

# Agilent PolarSil GPC/SEC Columns

## User Manual



# Notices

## Document Information

Part Number 5994-6519EN  
Published October 2023

## Copyright

© Agilent Technologies, Inc. 2023

No part of this manual may be reproduced in any form or by any means (including electronic storage and retrieval or translation into a foreign language) without prior agreement and written consent from Agilent Technologies, Inc. as governed by United States and international copyright laws.

Agilent Technologies, Inc.  
2850 Centerville Road  
Wilmington, DE 19808, USA

## Warranty

The material contained in this document is provided "as is," and is subject to being changed, without notice, in future editions. Further, to the maximum extent permitted by applicable law, Agilent disclaims all warranties, either express or implied, with regard to this manual and any information contained herein, including but not limited to the implied warranties of merchantability and fitness for a particular purpose. Agilent shall not be liable for errors or for incidental or consequential damages in connection with the furnishing, use, or performance of this document or of any information contained herein. Should Agilent and the user have a separate written agreement with warranty terms covering the material in this document that conflict with these terms, the warranty terms in the separate agreement shall control.

## Technology Licenses

The hardware and/or software described in this document are furnished under a license and may be used or copied only in accordance with the terms of such license.

## Restricted Rights Legend

U.S. Government Restricted Rights. Software and technical data rights granted to the federal government include only those rights customarily provided to end user customers. Agilent provides this customary commercial license in Software and technical data pursuant to FAR 12.211 (Technical Data) and 12.212 (Computer Software) and, for the Department of Defense, DFARS 252.227-7015 (Technical Data - Commercial Items) and DFARS 227.7202-3 (Rights in Commercial Computer Software or Computer Software Documentation).

## Safety Notices

### CAUTION

A **CAUTION** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in damage to the product or loss of important data. Do not proceed beyond a **CAUTION** notice until the indicated conditions are fully understood and met.

### WARNING

A **WARNING** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in personal injury or death. Do not proceed beyond a **WARNING** notice until the indicated conditions are fully understood and met.

# Contents

<b>1</b>	<b>Introduction</b>	<b>4</b>
1.1	Specifications	5
1.2	Solvent compatibility	5
1.3	Modifiers	6
<b>2</b>	<b>Column Installation</b>	<b>7</b>
2.1	Preparing the GPC/SEC system	7
2.2	Eluent flow rate	7
2.3	Temperature	8
2.4	Tubing	8
2.5	Connectors	8
2.6	Connecting the column to the injector or the preceding column in a set	9
2.7	Connecting columns in series	9
2.8	Connecting the column to the detector	10
2.9	Testing column performance	10
<b>3</b>	<b>Usage Guidelines</b>	<b>13</b>
3.1	Eluent preparation	13
3.2	Transferring to different solvents	13
3.2.1	Changing to eluents of similar polarity index (e.g. column in DMAc, P.I. 5.0 to 8.0)	13
3.2.2	Changing to eluent of dissimilar polarity index (e.g. column in DMAc, P.I. <5.0 or > 8.0)	14
3.3	Sample preparation	14
<b>4</b>	<b>Care and Maintenance</b>	<b>16</b>
4.1	Storage	16
4.2	Troubleshooting	16
4.3	Column repair	17
4.3.1	Partially dried-out columns	17
4.3.2	Partially blocked columns	17
4.4	Changing frits	18
<b>5</b>	<b>Ordering Information</b>	<b>19</b>

# 1

## Introduction

Thank you for choosing an Agilent PolarSil column, a high-quality, high-performance product ready for immediate use. To prolong the lifetime and to optimize its performance, please read this user manual thoroughly, paying special attention to the warnings, hints, and tips sections.

Each column has a unique serial number and is quality checked for efficiency in the eluent in which the column is delivered. The test conditions and results are shown on the certificate of analysis (CoA), which is available on our website at:

**[www.agilent.com/en/ecertificates-of-performance](http://www.agilent.com/en/ecertificates-of-performance)**

For additional information about Agilent GPC/SEC products, see

**[www.agilent.com/en/product/gpc-sec-columns-standards](http://www.agilent.com/en/product/gpc-sec-columns-standards)**

## 1.1 Specifications

PolarSil columns are packed with highly porous polar modified silica particles and are supplied as standard in Dimethylacetamide (DMAc). The columns are tested for efficiency with butylated hydroxytoluene (BHT).

