

# LSC in Practice

## Viscosity of ULTIMA-Flo Cocktails

### Problem

To help determine the proper pump for use in a laboratory automation project, a researcher contacted our laboratories and requested viscosity data for ULTIMA-Flo™ M (PerkinElmer part number 6013579) and ULTIMA-Flo AP (PerkinElmer part number 6013599) for various ratios with water.

### Discussion and Recommendation

Water was added to both ULTIMA-Flo M and ULTIMA-Flo AP at ratios up to 1:1 and the viscosities were measured at 25°C. The results are shown below:

Cocktail	No Sample	3:1 With Water	2:1 With Water	1:1 With Water
ULTIMA-Flo AP	14 cSt	22 cSt	22 cSt	19 cSt
ULTIMA-Flo M	17 cSt	28 cSt	23 cSt	23 cSt

In the above table, cSt represents centiStokes and is equivalent to mPa.s (milliPascal seconds).