# Determination of Ammonia in Sodium Bicarbonate Using a Compact Ion Chromatography System

Carl Fisher, Thermo Fisher Scientific, Sunnyvale, CA, USA

## **Key Words**

Integrion, IonPac CS16, EGC 500, Suppressed Conductivity, Hemodialysis Solution, Ion Chromatography

#### Introduction

One of the main pharmaceutical uses of sodium bicarbonate is for hemodialysis, and the quality and purity requirements for hemodialysis-grade sodium bicarbonate are stringent. Ammonia is one impurity that is assayed in sodium bicarbonate because the presence of elevated levels of ammonia in the blood stream can have adverse effects. This application proof note demonstrates the determination of ammonia (as ammonium) in sodium bicarbonate, using the method published in Application Note 1073.¹ In this proof note, the method is performed using a Thermo Scientific™ Dionex™ Integrion™ ion chromatography system.

### Method

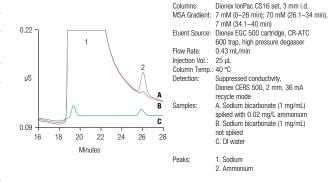
Method				
IC System:	Thermo Scientific Dionex Integrion IC system			
Columns:	Thermo Scientific Dionex IonPac CS16 Analytical (3 $\times$ 250 mm) Thermo Scientific Dionex IonPac CG16 Guard (3 $\times$ 50 mm)			
Eluent:	Methanesulfonic acid			
Gradient:	Retention Time [min]	Flow [mL/min]	Concentration [mM]	
	0.0	0.43	7.0	
	26.0	0.43	7.0	
	26.1	0.43	70.0	
	34.0	0.43	70.0	
	34.1	0.43	7.0	
	40.0	0.43	7.0	

Flow Rate:	0.43 mL/min		
Injection Volum	e: 25 μL		
Temperature:	40 °C		
Detection: Suppressed conductivity, Thermo Scientific™ Dionex™ CERS™ 500 (2 Electrolytically Regenerated Suppressor			

#### Reference

1. Thermo Scientific Application Note 1073: Determination of Ammonia in Sodium Bicarbonate. Sunnyvale, CA [Online] <a href="http://www.thermoscientific.com/content/dam/tfs/ATG/CMD/CMD%20Documents/AN-1073-Determination-Ammonia-Sodium-Bicarbonate-AN70815-E.pdf">http://www.thermoscientific.com/content/dam/tfs/ATG/CMD/CMD%20Documents/AN-1073-Determination-Ammonia-Sodium-Bicarbonate-AN70815-E.pdf</a> (accessed Jan. 8, 2016)





 $\label{lem:figure 1.Determination of ammonia in sodium bicarbonate.}$ 

For application support, visit the <u>AppsLab Library</u> 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. <a href="https://www.thermoscientific.com/appslab">www.thermoscientific.com/appslab</a>





©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.