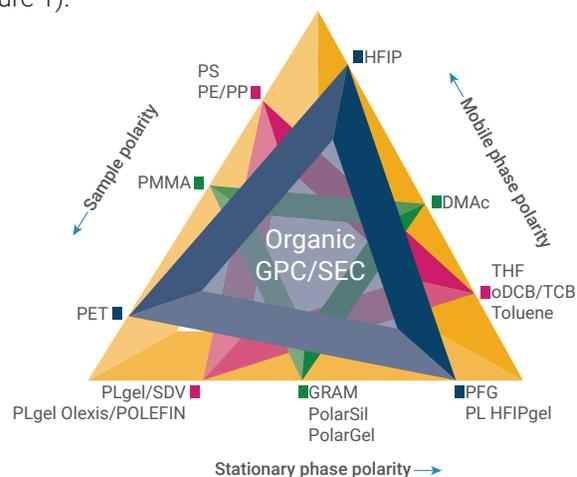
**Table 1. Operational specifications for Agilent PolarSil columns.**

Column Type	Maximum Operating Pressure per Column (bar/psi)	Maximum Operating Pressure for Three Column Set (bar/psi)	Maximum Operating Temperature	Efficiency Plates/m (BHT in DMAc)
PolarSil 3 $\mu$	100/1450	200/2901	70 °C	>50,000
PolarSil 5 $\mu$	60/870	120/1740	70 °C	>30,000

## 1.2 Solvent compatibility

PolarSil is compatible with a wide range of eluents and has optimum performance, when used with eluents that have a polarity index (P.I.) in the range 5.0 to 8.0. Transfer to compatible eluents is easy provided the guidelines in this user manual are followed.

For solvents outside this polarity range, other Agilent column materials are more appropriate to maintain a balanced polarity between, sample, eluent, and packing material for a size-only based separation based on the “magic triangle” concept (Figure 1).



**Figure 1. “Magic triangle” illustrating the fact that optimal GPC/SEC performance requires a balance between sample, stationary phase, and mobile phase polarity.**

However, it is possible to use PolarSil columns with eluents with a P.I. in the ranges 0 to 5.0, or 8.0 to 10.0. Use is restricted to neutral or acid conditions pH equal to or less than 8.0.



Certain solvents, although chemically compatible with the PolarSil packing material, are not recommended for use with PolarSil because they normally require temperatures above the recommended operating conditions as shown in Table 2.

**Table 2. The compatibility of solvents with Agilent PolarSil columns.**

Solvent	Polarity Index	Operating Temperature (°C)	Compatibility	Recommended Alternative
Acetone	5.1	25	Compatible	
Acetonitrile	5.8	25	Compatible	
Benzene	2.7	25	Possible	SDV, PLgel
Carbontetrachloride*	1.6	25	Possible	SDV, PLgel
Chloroform*	4.1	25	Possible	SDV, PLgel
Chloronaphthalene*		200	Unsuitable	POLEFIN, PLgel Olexis
Cyclohexane	0.2	25	Possible	SDV, PLgel
Decalin		150	Unsuitable	POLEFIN, PLgel Olexis
Dichloroethane*	3.5	25	Possible	SDV, PLgel
Dichloromethane*	3.1	25	Possible	SDV, PLgel
Dimethylacetamide	6.5	70	Compatible	
Dimethylformamide	6.4	70	Compatible	
Dimethylsulfoxide	7.2	70	Compatible	
Dioxane	4.8	25	Possible	SDV, PLgel
Ethanol	5.2	25	Compatible	
Ethyl Acetate	4.4	25	Possible	SDV, PLgel
Hexafluoroisopropanol*		40	Possible	PFG, PL HFIPgel
n-Methylpyrrolidone	6.7	70	Compatible	
m-Cresol		100	Unsuitable	POLEFIN, PLgel Olexis
Methanol	5.1	25	Compatible	
Methyl Ethyl Ketone	4.5	25	Possible	SDV, PLgel
o-Chlorophenol*		100	Unsuitable	POLEFIN, PLgel Olexis
o-Dichlorobenzene*	2.7	150	Unsuitable	POLEFIN, PLgel Olexis
Tetrahydrofuran	4.2	25	Possible	SDV, PLgel
Toluene	2.3	25	Possible	SDV, PLgel
Trichlorobenzene*		150	Unsuitable	POLEFIN, PLgel Olexis
Trichloroethane*		25	Possible	SDV, PLgel
Trifluoroethanol*		40	Possible	PFG, PL HFIPgel
Xylene	2.5	25	Possible	SDV, PLgel
Water	10.2	25	Possible	SUPREMA, PROTEEMA, PL aquagel-OH, AdvanceBio

