

Important notice on fittings for InfinityLab Poroshell and AdvanceBio PEEK-lined columns:



Dear Agilent Customer,

Thank you for your purchase of a PEEK-lined Agilent column.

Please note that care must be taken to avoid damage to PEEK-lined columns during installation. Combined compression and rotation may cause internal damage, so fittings without ferrules (such as PEEK finger-tight fittings) are not recommended.

Agilent offers fittings for bio-inert and stainless steel systems.

To choose the best fitting and capillary for your bio-inert instrument setup, please visit www.agilent.com/chem/bioinertfittings

To choose the best fitting and capillary for your stainless steel system, please visit www.agilent.com/chem/fittings

Less stress. More reliable connections.

With Agilent InfinityLab fittings, you can be sure that you're getting a perfect column connection, every time.



InfinityLab Quick Connect fittings are truly "finger-tight," reusable fittings for UHPLC, leak-free to 1300 bar. You don't need any tools. Simply use your fingers to close the lever for a perfect connection every time. This fitting is perfect for the column inlet.



With **InfinityLab Quick Turn fittings**, you will get either a finger-tight connection (leak-free to 400 bar), or a UHPLC connection (leak-free to 800 bar with mounting tool p/n 5043-0915, and 1300 bar after a quarter turn of a wrench). The spring-loaded design guarantees zero-dead-volume, and makes it ideal for connections at the column outlet and detector.

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.0760069444

© Agilent Technologies, Inc.,
2020

Printed in the USA, April 2020
p/n 830147-003