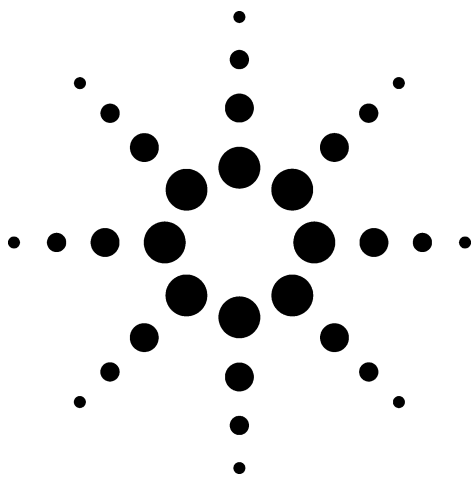




High-Speed Separation of Parabens

Application Food Analysis Robert Ricker

Parabens, or para-hydroxy benzoic acid alkyl esters are popular preservatives used in the cosmetic and food industry to battle microbe degradation. Analysis time of less than one minute is accomplished with a Rapid Resolution SB-C18, 4.6 x 30 mm column. Columns with particle sizes under 5 μm and dimensions up to ten-times smaller than traditional analytical-size columns are ideal for high-speed methods. If desired, increasing temperature cuts analysis time further.

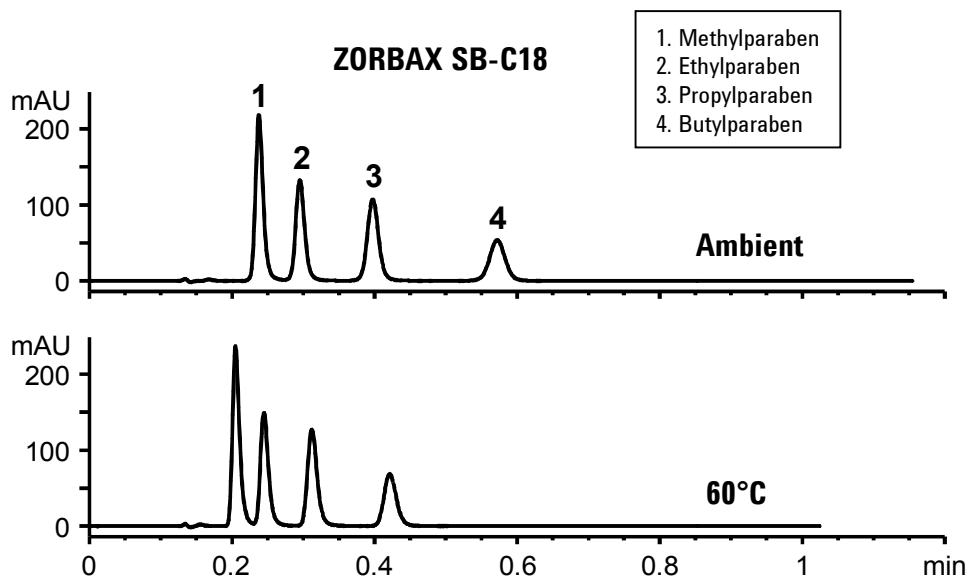
Extra-column volume is a crucial factor in chromatographic performance. Here, no modification to the modern instrument is necessary for optimal resolution.

Operating Conditions:

HPLC System:	Agilent 1100 with quaternary pump
Column:	ZORBAX StableBond-C18 Rapid-Resolution (3.5 μm) Cartridge-Column, 4.6 x 30 mm
Mobile Phase:	Agilent Part No. 833975-902 0.1% H_3PO_4 : ACN, (50:50)
Detection:	UV 254 nm with standard flow cell (13 μL)
Flow:	2 mL/ min.
Inj. Volume:	1 μL
Temperature:	top: ambient, bottom: 60°C

Highlights

- Reducing column length and particle size simultaneously can:
 - Reduce analysis time
 - Maintain resolution
 - Reduce solvent use
- Elevated operating temperature is effective in reducing run time.
- ZORBAX StableBond SB-C18 can operate at higher temperatures and lower pH than other commercial reversed-phase columns.



Agilent Technologies

*Robert Ricker is an application chemist
based at Agilent Technologies, Wilmington,
Delaware.*

For more information on our products and
services, visit our website at:
www.agilent.com/chem

Copyright© 2002 Agilent Technologies, Inc.
All Rights Reserved. Reproduction,
adaptation or translation without prior
written permission is prohibited, except as
allowed under the copyright laws.

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.

Printed in the USA
April 25, 2002
5988-6356EN



Agilent Technologies