/IC%20and%20RFIC%20Columns/AU-194-IC-Sulfate-Chloride-Denatured-Ethanol-AN71058-EN.pdf> (accessed Jan. 14, 2016)

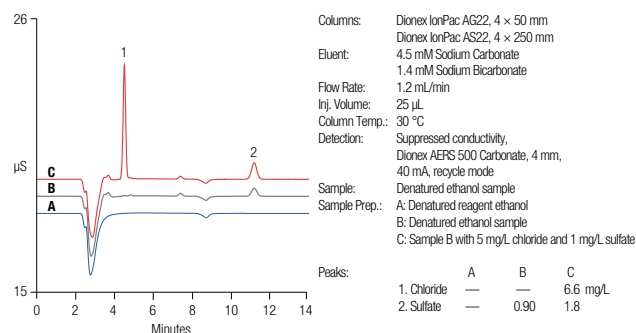


Figure 1. Separation of chloride and sulfate in denatured ethanol samples: A) Denatured reagent ethanol, B) Denatured ethanol sample, C) Sample B with 5 mg/L chloride and 1 mg/L sulfate.

For application support, visit the [AppsLab Library](#) where you can find detailed method information, chromatograms and related compound information. All the information needed to run, process and report the analysis is available in ready-to-use eWorkflows, which can be executed directly in your chromatography data system. www.thermoscientific.com/appslib



www.thermoscientific.com/integrion

©2016 Thermo Fisher Scientific Inc. All rights reserved. ISO is a trademark of the International Standards Organization. All other trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific 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.

AB71907-EN 0116S



Thermo Fisher Scientific,
Sunnyvale, CA USA is
ISO 9001 Certified.

Thermo
SCIENTIFIC

A Thermo Fisher Scientific Brand