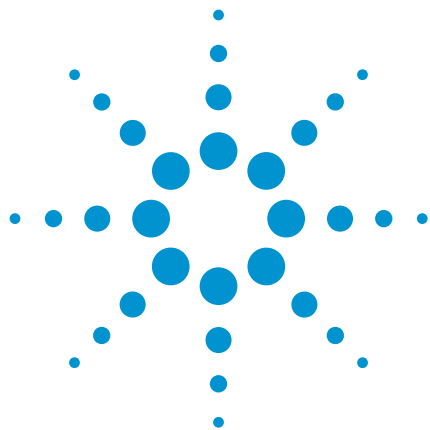




PEG Separations by Aqueous SEC

Agilent PL aquagel-OH 30 8 μm Columns

Technical Overview

Introduction

PL aquagel-OH 30 8 μm high performance columns are extremely hydrophilic and so neutral polymers such as polyethylene oxide/glycol (PEO/PEG) can be run with water as eluent.

Such a separation is illustrated in Figure 1 with the analysis of a range of Agilent PEG standards.

PL aquagel-OH columns are ideal for relatively low molecular weight separations by size exclusion chromatography. They combine high pore volume and high column efficiency for maximum resolution.



Agilent Technologies

Conditions

Columns: 2 x PL aquagel-OH 30 8 μ m,
300 x 7.5 mm (p/n PL1120-6830)

Eluent: Water

Flow Rate: 1 mL/min

Detection: RI

Peak Identification

1. 12,600

2. 4,120

3. 1,470

4. 440

5. 106

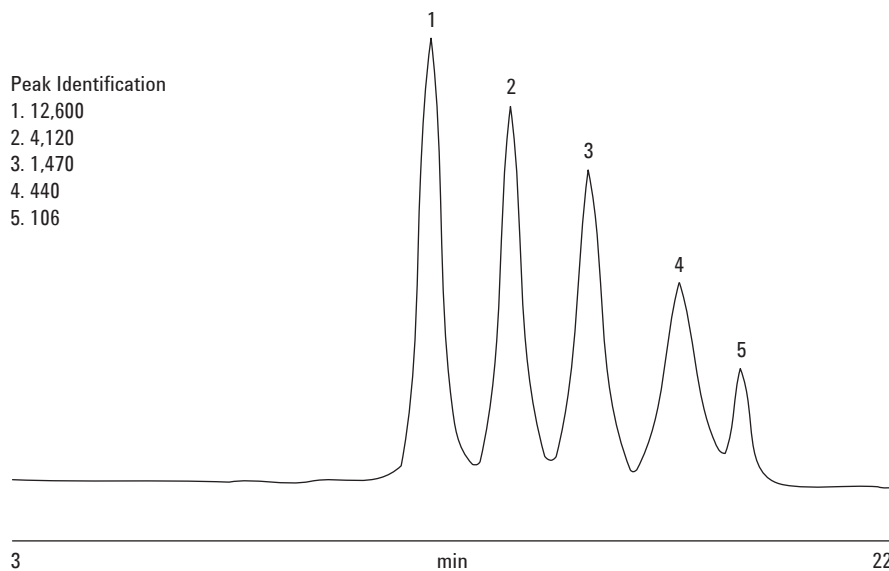


Figure 1. Good separation of a mixture of PEG standards using PL aquagel-OH columns

These data represent typical results. For further information, contact your local Agilent Sales Office.

www.agilent.com/chem

This information is subject to change without notice.

© Agilent Technologies, Inc. 2010

Published in UK, September 20, 2010

SI-01645



Agilent Technologies