

Agilent PLgel 5 μm MIXED-C Columns

GPC Calibrations and Hydrodynamic Volume

Technical Overview

Introduction

When the same column is calibrated with standards of different chemical composition, the shape of the calibration should be similar. The plots will differ according to the hydrodynamic volume of each polymer in the element used.



Agilent Technologies

This effect is shown here using different standards on a PLgel 5 μm MIXED-C column.

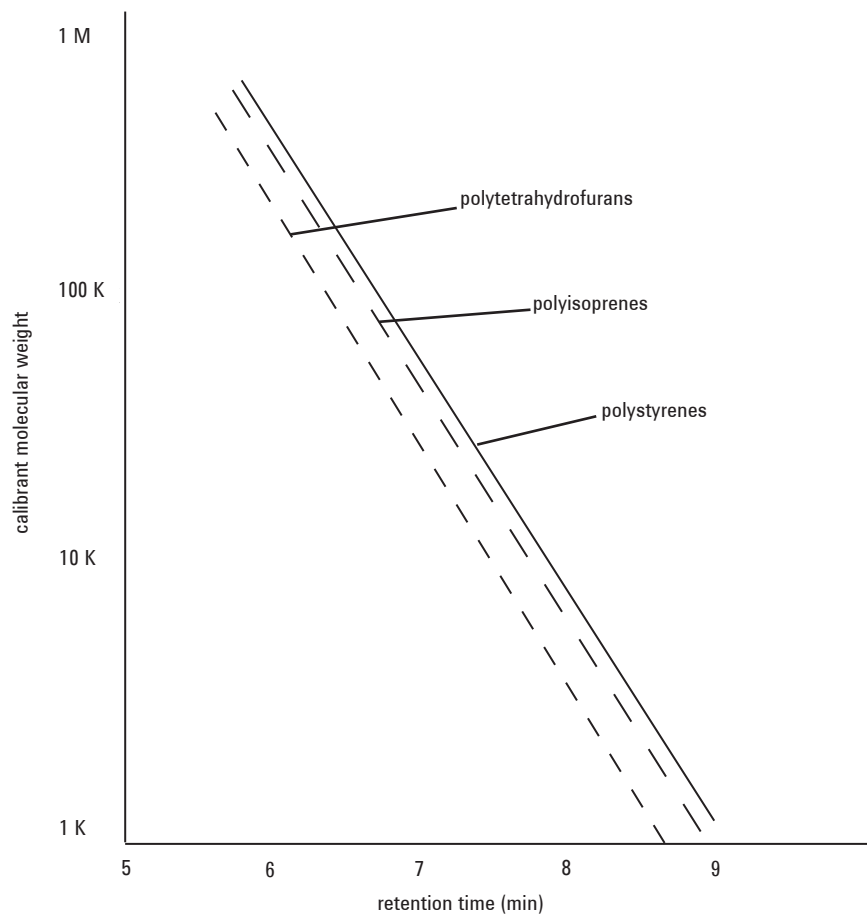


Figure 1. PLgel 5 μm MIXED-C, 300 x 7.5 mm (part number PL1110-6500)

www.agilent.com/chem

This information is subject to change without notice.

© Agilent Technologies, Inc. 2010

Published in UK, September 22, 2010

SI-1790



Agilent Technologies