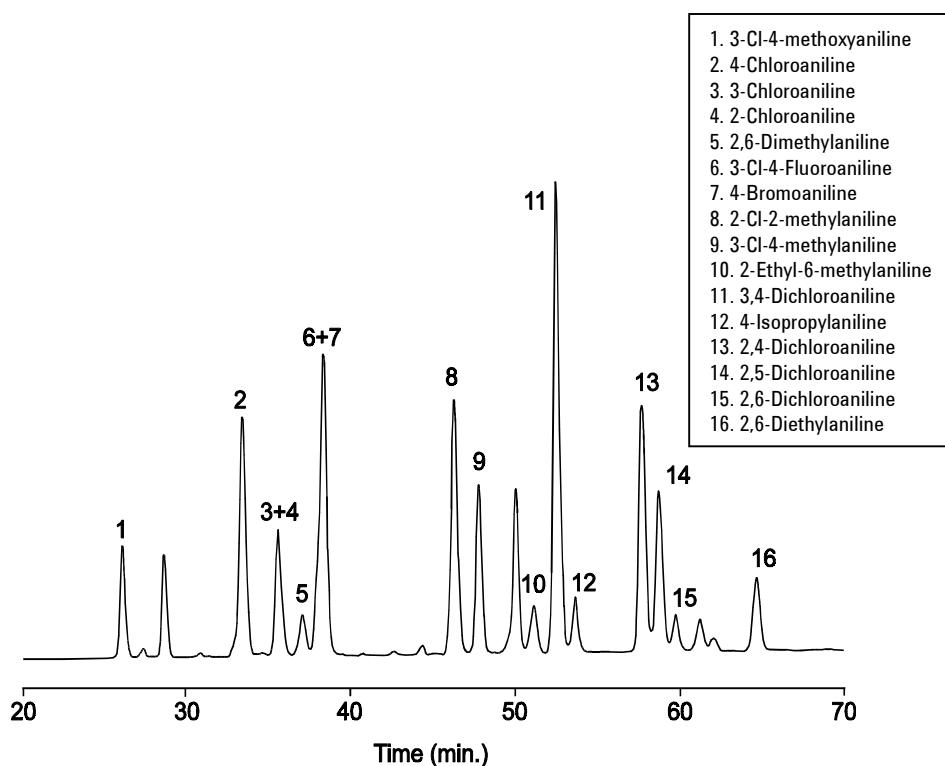




Reversed-Phase Separation 16 Substituted Anilines

Application
Environmental
Robert Ricker



Courtesy of Dr. Schlett, Gelsenwasser A.G.

Highlights

- Separation of 16 environmentally important substituted-anilines.
- Excellent peak shape is afforded by improved silanol interactions.
- ZORBAX StableBond ensures reproducible results time-after-time.

Conditions:

ZORBAX SB-C18 (3 x 250 mm) (Agilent P/N: 880975-902)

Mobile Phase: A=20mM Potassium Acetate, pH 6.5; B=ACN

t = 0 10% B, hold 2 min.

t = 2 10% to 45% B over 68 min.

t = 70 45% B, hold 15 min.

t = 85 45% to 90% B over 1 min.

t = 86 90% to 95% B over 7 min

Equilibration = 20 min., Posttime = 10 min

Injection 25µl, 0.35 mL/min, Ambient, Detect. UV (254 nm) Ref (460 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
5988-6343EN



Agilent Technologies