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.





10009 - Column Application Note **Characterization of Cellulose Triacetate**

Cellulose triacetate is prepared from cellulose with acetic acid anhydrid in acetic acid, catalyzed by strong acids like sulfuric acid. Cellulosetriacetate is mainly used for photographic films but also in many other applications. It is soluble in dichloromethane and acetic acid. Cellulosetriacetate resists against aliphatics, but decomposes in alcohols, acetone, acetic acid and bases.

Experimental Setup

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

Shodex-RI71 Detection:

Calibration: ReadyCal-Kit Poly(styrene)

Data processing: **PSS WinGPC**

Recommandations for Sample Concentration

narrow PDI

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

Suitable Columns

low molecular weights: P/N 201-0001 (set of 3) OR sda083003lis (1 linear) P/N 201-0002 (set of 2) OR sda083005lim (1 linear) medium molecular weights: P/N 201-0003 (set of 3) OR sda083005lxl (1 linear) P/N 202-0001 (set of 3) high molecular weights:

ultrahigh molecular weights:

Elugram and Calibration separation on PSS SDV

Molar Mass Distribution separation on PSS SDV





