

# GPC/SEC with SDV Columns and 2-Methyltetrahydrofuran

Using a mobile phase from a renewable source

## Authors

Jasmin Preis and  
Mathias Glaßner  
Agilent Technologies, Inc.

## Abstract

2-Methyltetrahydrofuran (2-methylTHF) is used as an eluent in GPC/SEC measurements with SDV columns. Reliable chromatography is displayed with polystyrene and polymethyl methacrylate standards as examples.

## Introduction

Typical eluents for GPC/SEC are nonpolar organic solvents such as tetrahydrofuran (THF) and toluene, which are based on petrochemicals. Recently, various solvents from renewable sources have become available. 2-MethylTHF is one example, which can be derived from biomass and is a potential sustainable replacement for THF.<sup>1</sup>

For an alternative solvent such as 2-methylTHF to be used as the mobile phase for GPC/SEC analysis, the following requirements need to be fulfilled: compatibility with the stationary phase is required, as well as good solubility of analytes. In addition, interaction-free chromatography and detection of the analytes with typical detectors such as refractive index (RI) detectors or UV-Vis detectors is essential.

## Experimental

**Table 1.** Instrument and sample conditions.

|                      | Conditions   |
|----------------------|--|
| Pump                 | Isocratic pump<br>Flow rate: 1 mL/min<br>Mobile phase: 2-methyltetrahydrofuran   |
| Injection System     | Autosampler<br>Injection volume: 20 $\mu$ L  |
| Columns              | SDV high MW combination:<br>SDV 5 $\mu$ m precolumn, 8 $\times$ 50 mm (p/n SDA080505)<br>SDV 5 $\mu$ m 1,000 $\text{\AA}$ , 8 $\times$ 300 mm (p/n SDA0830051e3)<br>SDV 5 $\mu$ m 100,000 $\text{\AA}$ , 8 $\times$ 300 mm (p/n SDA0830051e5)<br>SDV 5 $\mu$ m 1,000,000 $\text{\AA}$ , 8 $\times$ 300 mm (p/n SDA0830051e6) |
| Temperature          | 23 $^{\circ}$ C  |
| Sample Concentration | 1 mg/mL (0.5 mg/mL >1,000,000 Da)  |
| Calibration          | Agilent ReadyCal-Kit polystyrene (p/n PSS-PSKTR1)<br>Agilent ReadyCal-Kit polymethyl methacrylate (p/n PSS-MMKTR1)   |
| Detectors            | Refractive index (RI) detector   |
| Software             | Agilent WinGPC   |

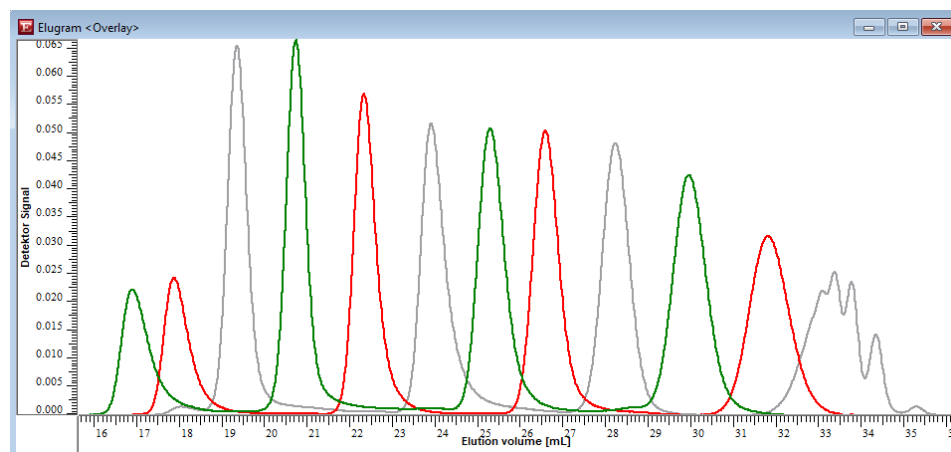
## Results and discussion

SDV columns are a typical stationary phase for nonpolar solvents, and these columns are fully compatible with 2-methylTHF. Solvent exchange between THF and 2-methylTHF is straightforward, without any loss of column efficiency when measuring plate count.

Polystyrene (PS) and polymethyl methacrylate (PMMA) reference materials show good solubility in 2-methylTHF, and GPC/SEC measurements are tested.

Using a RI detector leads to reasonable signal-to-noise ratios for both polymer types. Figure 1 depicts an overlay of three different PS mixtures with 12 standards in total covering a molecular weight range of Mp 474 to 2,520,000 Da.

The overlay of 12 different PMMA standards as three mixtures, covering a comparable molecular weight range from Mp 800 to 2,200,000 Da, is shown in Figure 2.



**Figure 1.** Overlay of three different PS mixtures (RI traces).

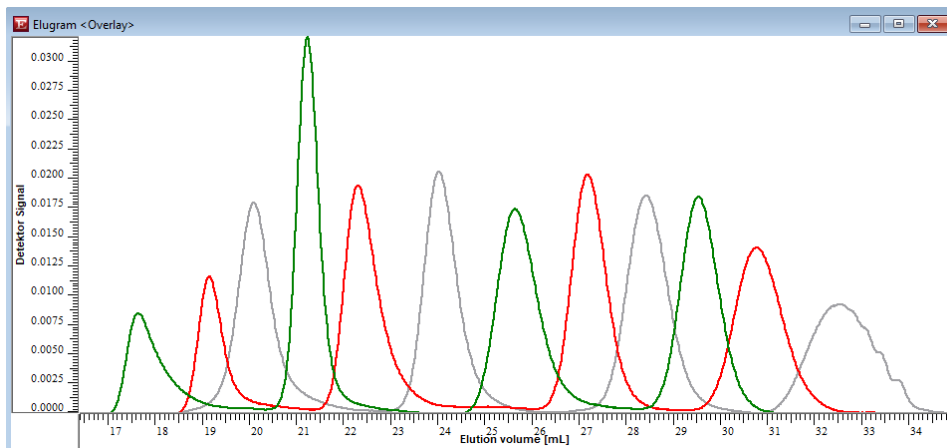


Figure 2. Overlay of three different PMMA mixtures (RI traces).

## Conclusion

2-MethylTHF, a sustainable replacement for THF, is feasible as a mobile phase in GPC/SEC for common nonpolar applications such as polystyrene or poly(meth)acrylate measurements. The use of SDV columns as stationary phase enables robust and reliable GPC/SEC measurements with RI detection.

## References

1. Pace, V. *et al.* 2-Methyltetrahydrofuran (2-MeTHF): A Biomass-Derived Solvent with Broad Application in Organic Chemistry. *Chem. Sus. Chem.* **2012**, 5(8), 1369–1379.