

# Preparative Gel Permeation Chromatography - Recycling

## Technical Overview

### Introduction

In preparative gel permeation chromatography, resolution of sample components can be improved by recycling the sample back through the column a number of times before fraction collection. This is a convenient way of minimizing cross-fraction contamination. The high column volume of the Agilent PLgel 10  $\mu\text{m}$  100  $\text{\AA}$ , 25  $\times$  300 mm prep column facilitates recycle capability (Figure 1). The improved resolution after three cycles is apparent in Figure 2.

### Conditions

Column	2 $\times$ Agilent PLgel 10 $\mu\text{m}$ 100 $\text{\AA}$ , 25 $\times$ 300 mm (p/n PL1210-6120)
Eluent	THF
Flow rate	9.0 mL/min
Detector	VWD, 254 nm
System	Agilent 1260 Infinity GPC-SEC Analysis System



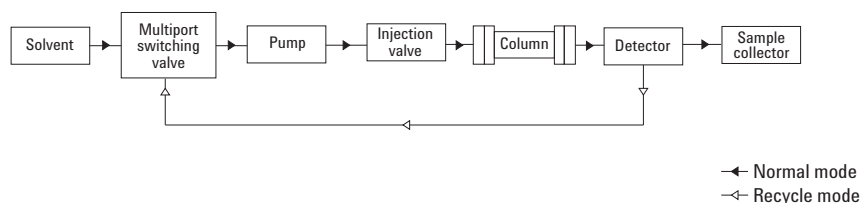


Figure 1. Schematic diagram of a preparative GPC system set up for recycling before fraction collection.

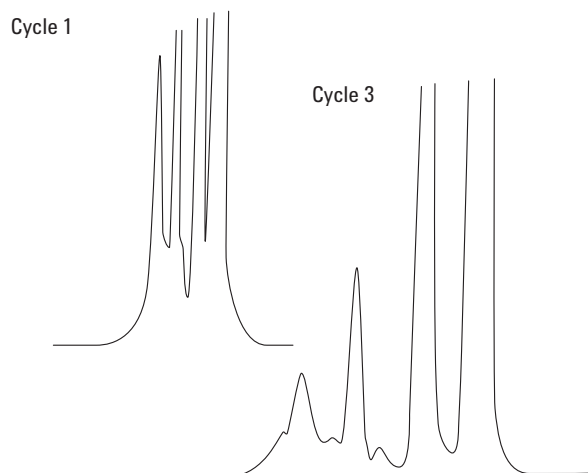


Figure 2. Improved resolution of sample components after three cycles on a set of preparative Agilent PLgel 10  $\mu\text{m}$  columns

## GPC/SEC Columns and Calibrants from Agilent

Agilent offers a comprehensive portfolio of GPC/SEC columns and calibrants for high-performance separations based on molecular size in solution. Agilent delivers leading solutions for characterizing and separating polymers by GPC/SEC, and manufactures all components for accurate polymer analysis.

Look at the Agilent Literature Library on [www.agilent.com/chem/gpc-sec](http://www.agilent.com/chem/gpc-sec) for a comprehensive range of application notes and technical overviews to help you get the best from your Agilent GPC/SEC columns and instruments.

[www.agilent.com/chem](http://www.agilent.com/chem)

Agilent shall not be liable for errors contained herein or for incidental or consequential damages in connection with the furnishing, performance, or use of this material.

Information, descriptions, and specifications in this publication are subject to change without notice.

© Agilent Technologies, Inc., 2015  
 Printed in the USA  
 April 30, 2015  
 5990-8427EN