

Dear Agilent Customer,

Thank you for purchasing Agilent products.

To reduce our environmental footprint, we have moved our ZORBAX Reliance Cartridge Guard Columns data sheet online:

www.agilent.com/chem/reliance-guards-user-guide

For more information

For more information on our products, visit
www.agilent.com

For technical support in the US and Canada, call
1-800-227-9770, or call your local Agilent sales office.

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.

DE.0458680556

© Agilent Technologies, Inc., 2020
Printed in the USA, April 2020
p/n 830147-002



[Go to data sheet](http://www.agilent.com/chem/reliance-guards-user-guide)

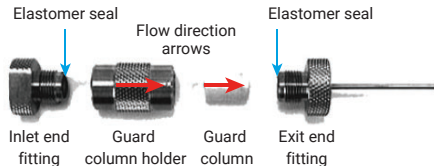


Agilent

| Trusted Answers

Agilent ZORBAX Cartridge Guard Columns

Assembly guidelines



First, loosely screw the inlet end fitting into the inlet end of the guard column holder, leaving one or two threads showing. Place the assembly on the benchtop, open end up. Put the polymeric guard cartridge into the holder with the arrow on the cartridge pointing up. With the guard column holder still vertical on the benchtop, screw the guard column exit end fitting into it until it is finger tight. Hand tighten the full assembly by gripping both ends and turning them clockwise until tight. To connect the guard column assembly to the analytical column, use the integral column connector provided (0.010 inch id $\times$ 30 mm length).

Alternatively, you can use a different inlet end fitting (included) on the exit end of the guard column holder. This allows you to use other connecting tubing, if desired.

Operational guidelines

- Do not use a mobile-phase pH outside of the range of the analytical column being used.
- Only finger-tighten the cartridge holder to avoid damaging the polymeric cartridge.
- HPLC guards may be used for temperatures up to 80 °C. Do not exceed the maximum operating temperature of the analytical column.
- Routinely check for the presence and integrity of the perfluoro-elastomer seals on the end fittings before every assembly operation. Replace when necessary.
- Minimize the system dead volume to maintain efficiency.
- All cartridges are shipped dry. Before use, flush with methanol to wet packing, then flush with mobile phase to ensure the highest column performance. Use the minimum volumes:

Cartridge id (mm)	MeOH Wetting Vol. (mL)	Mobile Phase Flush (mL)
2.1	0.5	2
3	2	6
4.6	3	10
9.4	10	20