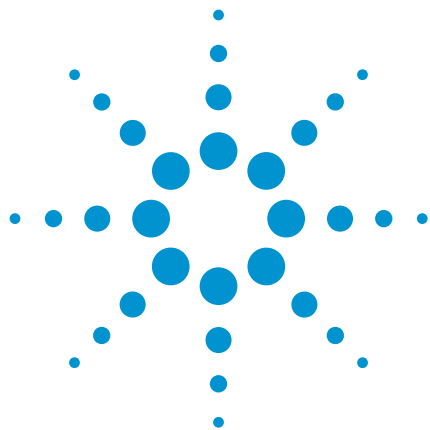




Easy to Use Pre-prepared Calibration Kits

Agilent EasiCal Pre-prepared Calibration Kits

Technical Overview

Introduction

EasiCal PS-1 pre-prepared standards can be used to calibrate Agilent PLgel 10 μm MIXED-B columns. All of the standards are resolved as the column exhibits a linear calibration over the full range of molecular weight covered.



Agilent Technologies

EasiCal polystyrene standards make column calibration easy and convenient. The EasiCal system consists of two different combs, both with ten detachable spatulas supporting a mixture of five polymer standards. The thin film of polymer (approx 5 mg) on the tip of the PTFE spatulas rapidly dissolves when immersed in eluent to provide two GPC/SEC calibration solutions. There are two polystyrene EasiCal versions: PS-1 (Mw range 580-7,500,000) and PS-2 (Mw range 580-400,000). They can be used with any GPC system or column.

Conditions

Columns: 3 x PLgel 10 µm MIXED-B,
300 x 7.5 mm (p/n PL1110-6100)

Eluent: THF

Flow Rate: 1.0 mL/min

Detection: UV, 254 nm

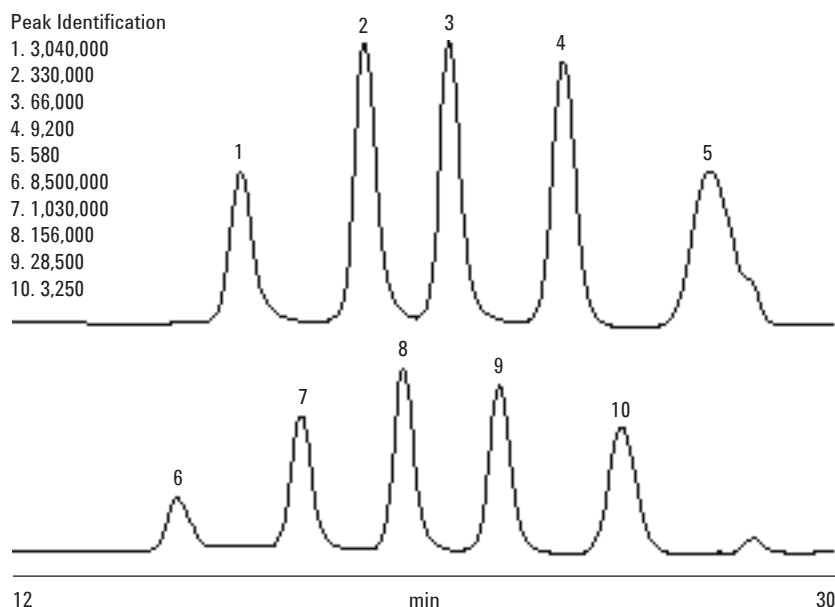


Figure 1. PLgel 10 µm MIXED-B column calibration using EasiCal PS-1 pre-prepared calibrants

www.agilent.com/chem

This information is subject to change without notice.

© Agilent Technologies, Inc. 2010

Published in UK, September 24, 2010

SI-01759



Agilent Technologies