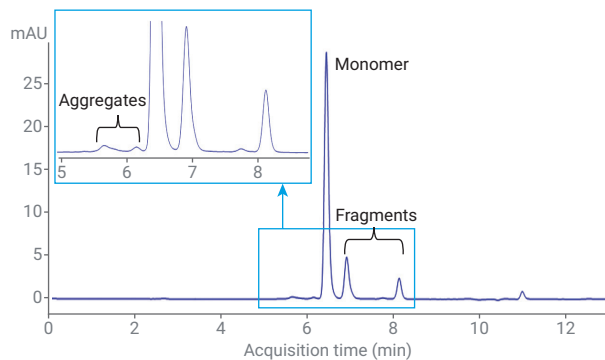


- Prepare your samples in the mobile phase, and ensure that they dissolve completely. Filter or centrifuge samples before injection.

**Note:** To maximize the lifetime of your column, we recommend using an Agilent Advance Bio SEC guard column.

#### Recommended starting conditions

| Parameter          | Value   |
|--------------------|---|
| Column             | AdvanceBio SEC 200 Å, 4.6 × 300 mm, 1.9 μm (p/n PL1580-5201)  |
| Instrument         | Agilent 1260 Infinity II Bio-inert LC System  |
| Flow Rate          | 0.35 mL/min   |
| Mobile Phase       | 150 mM sodium phosphate, pH 7.0   |
| Wavelength         | 280 nm  |
| Column Temperature | 25 °C   |
| Sample             | A stressed mAb (1 μg, injected onto column)<br>mAb sample stressed in 100 mM sodium bicarbonate pH 9 and incubated overnight at 40 °C |



## Column care and cleaning

### Column care

An increase in backpressure and decrease in performance may occur over time. If the pressure has increased, first identify if this increase is due to a guard column that may need replacing. If the increase in pressure is in a system component, such as tubing or a filter, replace the component and retest.

### Column cleaning instructions

It may be possible to restore performance of the columns using one of the following cleaning solutions:

- **For strongly adsorbed contaminants:** high salt concentration at low pH (for example, 0.5 M Na<sub>2</sub>SO<sub>4</sub>, pH 3) or 0.5 M guanidine hydrochloride
- **Organic solvent for hydrophobic materials:** up to 50% methanol, ethanol, or isopropanol
- **Acidic reagents for basic contaminants:** 0.1% TFA, formic acid, or acetic acid in 15% acetonitrile

Always flush the column in the direction of the flow arrow, and adjust the flow rate to keep the pressure below 400 bar. Rinse with at least five column volumes of ultrapure water before and after flushing with at least 20 column volumes of the cleaning solution.

**It is not recommended to use all three cleaning buffers sequentially. Choose the most appropriate buffer for your probable contaminant. Take care to avoid precipitation of buffer salts, and avoid overpressuring the column due to mobile phase viscosity differences.**

### Recommended storage

- **Shorter-term storage (less than two weeks):** store the column in the mobile phase.
- **Extended storage (longer than two weeks):** store in 100 mM sodium phosphate, pH 6.7 with 0.02% NaN<sub>3</sub> or 20% methanol in water for a minimum of 10 column volumes. Lower flow rates (0.1 mL/min) are recommended for 20% methanol flush to avoid over-pressuring the column due to high viscosity.
- Storage temperature is room temperature.

### Ordering details

| Description                                   | Part Number   |
|---|---------------|
| AdvanceBio SEC 120Å 1.9 μm, 4.6 × 300 mm      | PL1580-5250   |
| AdvanceBio SEC 120Å 1.9 μm, 4.6 × 150 mm      | PL1580-3250   |
| AdvanceBio SEC 120Å 1.9 μm guard, 4.6 × 30 mm | PL1580-1250   |
| AdvanceBio SEC 200Å 1.9 μm, 4.6 × 300 mm      | PL1580-5201   |
| AdvanceBio SEC 200Å 1.9 μm, 4.6 × 150 mm      | PL1580-3201   |
| AdvanceBio SEC 200Å 1.9 μm guard, 4.6 × 30 mm | PL1580-1201   |
| AdvanceBio SEC 120Å 1.9 μm, 2.1 × 150 mm PEEK | PL1980-3250PK |
| AdvanceBio SEC 120Å 1.9 μm, 2.1 × 50 mm PEEK  | PL1980-1250PK |
| AdvanceBio SEC 200Å 1.9 μm, 2.1 × 150 mm PEEK | PL1980-3201PK |
| AdvanceBio SEC 200Å 1.9 μm, 2.1 × 50 mm PEEK  | PL1980-1201PK |

