



Successful GPC Calibration Using Polyethylene Oxide

Agilent PLgel Individual Pore Size Columns

Technical Overview

Introduction

As part of the PLgel family, the individual pore size GPC columns offer high resolution over a specific molecular weight range. The linear portion of the calibration curve, where the slope is at its shallowest, defines the molecular weight region over which optimum resolution will be achieved.



Agilent Technologies

Calibration curves show the successful use of polyethylene oxide standards for GPC calibration in DMF on PLgel 5 μm 500 \AA columns. The slight difference between the calibration curves is due to differences in hydrodynamic volume for each polymer/solvent system.

Conditions

Column: PLgel 5 μm 500 \AA , 300 x 7.5 mm
(part number PL1110-6525)

Flow Rate: 1.0 mL/min

Detection: RI

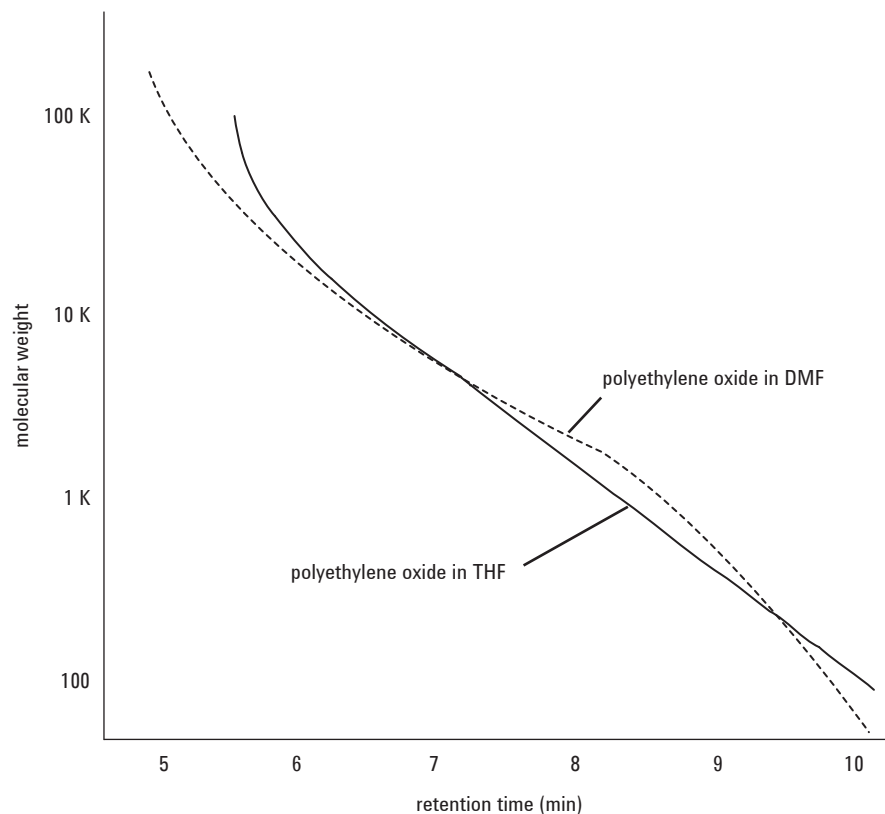


Figure 1. Calibration curves show the successful use of polyethylene oxide standards for GPC calibration in DMF on PLgel 5 μm 500 \AA columns

www.agilent.com/chem

This information is subject to change without notice.

© Agilent Technologies, Inc. 2010

Published in UK, September 23, 2010

SI-01753



Agilent Technologies