\* When using chlorinated or fluorinated solvents, use PEEK/titanium frits to avoid corrosion problems. PolarSil columns are equipped with PEEK/titanium frits upon delivery.

Mixed solvent systems can be used with PolarSil columns provided the recommended Polarity Index range is not exceeded.

### 1.3 Modifiers

It is possible to use small concentrations of modifiers to improve separation performance and reduce tailing. Salts such as LiBr, LiCl, and NH<sub>4</sub>Cl in concentrations of 10 to 100 mmol/L can be added to minimize intermolecular interactions and enable proper solvation. Similarly, up to 50 mmol/L acid modifiers such as trifluoroacetic acid or acetic acid can be used to improve performance when analyzing samples with acidic groups.

## 2

# Column Installation

When using the column for the first time, it is recommended to use the eluent that the column was originally supplied in and to perform a system plate count noting the backpressure due to the column. However, this is not essential.

### 2.1 Preparing the GPC/SEC system

Remove any existing columns and store according to the manufacturer's instructions. Connect the injector directly to the detector with appropriate tubing and connectors (Section 2.5). Transfer the GPC/SEC system from the current eluent to that required, paying attention to the miscibility and compatibility with the hardware as per the manufacturer's instructions.

### 2.2 Eluent flow rate

The optimum eluent flow rates as well as the volume of the solvent contained in the column (the "column volume") depend on the column dimensions (Table 3).

**Table 3. Recommended flow rates.**

Column Dimensions	Typical Flow Rate (mL/min)	Optimum Flow Rate (mL/min)	Column Volume (mL)
30 mm x 4.6 mm	0.1 to 0.7	0.33	0.5
250 mm x 4.6 mm	0.1 to 0.7	0.33	4.2
50 mm x 8.0 mm	0.3 to 2.0	1.0	2.5
300 mm x 8.0 mm	0.3 to 2.0	1.0	15

Avoid subjecting the column to any sudden shocks, such as sudden changes in flow rate. When applying flow to the column, set the flow rate to 0.0 mL/min and turn on the pump. Always increase the flow rate in small increments, e.g., 25% of the optimum flow rate every 15 seconds until the desired operating flow rate is reached.



Even if the flow is in the allowed range, do not exceed the maximum pressure recommended for the column.

## 2.3 Temperature

When using viscous eluents, increasing the temperature decreases the eluent viscosity and therefore the column pressure and is also beneficial to the separation. In such cases, use a low flow rate initially and heat the column at a rate no greater than 2 °C/min, ensuring that the final operating temperature is at least 10 °C below the boiling point of the eluent. When at temperature, the flow can be increased to the required rate. To cool the columns down, set a low flow rate and turn off the heat to allow the columns to cool down naturally. The maximum operational temperature is 70 °C.

## 2.4 Tubing

For best results, use 1/16 in OD stainless steel tubing for connections between columns, injectors, and detectors. The ID of the tubing should be 0.007 in (0.18 mm) for microcolumns (4.6 mm ID) and 0.010 in (0.25 mm) for analytical columns (8.0 mm ID). For optimum performance, use short lengths of machine-cut capillary tubing thus ensuring that the tubing fits flush inside the fitting, minimizing dead volume and system dispersion.



Plastic tubing, such as PEEK, should be used with caution, checking the temperature and pressure ratings and compatibility with the eluents being used.

