## **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.





## # 10307 - Column Application Note Characterization of Poly(vinyl alcohol)

Poly(vinyl alcohol) (PVOH, PVA or PVAL) is a water-soluble synthetic polymer. It has excellent film forming, emulsifying, and adhesive properties. It is also resistant to oil, grease and solvent. Unlike most vinyl polymers, PVA is not prepared by polymerization of the corresponding monomer. The monomer, vinyl alcohol, almost exclusively exists as the tautomeric form, acetaldehyde. PVA instead is prepared by partial or complete hydrolysis of polyvinyl acetate to remove acetate groups.

**Experimental Setup** 

Mobile Phase: Water Sodium chloride 0.1M Methanol (10-30%)

Stationary Phase: **PSS SUPREMA** 

Flow rate [mL/min]: 1,00 Temperature [°C]: 25

Detection: Shodex-RI71 Calibration: Kit Pullulan **PSS WinGPC** Data processing:



narrow PDI

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

20 Injection volume [µL]:

**Suitable Columns** 

low molecular weights:

precolumn, 100 Å, 100 Å precolumn, 30 Å, 1 000 Å, 1 000 Å precolumn, 100 Å, 3 000 Å, 3 000 Å medium molecular weights: high molecular weights:

ultrahigh molecular weights:

separation on PSS SUPREMA

## Overlay of different molar masses (samples with a broad molar mass distribution)



