

# SEPARATION CHALLENGE ACCEPTED

## Quasar Liquid Chromatography Columns

Whatever your separation challenge, your choice of liquid chromatography (LC) column can make all the difference. Our Quasar™ portfolio of LC columns allows you to achieve rugged and reproducible results – batch to batch and column to column – with an all-encompassing, flexible solution that meets the diverse, changing needs of analysis.

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.

For flexibility, we provide a wide range of column sizes, including shorter columns packed with smaller particle sizes for shorter run times and better productivity. Plus, our scalable columns facilitate easy method transfer between HPLC and UHPLC technology platforms – and the smaller particle sizes means optimized sensitivity for those applications. Whatever your separation need, we have a chemistry or dimension to fill it.

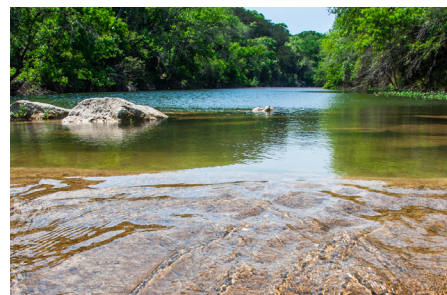
### Our wide range of column sizes has its benefits, including:

- 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



# Choose the Columns that are Right For You

With Quasar LC columns, you now have the tools to confidently say: **separation challenge accepted.**



## Quasar HPLC and UHPLC Columns

- High sensitivity for mass spectrometry (MS) applications
- High efficiency for complex separations
- Supports both high- and low-throughput environments
- Increases productivity and reduce run times
- Excellent pH stability across commonly used mobile phase buffers
- High-sample loading capacity

## Quasar SPP Columns

- Next-generation superficially porous particle (SPP) phases that promise productivity
- Faster run times and method development
- No specialized filtrations of sample and mobile phase
- Optimized low-band spreading
- Lower back pressures compared to sub 2 $\mu$ m columns with robust operation

For more information about Quasar Columns, visit [www.perkinelmer.com/quasarlc](http://www.perkinelmer.com/quasarlc)

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



For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright © 2019, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

014465\_05

PKI