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

For Research Use Only. Not for use in diagnostic procedures.

DE.2752893519

This information is subject to change without notice.

© Agilent Technologies, Inc. 2019, 2020

Printed in the USA, July 1, 2020

5994-0739EN

## Agilent AdvanceBio SEC 1.9 μm Column User guide

Agilent size exclusion columns optimized for separation and characterization of size variants are offered in two pore sizes: Advance Bio SEC 120Å is best suited for analysis of peptides and small therapeutic proteins; AdvanceBio SEC 200Å was designed for characterizing mAbs and ADCs. In addition to 4.6 mm id stainless steel columns, AdvanceBio SEC 1.9 μm is also available in PEEK-lined stainless steel columns in 2.1 mm id. These bioinert columns provide a metal-free flow path and are particularly suited to SEC-MS applications.



## Getting started

A column performance report, including a column-specific quality control (QC) test chromatogram and a batch-specific protein separation, is enclosed with every Agilent AdvanceBio SEC 200 Å 1.9 µm column. The QC test system has been modified from a standard system to minimize dead volume, so it may vary from the system used in your lab. This modification enables better evaluation of the column efficiency, and ensures a more consistent product. An optimized LC system will generate similar results to the chromatogram in the column performance report. For the best chromatographic results, it is recommended that a low dispersion LC is used. You can optimize your LC for maximum resolution by minimizing tubing internal diameter (id) and length between the sample injector and the column and between the column and detector. Low volume micro UV detector flow cells may be necessary with 2.1 mm id columns.

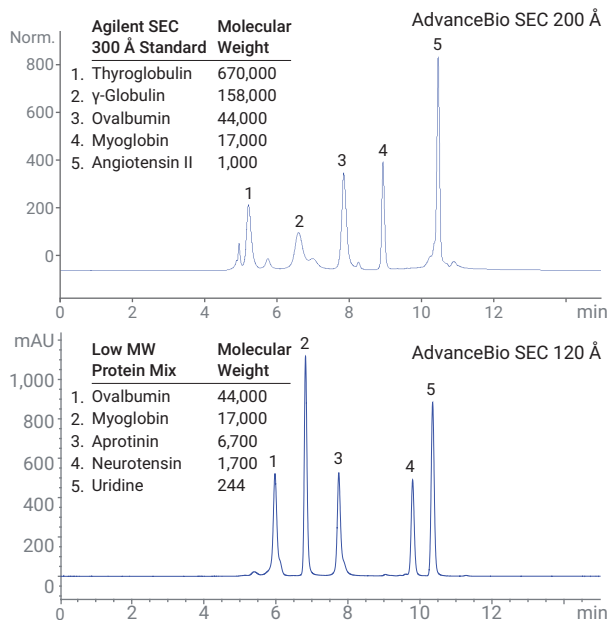
Ensuring proper column connection is important. Agilent recommends InfinityLab Quick Connect LC fittings (p/n 5067-5966), or Bio-inert UHP-FF fittings (p/n 5067-5695), particularly for use with PEEK-lined 2.1 mm id columns to avoid damage.

If you have specific questions, contact Agilent Technical Support at [www.agilent.com/chem/techsupport](http://www.agilent.com/chem/techsupport).

To monitor column and instrument performance, Agilent recommends running a standard test mixture regularly, such as the AdvanceBio SEC standards.

