Errata Notice

This document contains references to PSS or Polymer Standards Service. Please note that PSS is now Agilent. This document will be republished as an Agilent document in the future.





10035 - Column Application Note Characterization of Poly(isobutylene)

Poly(isobutylene) (PIB) can be made from the monomer isobutylene by cationic polymerization. Polyisobutylene is an elastomer and gas impermeable. The first major application of PIB/ butyl rubber (copolymer: 98% isobutylene, 2% of isoprene) was tire inner tubes. PIB (in the form of polyisobutylene succinimide) is used as an additive in lubricating oils and motor

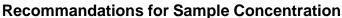
fuels.

Experimental Setup

Mobile Phase: Tetrahydrofuran Stationary Phase: PSS SDV Flow rate [mL/min]: 1,00 Temperature [°C]: 25

Detection: Shodex-RI71

Calibration: Kit Poly(isobutylene)
Data processing: PSS WinGPC



narrow PDI

M 100 Da - 10 000 Da: 2 g/L M 10 000 Da - 1 000 000 Da: 1-2 g/L

M > 1 000 000 Da: 0.5 g/L or less broad PDI (>1.5)

all molar masses: 3.0 - 5.0 g/L

Injection volume [µL]: 20

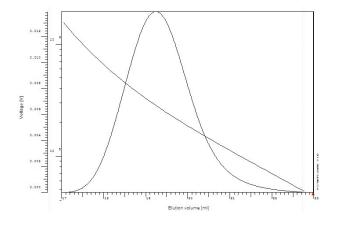


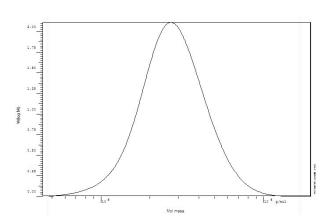
low molecular weights: P/N 201-0001 (set of 3) OR sda083003lis (1 linear) medium molecular weights: P/N 201-0002 (set of 2) OR sda083005lim (1 linear) high molecular weights: P/N 201-0003 (set of 3) OR sda083005lxl (1 linear)

ultrahigh molecular weights: P/N 202-0001 (set of 3)

Elugram and Calibration separation on PSS SDV

Molar Mass Distribution separation on PSS SDV







PSS Polymer Standards Service GmbH In der Dalheimer Wiese 5 55120 Mainz | Germany
 Phone
 +49 6131 96239-0

 Fax
 +49 6131 96239-11

 E-Mail
 info@pss-polymer.com

 Web
 www.pss-polymer.com

Polymer Standards Service-USA, Inc. 160 Old Farm Rd, Suite A Amherst | MA 01002 | USA Phone +1 413 835-0265
Fax +1 413 835-0354
E-Mail pssusa@pss-polymer.com
Web www.pss-polymer.com

DE54721197

5994-6277EN July 1, 2023