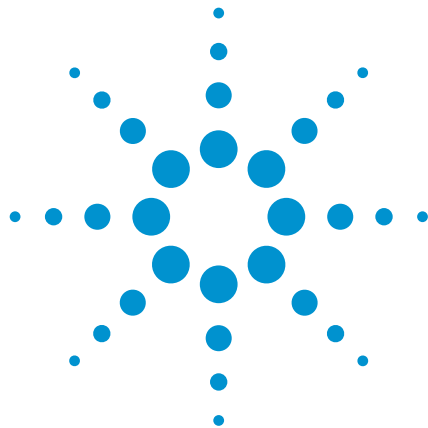




Efficient Analysis of Condensation Resins

Agilent PLgel 5 μ m MIXED-D Columns

Technical Overview

Introduction

The PLgel 5 μ m MIXED-D column is specifically designed for the analysis of polymers, paints and resin systems where material above 400,000 MW is unlikely to be present. High pore volume, combined with the 5 μ m efficiency, provides excellent resolution for low MW polymers and oligomers. Two, or even three, PLgel 5 μ m MIXED-D columns are the perfect replacement for the popular $10^4/500 \text{ \AA}$ or $10^4/10^3/500 \text{ \AA}/100 \text{ \AA}$ column combinations.



Agilent Technologies

A variety of resins, produced by condensation polymerization, can be analyzed using three columns.

Column: 3 x PLgel 5 μ m MIXED-D, 300 x 7.5 mm
(part number PL1110-6504)
Eluent: THF
Flow Rate: 1.0 mL/min
Detection: RI

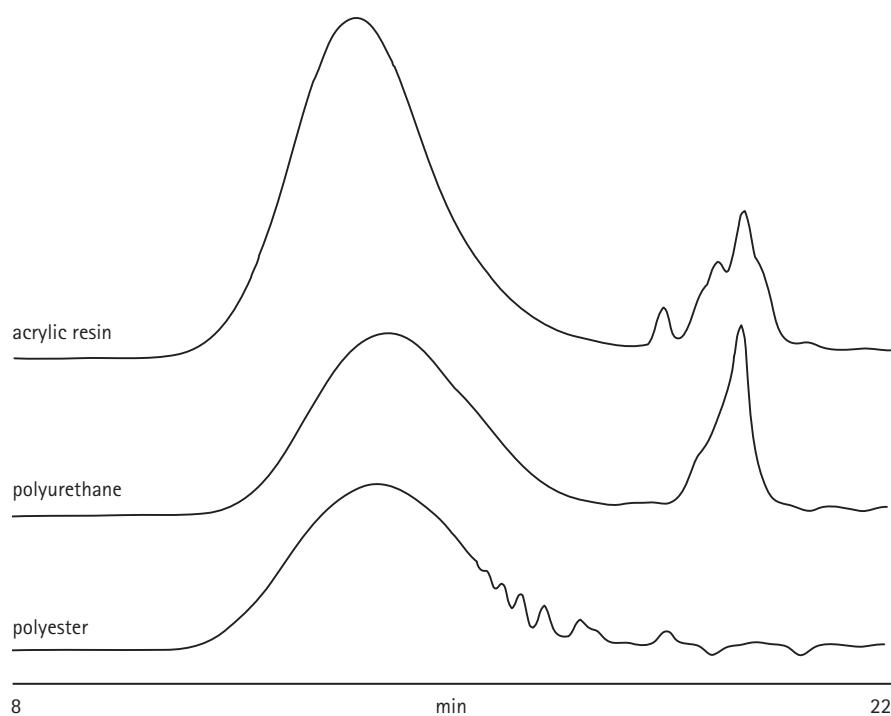


Figure 1. Analysis of resins using PLgel 5 μ m MIXED-D columns

www.agilent.com/chem

This information is subject to change without notice.

© Agilent Technologies, Inc. 2010

Published in UK, September 21, 2010

SI-01828



Agilent Technologies