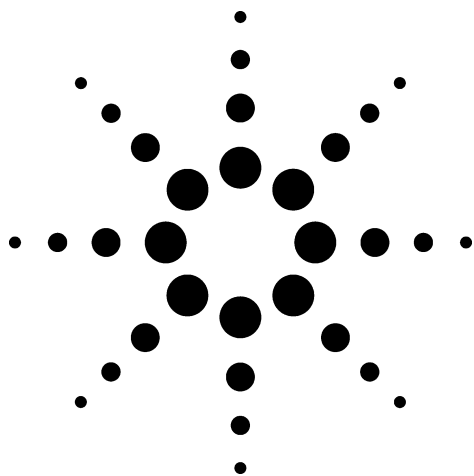


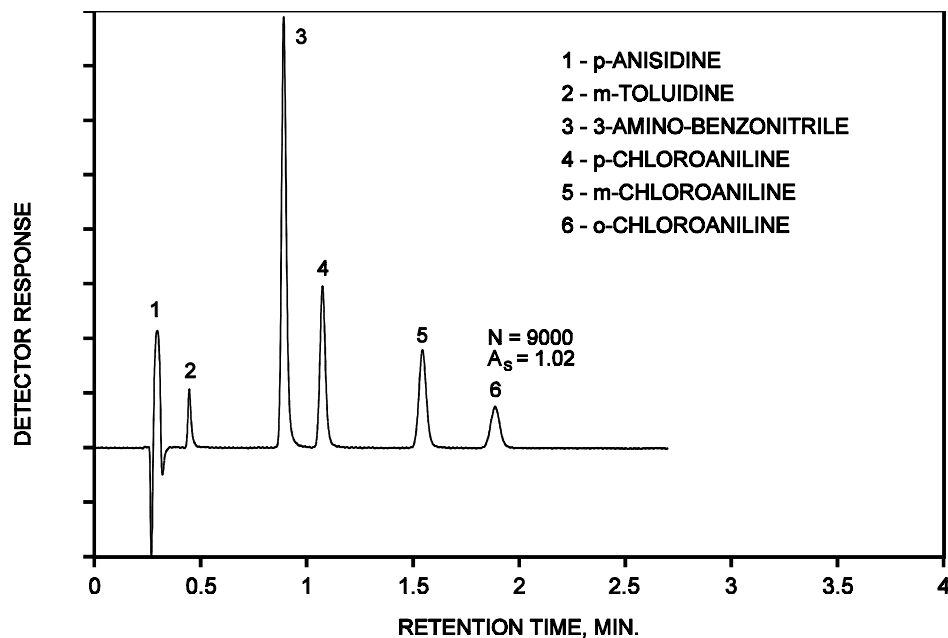


Rapid Separation of Substituted Anilines

Application

Organic Chemicals

Robert Ricker



Highlights

- Isocratic separation achieved in 2 minutes on column packed with 3.5µm ZORBAX SB-C8.
- Excellent peak efficiencies and peak shapes at 3.0 ml/min. with modest back pressure at 60°C.

Conditions: Instrument: HP-1050
ZORBAX SB-C8, 4.6 x 75 mm (Agilent P/N: 866953-906)
Mobile Phase: 20% ACN/80% 25 mM phosphate buffer, pH 2.5
Flow Rate: 3.0 mL/min.; Temp.: 60°C; Pressure: 111bar; UV 254 nm



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
5988-6367EN



Agilent Technologies