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.

10271 - Column Applic



10271 - Column Application Note Characterization of Gelatin

Gelatin is a polypeptide that is produced from hydrolysis of collagenes in skin and bones. It is obtained from acidic or from basic hydrolysis and contains polyaminoacids. Gelatin is used in food, beverage and in pharmaceutical industry as stabilizer for tablets and many other applications.

Experimental Setup

Mobile Phase: Stationary Phase: Flow rate [mL/min]: Temperature [°C]: Detection: Calibration: Data processing: Phosphate buffer pH 6.6 Sodium chloride 0.3M PSS PROTEEMA 1,00 25 GPC1100 Refractive index Kit Pullulan PSS WinGPC

Recommandations for Sample Concentration

narrow PDI M 100 Da - 10 000 Da: M 10 000 Da - 1 000 000 Da: M > 1 000 000 Da: broad PDI (>1.5) all molar masses: Injection volume [µL]:

Suitable Columns

low molecular weights: medium molecular weights: high molecular weights: ultrahigh molecular weights: 2 g/L 1-2 g/L 0.5 g/L or less

3.0 - 5.0 g/L 20

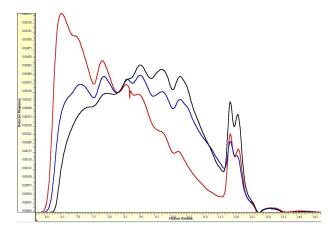


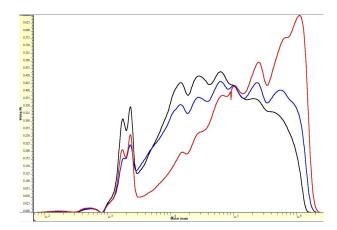
P/N pra080505 and 2x pra0830051e2 P/N pra080505 and pra0830053e2 P/N pra080505 and 2x pra0830051e3

Elugram

separation on PSS PROTEEMA

Molar Mass Distribution separation on PSS PROTEEMA







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

RA45128.3774768518

5994-6269EN August 1, 2023