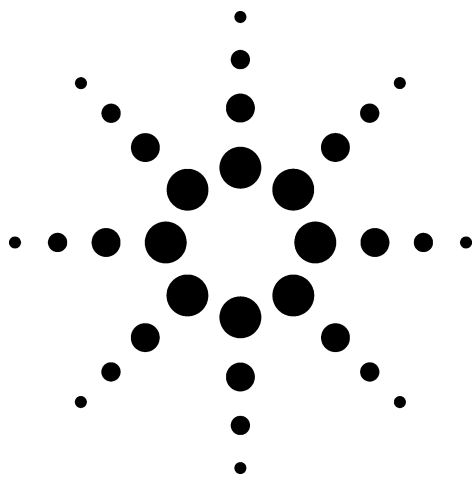


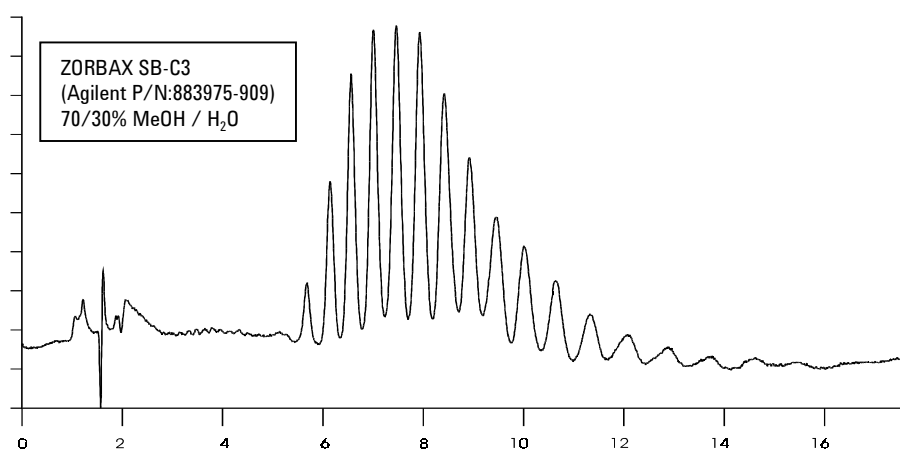


Decreasing Run Time by Changing Bonded Phase: Triton X-114

Application

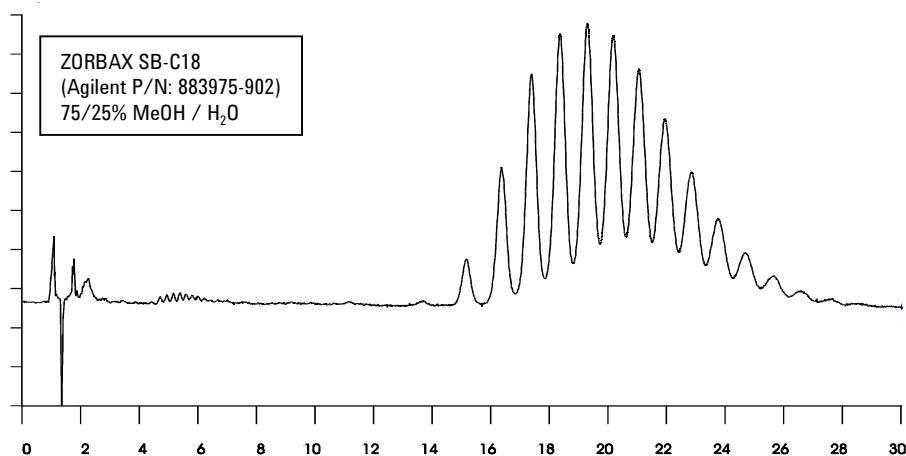
Polymers

Robert Ricker



Highlights

- Analysis time reduced to one-half without loss in resolution.
- Chromatograms show the advantage of having bonded phases with different retentivities.



Conditions:
ZORBAX SB-C3, SB-C18 (4.6 x 150 mm) Agilent P/N: (883975-909), (883975-902)
Mobile Phase: MeOH and H₂O (as indicated)
1 mL/min, 50°C, Detect. UV(225 nm)



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
5988-6377EN



Agilent Technologies