## 2.5 Connectors

Different column types use similar, but subtly different, fittings. The performance of the system can be negatively influenced, or, in the worst case, damage can occur if care is not taken. For best results, use standard stainless steel 1/16 in nuts and one-piece ferrule compression fittings.

### WARNING

The distance from the ferrule to the end of the capillary (stop depth or seating depth) must be 1 mm (Figure 2). Different column types have differing lengths of tubing protruding from the ferrule.

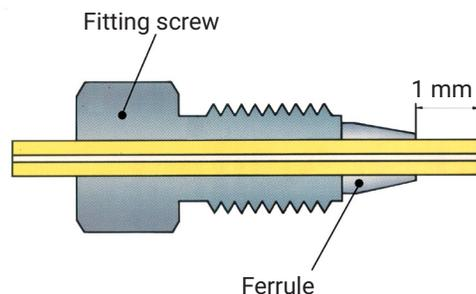


Figure 2. Schematic of the capillary connection.

Check the compatibility of the two connectors between the column and the injector and detector. For best practice, use new connectors with the Agilent column. The old connectors, together with the columns being replaced, can be put into storage if these columns will be reused.

Plastic connectors and ferrules, such as PEEK, should be used with caution, checking the temperature and pressure ratings as well as the compatibility with the eluents being used. However, these connectors offer some practical advantages in that the ferrule does not permanently grip the tubing. The connector position can be easily adjusted to accommodate the different lengths of protruding tubing used by the different manufacturers, forming a leak-tight connection with little dead volume.

## 2.6 Connecting the column to the injector or the preceding column in a set

Please use care when you install and test your new column. Remove the end plugs from the column and set aside for use when storing the column. When making the connections, place the wrenches on the fitting screw and column end fitting. Never use the flats that are machined on the column body.

Noting the direction of flow displayed on the column, connect the inlet fitting with the injector or preceding column using the appropriate nuts and new ferrules at each end of the tubing. Ensure that the capillary is as far as possible into the injector or column head before tightening the nut. Do not overtighten. The nut should be tightened enough so that the ferrule bites into the capillary tubing and no leak occurs. Excess force should not be required.

To prevent trapped air being pumped into the column, unscrew the nut at the column inlet and slowly pump a few drops of eluent to the waste before reconnecting the column.

## 2.7 Connecting columns in series

For connections between columns use the column connector delivered with each Agilent column. The precolumn should be connected to the injector and the remaining pore sizes connected in sequence of increasing pore size so that the largest porosity is closest to the detector. Only combine columns of the same particle size. To avoid porosity mismatch, we advise using the recommended column combinations shown in Table 4. Do not combine linear/mixed bed/multipore columns with single porosity columns. Each time a column is added, the procedure "Connecting the column to the injector or the preceding column in a set" should be followed.

Table 4. Recommended analytical column sets.

Separation Range [Da]	Description
100 to 40,000	1 x PolarSil precolumn, 5 $\mu\text{m}$ , 8 x 50 mm (part number PSA080505) 3 x PolarSil analytical column, 5 $\mu\text{m}$ , 100 $\text{\AA}$ , 8 x 300 mm (part number PSA0830051E2)
100 to 1,000,000	1 x PolarSil precolumn, 5 $\mu\text{m}$ , 8 x 50 mm (part number PSA080505) 3 x PolarSil analytical column, 5 $\mu\text{m}$ , linear M, 8 x 300 mm (part number PSA083005LIM)

## 2.8 Connecting the column to the detector

Connect the outlet fitting of the column to a long piece of 1/16 in tubing to take the eluent to the waste. Pump three to four column volumes of eluent through the column to the waste at the recommended operational flow rate. Check for any leaks. Stop the flow, remove the waste tubing from the outlet, and connect to the detector using the appropriate nuts and new ferrules at each end of the tubing. Apply the desired flow rate, again checking for leaks. When the detector has stabilized, the column is ready for use.

## 2.9 Testing column performance

To verify correct installation, the column can be tested by measuring the plate count/meter and the resolution and comparing with the CoA, noting that the measurements are a reflection of the dispersion in the whole system and not just the column alone. See Sections 2.2 and 3.3 for optimum flow rates, concentrations, and injection volumes.