The AdvanceBio SEC 200 Å 1.9 µm columns are recommended for use with UV, DAD, LS detectors, and for SEC-MS applications under denaturing and native mode.

| Parameter          | Value  |
|--------------------|--|
| Columns            | AdvanceBio SEC 200Å, 4.6 × 300 mm, 1.9 µm (PL1580-5201)<br>AdvanceBio SEC 120Å, 4.6 × 300 mm, 1.9 µm (PL1580-5250)   |
| Flow Rate          | 0.35 mL/min  |
| Mobile Phase       | 150 mM sodium phosphate, pH 7.0  |
| Wavelength         | 220 nm   |
| Column Temperature | Ambient  |
| Samples            | AdvanceBio SEC 300 Å Protein Standard (p/n 5190-9417);<br>Low MW Protein Mix (Ovalbumin 5.0 mg,<br>Myoglobin 5.1 mg, Aprotinin 5.1 mg, Neurotensin 2.3 mg,<br>and Uridine 2.4 mg) in 1 mL mobile phase |



## Important safety considerations

- All points of connection in an LC system are potential sources of leaks. Users of liquid chromatographic instruments should be aware of the potential toxicity or flammability of their mobile phases.
- Do not remove the column end fitting.

## Using your column

### Column specifications

| Parameter                                   | Value  |
|---|--|
| Shipping Solvent/Long Term Storage Solution | pH 6.7 100 mM sodium phosphate buffer with 0.02 % NaN <sub>3</sub>   |
| Working Flow Rate                           | for 4.6 × 150 mm 0.1 to 0.7 mL/min<br>for 4.6 × 300 mm 0.1 to 0.5 mL/min<br>for 2.1 mm id columns 0.05 to 0.10 mL/min  |
| Maximum Pressure                            | 620 bar (9,000 psi)  |
| pH Stability                                | 2 to 8   |
| Salt Concentration                          | ≤0.5 M   |
| Mobile Phase Compatibility                  | Compatible with all the SEC mobile phases for UV. Phosphate buffer, pH 7.0 with different salt concentrations and denaturing and native mode SEC-MS mobile phases. |
| Operating Temperature                       | 20 to 40 °C (recommended),<br>80 °C (maximum)  |

**Note:** Working at extremes of the operating parameters may reduce column lifetime.

### Installation

- Remove both end plugs, and ensure that your system's flow direction matches the arrow on the column. Do not use the column with the flow in the reverse direction.
- Use an Agilent InfinityLab Quick Connect LC fitting (p/n 5067-5966) to quickly connect the column to your LC

instrument, or Bio-inert UHP-FF fittings (p/n 5067-5695) for use with PEEK-lined 2.1 mm ID columns.

### Column conditioning

- The columns are shipped in 100 mM sodium phosphate buffer, pH 6.7, containing 0.02% NaN<sub>3</sub>, and must first be flushed into the mobile phase required for your separation. Ramp the flow rate slowly from 0.0 mL/min up to the intended operating flow rate over a period of several minutes. If possible, the maximum flow gradient should be set at 0.1 mL/min/min. Equilibrate the column by flushing for a minimum of 10 column volumes, or until the baseline is stable.

### Instructions for use

- Columns are compatible with commonly used aqueous buffers, including 150 mM sodium phosphate at pH 7.0 with or without the addition of other salts. Salt concentration should not be more than 0.5 M. It is recommended that the percentage of organic solvent be less than 50%. For native mode SEC-MS, ammonium acetate is recommended. Flush the column extensively before connecting to your MS detector in case nonvolatile mobile phase salts are still present. When changing eluents, always take into account the viscosity and risk of salt precipitation. If you are unsure, flush the column first with high-purity water before introducing a new eluent.
- Mix your buffers freshly using high-purity components and ultrahigh purity water such as Milli-Q or Nanopure. Filter buffers through a 0.2 or 0.45 µm filter, and degas before use. Filtering will remove particulates, and help reduce the risk of bacterial growth, which will otherwise damage the column and your LC system.