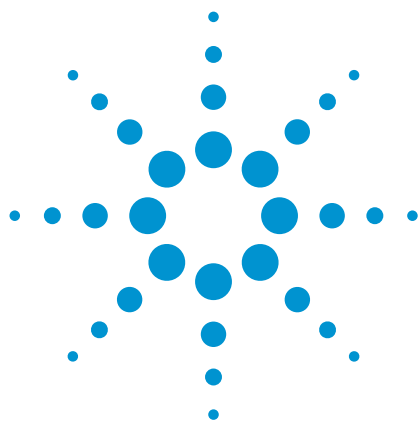




USP Analysis of Sugar Alcohols on an Agilent Hi-Plex Ca Column – Mobile Phase Effects

Application Note

Pharmaceutical

Authors

Stephen Ball, Linda Lloyd
Agilent Technologies, Inc.

Introduction

Sugar alcohols, or polyols, are hydrogenated carbohydrates commonly used to replace sucrose in foods. They are often used with high-intensity artificial sweeteners to counter their low sweetness.

The separation of seven sugar alcohols on an Agilent Hi-Plex Ca column can be altered by introducing acetonitrile into the mobile phase.



Agilent Technologies

Experimental

Conditions

Column	Agilent Hi-Plex Ca USP L19, 4.0 × 250 mm, 8 µm (p/n PL1570-5810)
Mobile phase	100% DI H ₂ O (initially)
Flow rate	0.15 mL/min
Injection volume	10 µL
Temperature	90 °C
Detector	RI

Sample Preparation

The seven sugar alcohols — iso-erythritol, adonitol, arabitol, mannitol, xylitol, dulcitol, and sorbitol — are made up to a concentration of 10 mg/mL in water. See Figure 1.

When pure water is used for the mobile phase, several of the sugar alcohols in the sample either partially or completely co-elute. Modifying operating temperature or flow rate is very unlikely to give a good separation between these compounds.

Introducing acetonitrile into the mobile phase has a significant effect on the selectivity of the Agilent Hi-Plex Ca material and results in a doubling of the retention time. As a result, the mobile phase and flow rate conditions need to be modified as follows:

Conditions

Mobile phase	30:70 acetonitrile:100% DI H ₂ O
Flow rate	0.20 mL/min
Temperature	90 °C

The same quantity of test solution is injected. See Figure 2.

Conclusion

As can be seen by comparing the two chromatograms, using 30% acetonitrile gives extra retention for the sugar alcohols and, as a result, increases the resolution between them. All seven sugar alcohols are now either partially or completely separated. It also gives a change in the elution order.

For More Information

These data represent typical results. For more information on our products and services, visit our Web site at www.agilent.com/chem.

Peak identification

1. Adonitol
2. Iso-erythritol
3. Mannitol, arabitol
4. Dulcitol
5. Sorbitol, xylitol

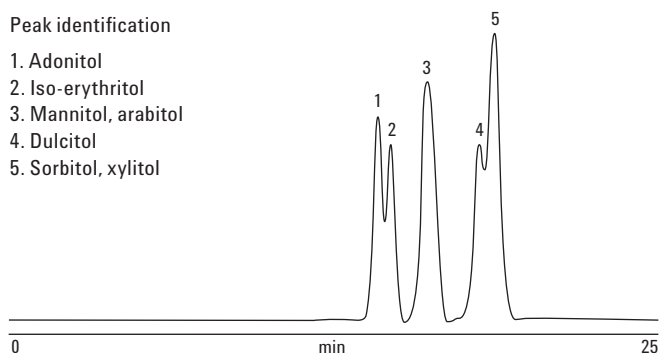


Figure 1. Raw data chromatogram of seven sugar alcohols on an Agilent Hi-Plex Ca USP L19 column.

Peak identification

1. Iso-erythritol
2. Adonitol
3. Arabitol
4. Mannitol
5. Xylitol
6. Dulcitol
7. Sorbitol

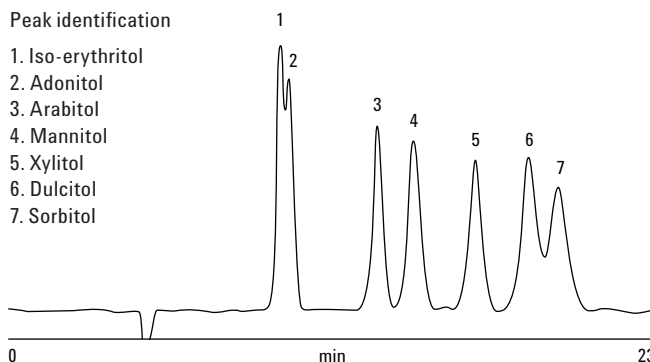


Figure 2. Raw data chromatogram of seven sugar alcohols on an Agilent Hi-Plex Ca USP L19 column after the introduction of acetonitrile into the mobile phase.

www.agilent.com/chem

Agilent shall not be liable for errors contained herein or for incidental or consequential damages in connection with the furnishing, performance, or use of this material.

Information, descriptions, and specifications in this publication are subject to change without notice. This publication was originally published in 2008.

© Agilent Technologies, Inc., 2013
Published in USA, January 11, 2013
SI-01672



Agilent Technologies