To measure plate count, inject a suitable low molecular test probe under test conditions as detailed on the CoA.

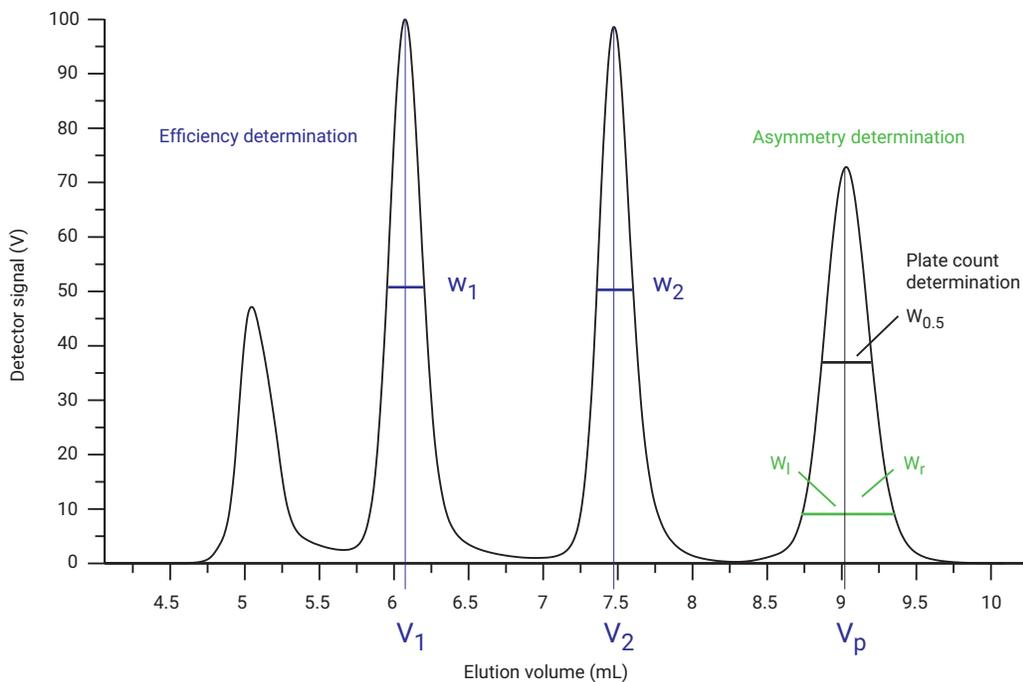


Figure 3. Variables used for determination of plate count, asymmetry, and resolution.

Determination of theoretical plates  $N$ :

$$N = \left( \frac{V_p}{\sigma} \right)^2 = 5.54 \left( \frac{V_p}{W_{0.5}} \right)^2$$

Where:

$V_p$  = Elution volume of test probe in mL

$\sigma$  = Peak dispersion

$W_{0.5}$  = Width of peak at  $\frac{1}{2}$  height in mL =  $2.35 \sigma$

$L$  = Column length in meters:

$$\frac{\text{Plates}}{\text{Meter}} = \frac{N}{L}$$

Determination of peak asymmetry  $A$ :

$$A = \frac{W_l}{W_r}$$

#### NOTE

The definition of asymmetry is as per ISO 13885-1. Sometimes, asymmetry is defined as the inverse of this formula.

A better way to determine the separation power for polymers is the calculation of the resolution factor. To do this, a mixture of polymer standards with narrow distribution of molar masses is injected:

$$\text{Resolution} = R_s = \frac{V_2 - V_1}{2(\sigma_1 + \sigma_2)} = \frac{V_2 - V_1}{0.86(w_1 + w_2)}$$

For comparison, it is better and easier to calculate the specific resolution:

$$\text{Specific resolution} = R_{sp} = \frac{0.579}{\sigma \cdot D} = \frac{R_s}{\lg\left(\frac{M_1}{M_2}\right)}$$

Where:

$M_1$ ,  $V_1$  and  $w_1$  are the molecular weight, elution volume, and peak width at ½ height of peak 1.

