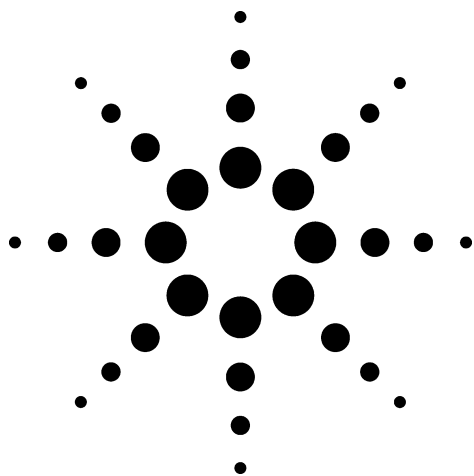
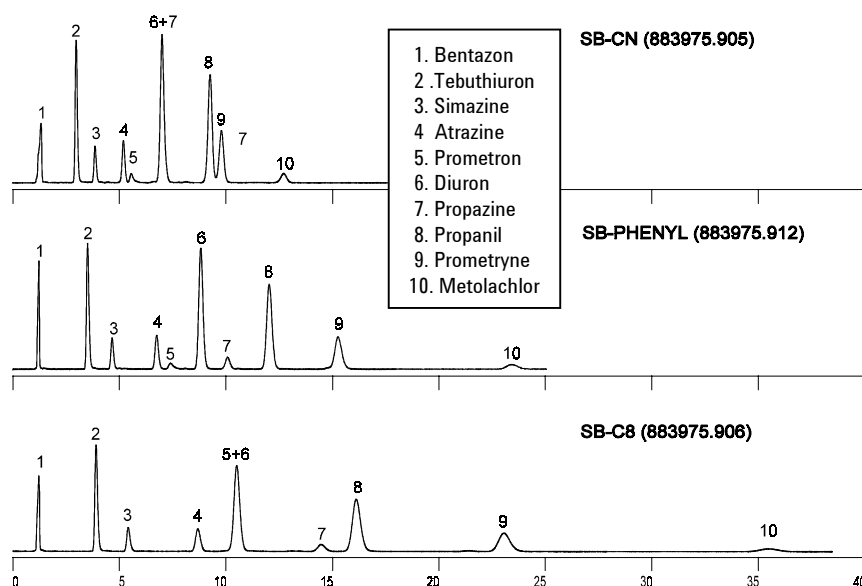




# Herbicides

## Examples of Bonded-Phase Selectivity Differences

Application  
Environmental  
Robert Ricker



Conditions:  
ZORBAX StableBond® Columns (4.6 x 150 mm) (Agilent P/N: above)  
Mobile Phase: 35% ACN, 65% H<sub>2</sub>O, 1mL/min

### Highlights

- Useful differences in band spacing can often be obtained by varying column type, keeping mobile-phase conditions constant.
- For the same mobile phase, stationary-phase strength (retentivity) is in the order C8>Phenyl>CN for these sterically protected bonded phases.
- Long-term stability and reproducibility are provided by ZORBAX StableBond columns.
- A wide range of pesticides can be separated with excellent peak efficiency and peak shape.



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](http://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  
5988-6340EN



**Agilent Technologies**