

## Preparing the High Mass Range Calibration Solution

The high mass range calibrant is a solution of 3.5 mg/μL polypropylene glycol (PPG) in a solvent of 65:35 methanol/10 mM sodium acetate.

The high mass range calibration procedure is designed to work with a PPG that has an average molecular weight of approximately 2700 ( $M_n \sim 2700$ ), which is Aldrich product number 202347. PPG 2700 is a viscous liquid. To order this compound from Sigma-Aldrich, write or call:

Sigma Chemical Company  
P. O. Box 14508  
St. Louis, Missouri U.S. 63178-9916  
(800) 325-3010 (U.S.)  
(905) 829-9500 (Canada)  
(314) 771-3750 (outside the U.S. or Canada)  
[www.sigmaaldrich.com](http://www.sigmaaldrich.com)

## B Sample Formulations for the Velos Pro Mass Spectrometer

Preparing the High Mass Range Calibration Solution

### ❖ To prepare the sodium acetate stock solution

Dissolve 0.082 gm of sodium acetate in 10 mL of water in a clean 20 mL glass vial and label the container **Sodium Acetate Stock Solution**.

### ❖ To prepare the PPG stock solution

1. Dissolve 0.7 gm of PPG 2700 in 7 mL of methanol in a clean 20 mL glass vial.

**Tip** Because PPG 2700 is a viscous liquid, use a glass pipette to transfer 0.7 gm of the liquid into a weigh boat, or weigh the liquid directly into a minimum 20 mL glass vial.

2. Add 2.3 mL of water to the vial.
3. Add 0.7 mL of the sodium acetate stock solution and label the container **PPG 2700 Stock Solution (70 µg/µL)**.

### ❖ To prepare the calibration solution

1. Pipette 6.65 mL of methanol into a clean 20 mL glass vial.
2. Add 2.15 mL of water to the vial.
3. Add 700 µL of the sodium acetate stock solution to the vial.
4. Add 500 µL of the PPG 2700 stock solution and label the container **Velos Pro PPG 2700 Calibration Solution (3.5 µg/µL)**.