$M_2$ ,  $V_2$  and  $w_2$  are the molecular weight, elution volume, and peak width at ½ height of peak 2.

$$D = -\frac{d \lg M}{dV} = \text{slope of the calibration curve}$$

$\sigma$  = Peak dispersion obtained from plate count measurements

#### NOTE

When first installing a column set, it is useful to perform plate count/meter and resolution measurements as well as noting the backpressure in the chosen eluent. These parameters can be used to monitor ongoing column performance and to diagnose problems.

### 3.1 Eluent preparation

Use high-quality HPLC grade eluents, free from large particles, filtered using a 0.45 µm filter if required. Degas the eluent thoroughly, preferably using an online degasser.

#### NOTE

When using PolarSil columns with laser light scattering detectors, extra steps must be undertaken to obtain optimum low noise baselines. Filtering solvents under vacuum through a 0.2 µm filter and avoiding the use of salts and other additives in the eluent can improve performance.

### 3.2 Transferring to different solvents

Only use PolarSil columns with compatible eluents. If the column must be used with another eluent, it can be changed by careful solvent exchange without loss of resolution. Differences in plate counts that can occur in these cases are often due to viscosity differences of the eluents. Before changing eluents, ensure that both the initial eluent and the desired eluent are completely miscible. If they are not miscible, seek an intermediate that is miscible with both e.g., acetone.



Do not exceed the recommended maximum column pressure while changing solvents. Frequent transfer of columns between eluents decreases the lifetime. It is recommended, if possible, to have columns dedicated to certain eluents.

#### 3.2.1 Changing to eluents of similar polarity index (e.g., column in DMAc, P.I. 5.0 – 8.0)

1. Prepare the GPC/SEC system with the new eluent. Connect the column/column set to the injector and attach some 1/16 in tubing to the outlet directed to the waste.
2. Pump approximately two column volumes of the eluent at a 10th of the optimum flow rate through the column(s) and then slowly increase the flow rate to half the optimum flow rate and pump a further eight column volumes, e.g., for 300 mm x 8.0 mm columns, start at 0.1 mL/min and increase to 0.5 mL/min.
3. Stop flow and connect the column(s) to the detector.
4. Start flow at 0.1 mL/min. Adjust column temperature to the required value (Section 2.3) and when at temperature, increase the flow rate step by step to the desired value.
5. Ensure that all detector lines are completely purged and that the system is thoroughly equilibrated before making measurements.

### 3.2.2 Changing to eluents of dissimilar polarity index (e.g., column in DMAc, P.I. <5.0 or > 8.0)

1. Prepare a 1:1 mixture of the now used eluent with the new eluent.
2. Follow steps as described in Section 3.2.1.1 and 3.2.1.2.
3. Now prepare the GPC/SEC system with 100% of the new eluent.
4. Follow the steps in Section 3.2.1.1 to 3.2.1.5

If unsure of the polarity difference, follow the procedure in Section 3.2.2.

### 3.3 Sample preparation

Samples and polymer standards used for column calibration should be prepared in the same eluent as the GPC/SEC system. To minimize so called “solvent peaks” when using RI detection, use eluent taken directly from the solvent reservoir of the GPC/SEC system.

#### Sample injection volume and concentration

The total amount that can be injected without deterioration in separation performance depends on the total column length and is calculated from the product of injection volume and the sample concentration. A modern approach is to keep the injection volume small to minimize dispersion caused by the loop itself. If required to improve detection, vary the injected mass. The information provided in Table 5 is for guidance only. Large pressure surges related to the injection must be avoided. It may be that, for viscous solvents, viscous samples (typically very high molecular weights), and smaller particle size GPC/SEC packings, steps must be taken to reduce the pressure surge. These steps include, but are not restricted to, using lower concentrations, larger injection volumes, lower flow rates, and increasing column temperature.

Table 5. Sample quantity recommendations for optimal GPC/SEC separation performance.

