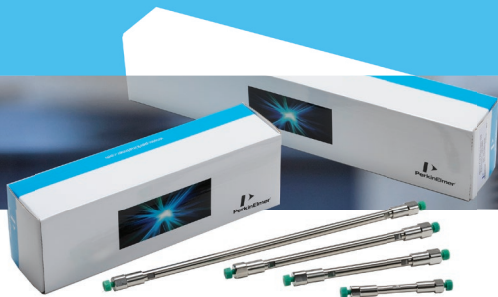


SEPARATION CHALLENGE ACCEPTED



Quasar™ Liquid Chromatography Columns



QUASAR LC COLUMNS FOR YOUR NEW LC 300 SYSTEM

Whatever your separation challenge, your choice of liquid chromatography (LC) column can make all the difference. Our Quasar portfolio of LC columns combined with the improved workflow and usability of the LC 300 system allows you to achieve rugged and reproducible results.

Ultrapure silica-based Quasar columns deliver a comprehensive range of chemistries, together with state-of-the-art, optimized bonding technology to give you a versatile, high-performing analytical solution for your increasingly complex samples.

Quasar HPLC and UHPLC Columns

- High sensitivity for trace analyte identification
- High efficiency for complex separations
- Excellent pH stability across commonly used mobile phase buffers

Quasar SPP Columns

- Next-generation superficially porous particle (SPP) phases that promise productivity
- Faster run times and method development
- Offers robust operation with lower back pressures when compared to sub 2 μm columns

Our Quasar columns offers benefits such as:

- Longevity to withstand high-throughput environments
- High-efficiency separations to facilitate trace-level analysis
- Good resolution and peak shapes for more effective compound separation
- Analysis of increasing polar and complex analytes

LC 300 HPLC and UHPLC Systems

- Robust and precise autosampler with integrated column oven and preheater
- Choice of 10K or 18K psi pump and autosampler
- Portfolio of five available detectors (PDA, MWD, UV/Vis, RI and FL), which can be used individually or in sequence
- Efficient and powerful instrument control with Simplicity™ Chrom CDS

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com

For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2020, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

NEW INSTRUMENT OFFER

FREE OF CHARGE

Quasar LC column with
your new LC 300 system

Please contact
your sales representative
for details.

* Terms and Conditions: No other PerkinElmer products qualify. This offer cannot be applied to prior orders or purchases. Limit one offer per instrument. Other restrictions may apply. Void where prohibited.

For a complete listing of our Quasar columns portfolio or more details, please call your local PerkinElmer representative or visit www.perkinelmer.com/quasar

