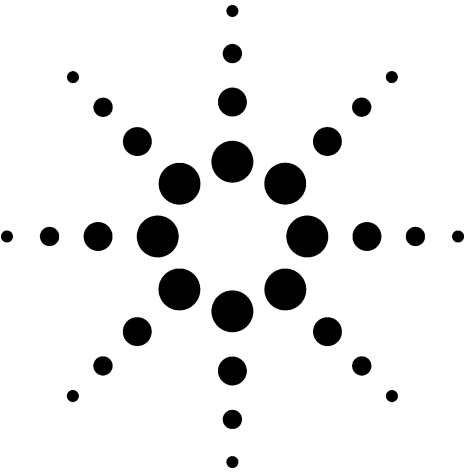




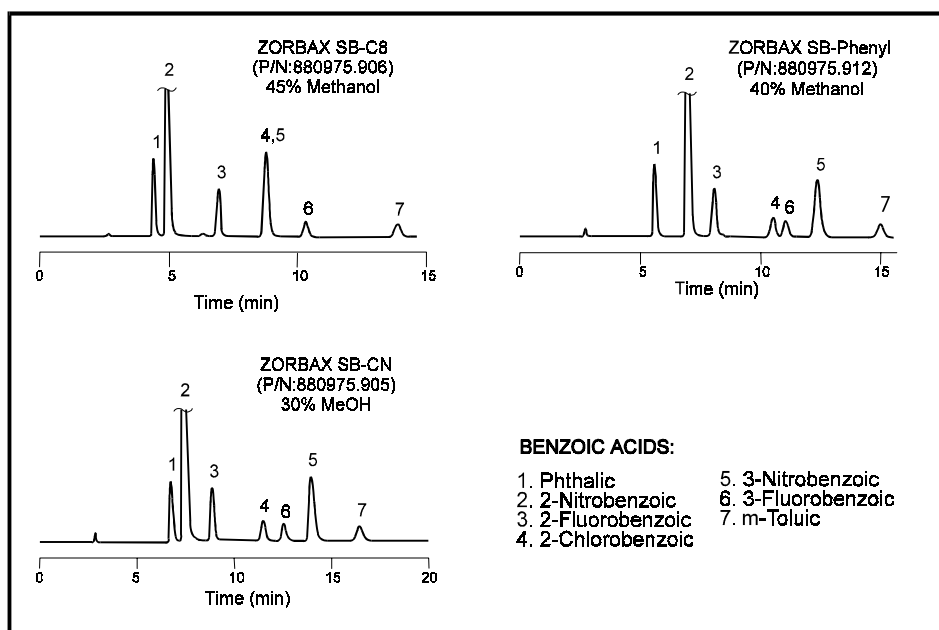
Benzoic Acids

Examples of Bonded-Phase Selectivity Differences

Application
Organic Chemicals
Robert Ricker

Highlights

- Significant selectivity differences are obtained by changing columns rather than mobile phase.
- Allows use of same series of high stability columns, rather than switching to columns from other manufacturers.
- For this sample mixture of Benzoic Acids, ZORBAX SB-CN provides the best overall resolution (minimum resolution is $R_s = 1.90$ for peak pair 4 and 6).



Conditions:
ZORBAX StableBond Columns (4.6 x 250 mm) (Agilent P/N: above)
Mobile Phase: 30-45% methanol (above) in 25mM Na Phosphate, pH 2.5
1.0 mL/min; 35°C



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
5988-6368EN



Agilent Technologies