Sample, Molecular Weight	Concentration (g/L)	Flow Rate	Injection Volume (μL) Column Set	
			Column Dimensions mm (L × ID)	
			250 × 4.6	300 × 8.0
Plate Count	1	Optimum	5	20
Narrow Stds. < 1 M Da	1	Optimum	5	20
Narrow Stds. > 1 M < 3 M Da	<0.5	Optimum	5	20
Narrow Stds. > 3 M Da	<0.5	50% optimum	5	20
Broad Stds., Samples < 1 M Da	1 to 3	Optimum	10	50
Broad Stds., Samples > 1 M Da	<1	50% optimum	10	50

Always ensure that samples are completely dissolved before injecting on the GPC/SEC system. High molecular weight materials require several hours for complete dissolution and are best left overnight before injection.

Sample filtration (0.5 μm or, for samples above 1 M Da, 2.0 μm) should be used after complete dissolution of the sample, to prevent blockage of the columns. However, it should be noted that filtration may potentially remove some of the sample as well as unwanted particulates.



To avoid damage of the primary column system through adsorption of sample impurities or contaminants within your samples, we strongly recommend the use of a precolumn.

## 4

# Care and Maintenance

## 4.1 Storage

If the columns are to be used again within one week, then it is not necessary to put the columns into storage provided that the eluent used does not degrade and there is no chance of salt precipitation. If feasible, the GPC/SEC system should either be put into recycle mode or be run at a very low flow rate.



Never leave the columns at elevated temperatures without flow or allow the columns to dry out. If recycle mode is used, replace with fresh eluent when the system is used again.

Replace any eluent that contains salt or acids with pure eluent before storage. All compatible eluents are suitable for storage, but the standard shipping solvent (Section 1.1) is recommended to avoid freezing at ambient temperatures. When removing the column(s) from the system, the end plugs must be replaced to prevent the column(s) drying out. If storing in volatile eluents, store in a cool place. Never let the eluent freeze.

## 4.2 Troubleshooting

Table 6. Troubleshooting guidance.

Problem	Potential Cause	Corrective Action	Recommendation
Increase in operating pressure	Blockage of solvent filter between pump and injector	Replace filter	Check pump seals and solvent quality
	Blocked or overtightened connecting tubing	Replace connector	Check sample solubility and filter samples
	Blocked inlet frit of column	Replace frit or replace column	Check solvent quality and sample solubility
	Adsorption of sample on column	Replace precolumn; undertake column regeneration	Alter analysis conditions to prevent adsorption
Decrease in operating pressure	Air in the system; leak	Check degassing of solvent; check connectors	Check if degasser is working; replace worn connectors
	Worn or badly made connections	Check and replace connectors if required	
Loss of plate count or resolution	Problem with injector	Check injector	
	Column frits blocked	Undertake column maintenance	
	Precolumn blocked/saturated	Replace precolumn	
	Analytical column damaged	Replace column	
Changing peak shape	Column damaged	Check plate counts	Replace or repair damaged column
	Sample adsorbing	Clean columns; replace precolumns	Find analysis conditions to prevent adsorption
No peak	Problem with injector	Check injector	
	Problem with detector	Check detector	
	Sample adsorbing	Clean columns; replace precolumns	Alter analysis conditions to prevent adsorption

When diagnosing problems, it is recommended to perform a series of logical experiments by removing components such as columns or connectors from the system one by one to identify which component is the cause of the problem.

### 4.3 Column repair

For best results, perform repairs on columns individually using pure eluents with no salts or acids. Always flush the columns to the waste.

The performance of a column may deteriorate due to blockage of the frits, sample adsorption, or damage to the packing bed causing an increase in backpressure.

#### 4.3.1 Partially dried-out columns

During storage, some of the solvent may evaporate through the threads (this typically happens during long storage times and with high storage temperatures and volatile eluents). It is good practice to keep columns with volatile eluents in a refrigerator when not in use, to prevent solvent evaporation. It is an indication of a partially evaporated column solvent if the expected pressure does not build up or the pump constantly readjusts the flow. In this case, immediately reduce the flow down to 0.1 mL/min (to prevent damaging the packing) and fill the column until no more bubbles appear at the column outlet and column backpressure is stable. Now add the next column using the previously mentioned installation procedure for installation until each column is installed in your system.

#### NOTE

If you want to install a complete set of columns whose solvent has been evaporated partially, you can connect all columns when solvent appears from the previous column. Then it can be pumped at 0.1 mL/min, unattended (for example overnight), with no negative effects.

