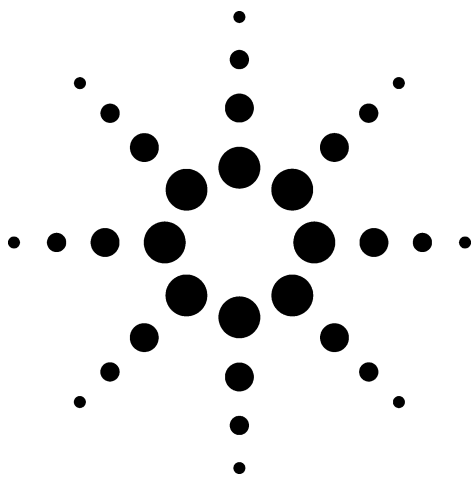




High-Speed Separation of Pesticide Mixture

Application
Environmental
Robert Ricker

Columns with particle sizes under 5 μm and columns with dimensions up to ten-times smaller than standard analytical sized columns are ideal for high-speed analyses. In this application, smaller particles and column dimensions are used to resolve six pesticides in less than two minutes.

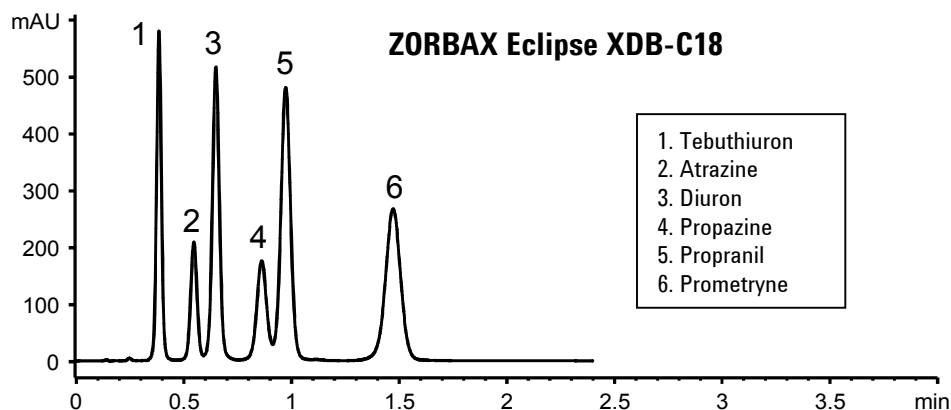
Extra-column volume is a crucial factor in chromatographic performance when using low volume columns. In this example, no modification to a modern instrument (plumbing or flow cell) is necessary.

Highlights

- Reducing column length and particle size simultaneously can:
 - Reduce analysis time
 - Maintain resolution
 - Reduce solvent use
- Higher flow rates are possible due to decreased pressure when using ZORBAX Rapid Resolution cartridge columns.

Operating Conditions:

HPLC System: Agilent 1100 with quaternary pump
Column: ZORBAX Eclipse XDB-C18 Rapid-Resolution (3.5 μm) Cartridge-Column, 4.6 x 30 mm
Agilent Part No. 931975-932
Mobile Phase: MeOH: water (60:40)
Detection: UV 254 nm with standard flow cell (13 μL)
Flow: 2 mL/ min.
Temperature: ambient



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
5988-6347EN



Agilent Technologies