

# Determination of High Sugar Concentrations in a Rice Wine Sample Using a Compact Ion Chromatography System

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

## Key Words

HPAE-PAD, Dionex Integrion RFIC System, Dionex CarboPac PA20 Column, Dionex EGC 500 KOH Eluent Generator, Alcohol

## Introduction

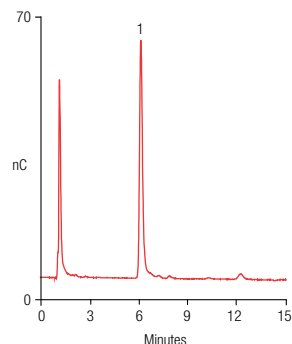
This application proof note demonstrates the determinations of glucose and fructose in a 100-fold diluted rice wine sample by HPAE-PAD. Typically, samples with g/L concentrations require greater than 10,000-fold dilutions to remain in the linear range of the very sensitive HPAE-PAD technique. However, in this proof note, the method is performed using a Thermo Scientific™ Dionex™ Integrion™ RFIC™ system equipped with a 0.4  $\mu$ L internal injection loop and the Thermo Scientific™ High Concentration Carbohydrate Analysis Kit<sup>1</sup> to extend the linearity from low mg/L to g/L concentrations. A wood hydrolysate application using this technique is available for reference.<sup>2</sup>

## Method

IC System:	Thermo Scientific Dionex Integrion RFIC system with column heater
Columns:	Thermo Scientific™ Dionex™ CarboPac™ PA20 Analytical (3 $\times$ 150 mm) Thermo Scientific Dionex CarboPac PA20 Guard (3 $\times$ 30 mm)
Eluent:	35 mM KOH with 100 mM KOH wash
Flow Rate:	0.5 mL/min
Injection Volume:	0.4 $\mu$ L
Temperature:	30 °C
Detection:	Pulsed amperometry, Au on PTFE, 62 mil gasket

## Reference

1. Thermo Scientific Product Specification 70749: Thermo Scientific High Concentration Carbohydrate Analysis Kit. Sunnyvale, CA [Online] <https://www.thermoscientific.com/content/dam/tfs/ATG/CMD/CMD%20Documents/Product%20Manuals%20&%20Specifications/PS-70749-High-Concentration-Carbohydrate-Analysis-Kit-PS70749-EN.pdf> (accessed Jan. 8, 2016)
2. Thermo Scientific Application Note 1089: Determination of Carbohydrates in Acid Hydrolysates of Wood. Sunnyvale, CA [Online] <http://www.thermoscientific.com/content/dam/tfs/ATG/CMD/CMD%20Documents/Application%20&%20Technical%20Notes/Chromatography/Ion%20Chromatography/AN-1089-Carbohydrates-Acid-Hydrolysates-Wood-AN70941-EN.pdf> (accessed Jan. 8, 2016)



Column: Dionex CarboPac PA20 guard,  
Dionex CarboPac PA20, 3 mm i.d.  
Eluent: 35 mM KOH, 100 mM KOH wash  
from 15 to 25 min  
Eluent Source: Dionex EGC 500 KOH cartridge,  
Dionex CR-ATC 500 trap column,  
high pressure degasser  
Flow Rate: 0.50 mL/min  
Inj. Volume: 0.4  $\mu$ L  
Column Temp.: 30 °C  
Detection: PAD, Au on PTFE, 62 mil gasket  
Waveform: 4-Potential Carbohydrate  
Ref. Electrode: Ag/AgCl  
Sample Prep.: 100x dilution with deionized water,  
filter, 0.2  $\mu$ m  
Peaks: 1. Glucose 222 mg/L

Figure 1. Glucose in a rice wine sample.

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](http://www.thermoscientific.com/appslib)



[www.thermoscientific.com/integrion](http://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.

AB71913-EN 0116S



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

**Thermo**  
SCIENTIFIC

A Thermo Fisher Scientific Brand