#### 4.3.2 Partially blocked columns

Columns become partially blocked when particulates get trapped in the frits or in the packing bed and/or sample components adsorb on the frits or packing material. Due to the varying chemical nature of possible samples and column history, there is no single treatment that is successful in all cases. It is often not possible to recover a column. Certain steps can be taken, either individually or combined to try to partially recover performance. These steps should be undertaken on individual columns to avoid potential contamination problems.

Reverse flush the column to the waste at a low flow rate (25% of the optimum flow rate) for at least four column volumes. Afterwards, reconnect the column in the correct direction. Slowly increase the flow rate to see if the backpressure generated on the column has reduced and if it can be operated under normal conditions. If the pressure is now reasonable, perform a plate count test and also, if desired, a resolution test. If the pressure remains high, either try the following steps or directly change the frits.

1. Increase the temperature of the column in combination with the procedure in Section 4.3.2 following the guidelines in Section 2.3. Flush for 10 column volumes.
2. Change the eluent and/or use modifiers to try to dissolve any adsorbed components in combination with the procedure in Section 4.3.2 and, if appropriate, step 1 in this section. Flush for 10 column volumes.

#### 4.4 Changing frits

This procedure can be used for changing both the inlet and outlet column frit. The inlet frit is particularly susceptible to blocking. The column should only be opened to change frits if a blockage could not be removed by the above measures.



**Figure 4. Components of the column head**  
**A) End fittings**  
**B) Fitting adapters**  
**C) Frits**

1. Remove the column from the system and replace the end plug in the end that is not being changed.
2. Using two wrenches, one placed on the flat of the column body nearest the end fitting (Figure 4) to be changed and the end fitting itself, slowly unscrew and remove the end fitting, the end fitting adapter, and finally slide off the frit and seal from the top of the column.
3. Inspect the top of the gel packing to ensure that there are no holes in the gel packing.
4. Inspect the end fitting adapter to ensure it is not damaged.
5. Rinse end fitting and fitting adapter to ensure that they are entirely clean and free of particles. Re-assemble the fitting adapter (flat end towards the frit) and a new frit in the end fitting.
6. Holding them at an angle of 45° to each other, screw the end fitting, the adapter, and new frit back onto the column until finger tight. Then use the wrenches to tighten by a further 60° turn.
7. Connect the column in reverse direction and flush for three column volumes to the waste, ensuring that gel particles do not elute due to a poorly seated frit and seal.
8. Connect the column in the correct direction and check the pressure and plate count.

## 5. Ordering Information

Table 7. Ordering information for spare parts.

Part Number	Description
299-2045	PEEK/titanium replacement frits for 8.0 mm ID column (pk of 2)
299-2046	PEEK/titanium replacement frits for 4.6 mm ID column (pk of 2)
PL1310-0008	Tubing ferrules, 1/16 in, 5/pk
PL1310-0007	Column connecting nuts, 1/16 in, 5/pk
PL1310-0048	Connecting tubing for inter-column connection, 100 mm length, ID 0.25 mm, 10/pk, for 8.0 mm and 20 mm ID columns, combine with ferrules (PL1310-0008) and nuts (PL1310-0007)
5021-1816	Connecting tubing for inter-column connection, 105 mm length, ID 0.17 mm, for 4.6 mm ID columns, combine with ferrules (PL1310-0008) and nuts (PL1310-0007)
5021-1818	Connector from column to injector/detector, 280 mm length, ID 0.17 mm, for 4.6 mm ID columns, combine with ferrules (PL1310-0008) and nuts (PL1310-0007)
5022-6508	Connector from column to injector/detector, 280 mm length, ID 0.25 mm, for 8.0 mm ID columns, combine with ferrules (PL1310-0008) and nuts (PL1310-0007)

Contact us:

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

Buy online:

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

Get social with Agilent:

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

Get answers to your technical questions and  
access resources in the Agilent Community:

**[community.agilent.com](http://community.agilent.com)**

DE07786567

This information is subject to change without notice.

© Agilent Technologies, Inc. 2023

Published in the USA, October 1, 2023

5994-6519EN

