



Agilent EasiCal Pre-prepared Calibration Kits

Fast and Convenient GPC Column Calibration

Technical Overview

Introduction

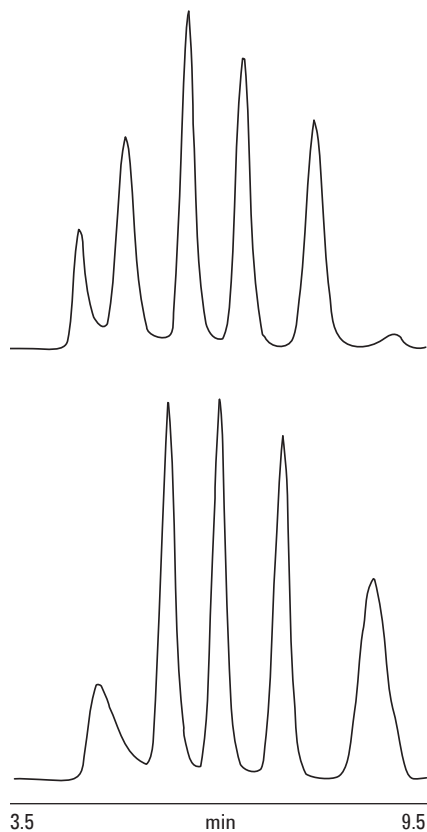
EasiCal polystyrene standards make column calibration easy and convenient. The EasiCal system consists of two different combs, both with ten detachable spatulas supporting a mixture of five polymer standards. The thin film of polymer (approx 5 mg) on the tip of the PTFE spatulas rapidly dissolves when 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 5820-400,000). EasiCal standards can be employed with any GPC system or column. In this example, they have been used with Agilent PLgel MIXED-C columns.



Agilent Technologies

Conditions

Eluent: THF
Flow Rate: 1.0 mL/min
Concentration: 0.05%
Detection: UV, 254 nm



**Figure 1. PLgel 5 μ m 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-1789



Agilent Technologies