

IC Assay for Lithium, Sodium, and Calcium in Lithium Carbonate Using a Compact Ion Chromatography System

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

Key Words

Integriion, IonPac CS16, EGC 500, Cationic Impurities

Introduction

Lithium carbonate is used to treat a number of mental health problems associated with chemical imbalances in the brain. Sodium and calcium are possible cationic impurities in lithium carbonate preparations. This application proof note shows the determination of lithium, sodium, and calcium in lithium carbonate, using the method published in Application Note 1090.¹ In this proof note, the method is performed using a Thermo Scientific™ Dionex™ Integriion™ ion chromatography system in combination with a Thermo Scientific™ Dionex™ IonPac™ CS16 cation-exchange column, an electrolytically generated MSA eluent, and suppressed conductivity detection.

Method

IC System: Thermo Scientific Dionex Integriion IC system

Columns: Thermo Scientific Dionex IonPac CS16 Analytical (3 × 250 mm)
Thermo Scientific Dionex IonPac CG16 Guard (3 × 50 mm)

Eluent: Methanesulfonic acid

Gradient:	Retention Time [min]	Flow [mL/min]	Concentration [mM]
	0.0	0.43	8.0
	15.0	0.43	8.0
	15.0	0.43	67.0
	20.0	0.43	67.0
	20.0	0.43	8.0
	25.0	0.43	8.0

Flow Rate: 0.43 mL/min

Injection Volume: 10 µL

Temperature: 40 °C

Detection: Suppressed conductivity, Thermo Scientific™ Dionex™ CERS™ 500 (2 mm)
Electrolytically Regenerated Suppressor

Reference

1. Thermo Scientific Application Note 1090: IC Assay for Lithium, Sodium, and Calcium in Lithium Carbonate. Sunnyvale, CA [Online] <http://www.thermoscientific.com/content/dam/tfs/ATG/CMD/CMD%20Documents/Application%20&%20Technical%20Notes/Chromatography/Ion%20Chromatography/AN-1090-IC-Lithium-Sodium-Calcium-Lithium-Carbonate-AN71001-EN.pdf> (accessed Jan. 8, 2016)

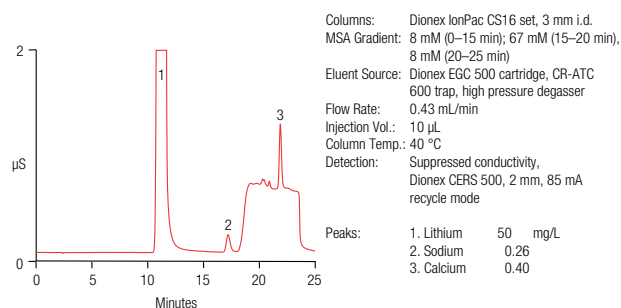
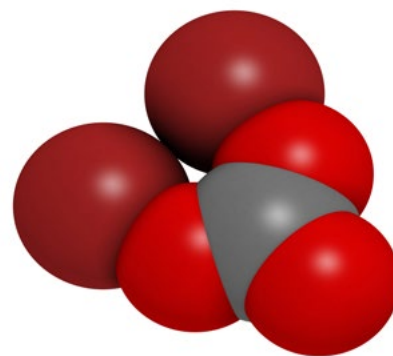


Figure 1. Lithium, sodium, and calcium in lithium carbonate.

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/appslab



www.thermoscientific.com/integriion

©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.

AB71919-EN 0116S



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

Thermo
SCIENTIFIC

A Thermo Fisher Scientific Brand