



Agilent PLgel 5 μ m MIXED-C Columns

Easy GPC Column Calibration Using Polystyrene Standards

Technical Overview

Introduction

Agilent EasiCal PS-1 pre-prepared standards can be used to calibrate PLgel 5 μ m MIXED-C columns. The 8,500,000 MW is excluded as the linear operating range of this column extends up to 2,000,000 MW.



Agilent Technologies

PLgel 5 μm MIXED-C columns are designed for rapid polymer analysis. With its linear calibration up to 2 million MW, this is the column of choice for highest resolution and accuracy in molecular weight distribution analyses. Rapid solvent change capability, excellent temperature stability and the high resolution of the PLgel 5 μm MIXED-C also provide ideal versatility for the modern R&D laboratory.

Conditions

Columns: 2 x PLgel 5 μm MIXED-C,
300 x 7.5 mm (part number
PL11110-6500)
Eluent: THF
Flow Rate: 1.0 mL/min
Detection: UV, 254 nm

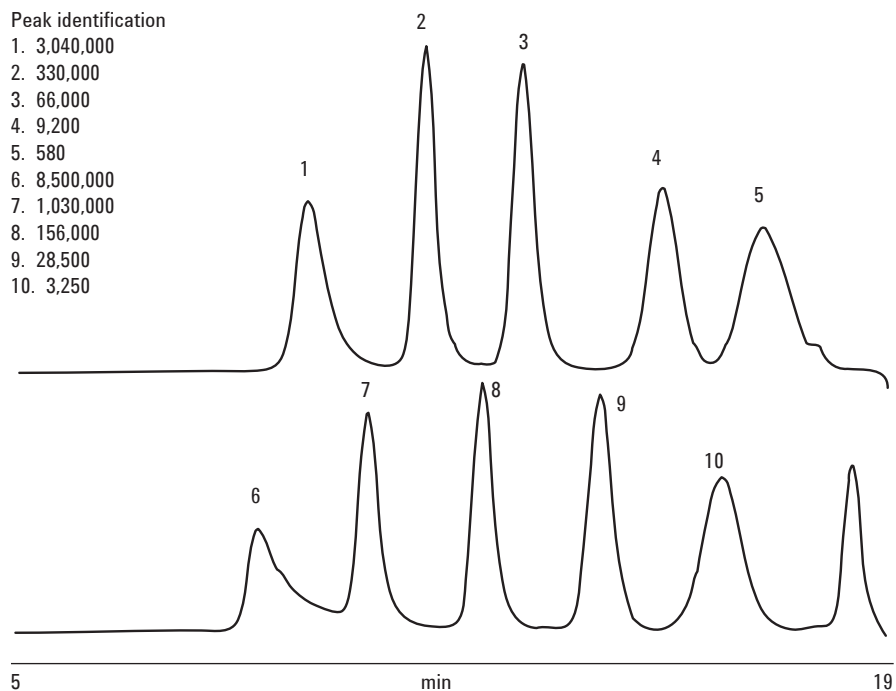


Figure 1. Calibration of PLgel 5 μm MIXED-C columns using EasiCal PS-1 standards

www.agilent.com/chem

This information is subject to change without notice.

© Agilent Technologies, Inc. 2010

Published in UK, September 22, 2010

SI-1788



Agilent Technologies