

Comparing Polycarbonate Samples with GPC

Agilent PLgel 5 μ m MIXED-C Columns

Technical Overview

Introduction

PLgel 5 μ m MIXED-C columns are designed for rapid polymer analysis. They are ideal for comparing batches of polycarbonate.



Agilent Technologies

With its linear calibration up to 2 million MW, PLgel 5 µm MIXED-C 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 the versatility essential for the modern R&D laboratory.

Conditions

Column: PLgel 5 µm MIXED-C,
300 x 7.5 mm (part number
PL1110-6500)
Eluent: Chloroform
Flow Rate: 1.0 mL/min
Detection: RI

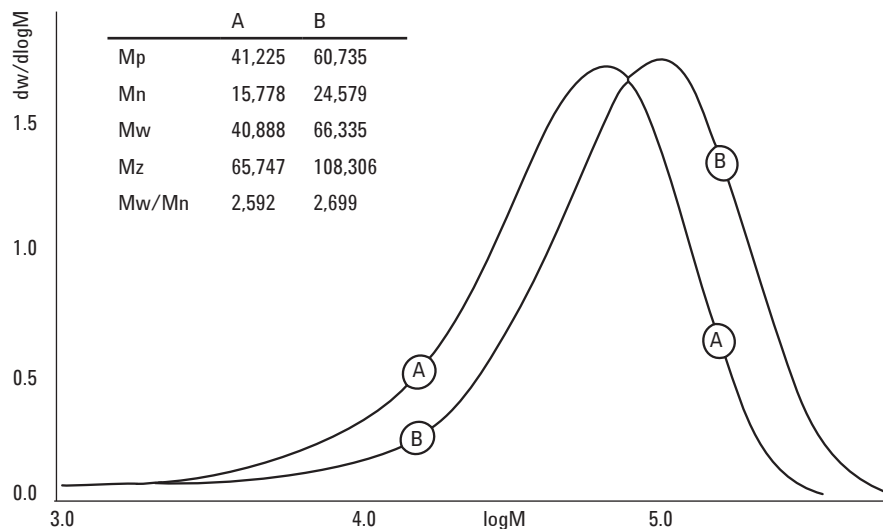


Figure 1. Differences in two batches of polycarbonate revealed by GPC with PLgel MIXED-C columns

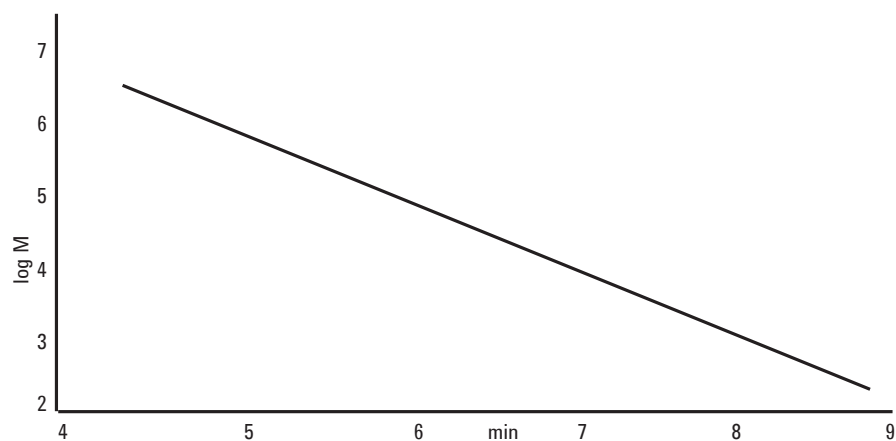


Figure 2. Polystyrene calibration curve

www.agilent.com/chem

This information is subject to change without notice.

© Agilent Technologies, Inc. 2010

Published in UK, September 22, 2010

SI-1792



Agilent Technologies