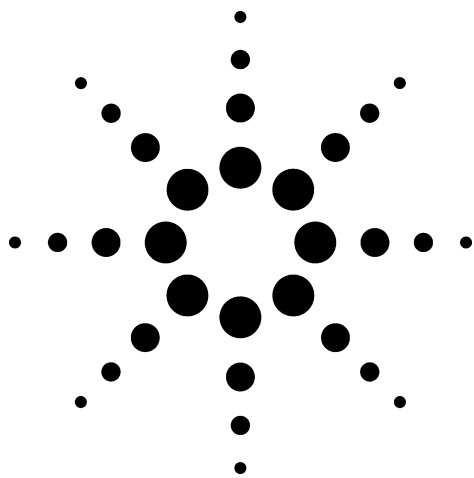
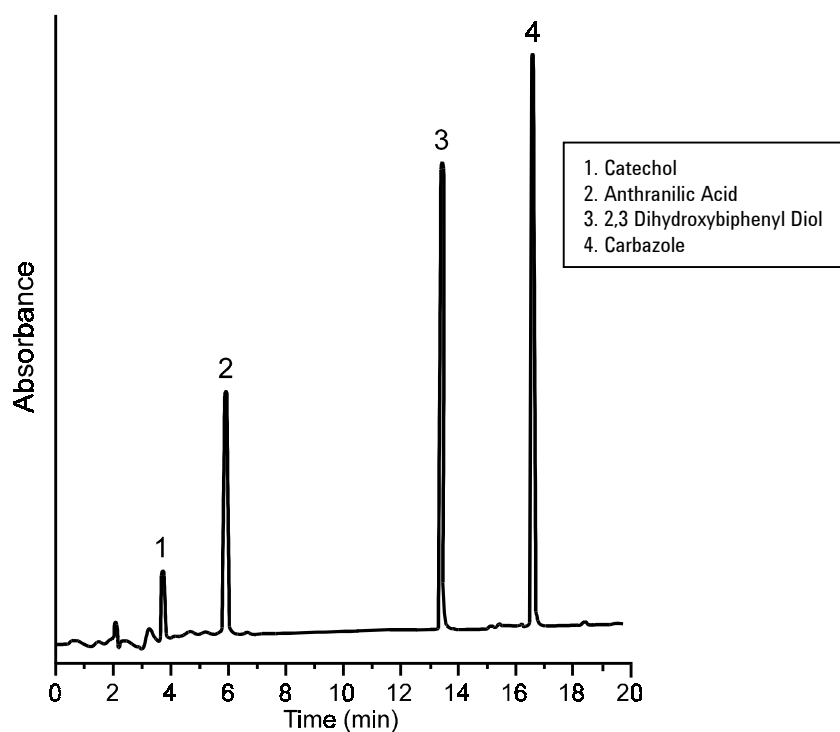




Separation of Catechol and Various Metabolites

Application
Environmental
Robert Ricker

Carbazole is an environmental pollutant of some concern. Sensitive detection of carbazole and related metabolites provides a means of following its breakdown in various sources. The postulated pathway is: **carbazole** → **2,3-aminobiphenyl diol** → **anthranilic acid** → **catechol**. (The 2nd intermediate is not commercially available and was replaced with 2,3-dihydroxybiphenyl diol, to act as a related marker.)



Courtesy of Phillip Gibbs, Energy Biosystems Corp. The Woodlands, TX;
Rich Willson, University of Houston, Dept. Chemical Engineering, Houston, TX

Conditions:
ZORBAX 300 SB-C18 (4.6 x 150 mm) (Agilent P/N: 883995-902)
Mobile Phase: Gradient 5-60% in 20 min., Wash, 95%B for 2 min.
A) H₂O with 0.1% TFA to pH3 with TEA; **B)** ACN with 0.085% TFA
Injection 15µL, 250-500 µg/mL each, 1 mL/min, 65°C, Detect. UV (233 nm)

Highlights

- The resolution of peaks 1 and 2, not previously obtained, was achieved using the ZORBAX 300SB-C18 at low pH.
- The steric protecting groups of the ZORBAX SB-C18 stabilize the bonded phase, permitting stable and reproducible chromatography at $\text{pH} \leq 1$.



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
5988-6344EN



Agilent Technologies