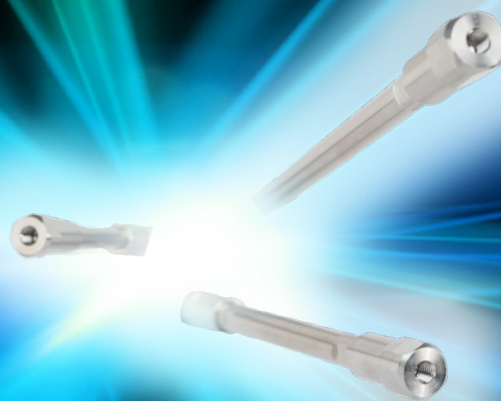


# WEBCAST: CHOOSING THE RIGHT C18 REVERSED PHASE HPLC COLUMNS



## Join Our Live Webcast

Whether you're testing food, water, or pharmaceuticals, success often depends on getting back to basics. The most commonly used LC reversed-phase alkyl bonded stationary phase is octadecyl carbon chain (C18)-bonded silica, which is denoted as USP classification L1. However, not all C18 columns are alike.

This webinar will provide details of a comparative study between twenty-four silica-based C18 phases and how C18 columns can be compared in terms of:

- Relative hydrophobicity
- Column efficiencies
- Silanol activity



C18 columns

Please join us and Kathryn Lawson-Wood to discuss the importance of choosing the right LC column for your analysis and how differences from one C18 column to another can result in differences in retention time, resolution and even selectivity.

### LIVE WEBCAST DETAILS:

**Date:** Friday, May 29, 2020

**Time:** 8:00 AM ET | 14:00 CET

**Duration:** 90 minutes



**Speaker:**

Kathryn Lawson-Wood,  
Consumables Product Specialist,  
PerkinElmer, Seer Green, UK

SEE DETAILS AND REGISTER NOW [www.perkinelmer.com/HPLCColumnsWebinar](http://www.perkinelmer.com/HPLCColumnsWebinar)

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