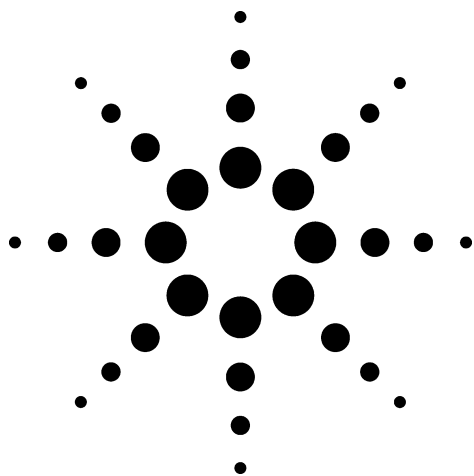
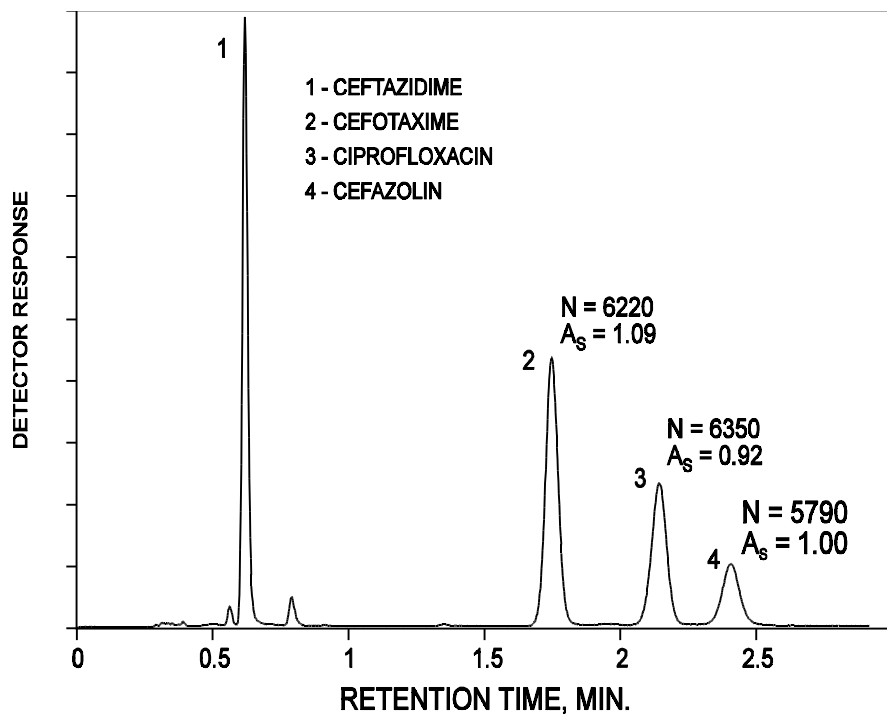




High-Speed Separation of Antibiotics Mixture

Application
Pharmaceutical
Robert Ricker



Highlights

- Excellent peak shape and efficiency from a 75-mm long column at 3.0 ml/min.
- SB-C8 bonded-phase provides reproducible and stable methods for pharmaceuticals.

Conditions:
ZORBAX SB-C8 (3.5µm); 4.6 x 75 mm (Agilent P/N: 866953-906)
Mobile Phase: 8.0% acetonitrile/92% 0.1% aqueous TFA
Flow Rate: 3.0mL/min.; Temperature: 60°C; UV detector, 260 nm
Sample: 1 µL containing 0.40, 0.36, 0.10 and 0.35 µg ea. of 1-4, resp.



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
5988-6379EN



Agilent Technologies