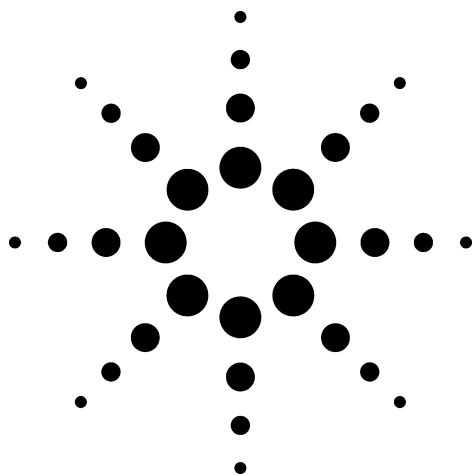
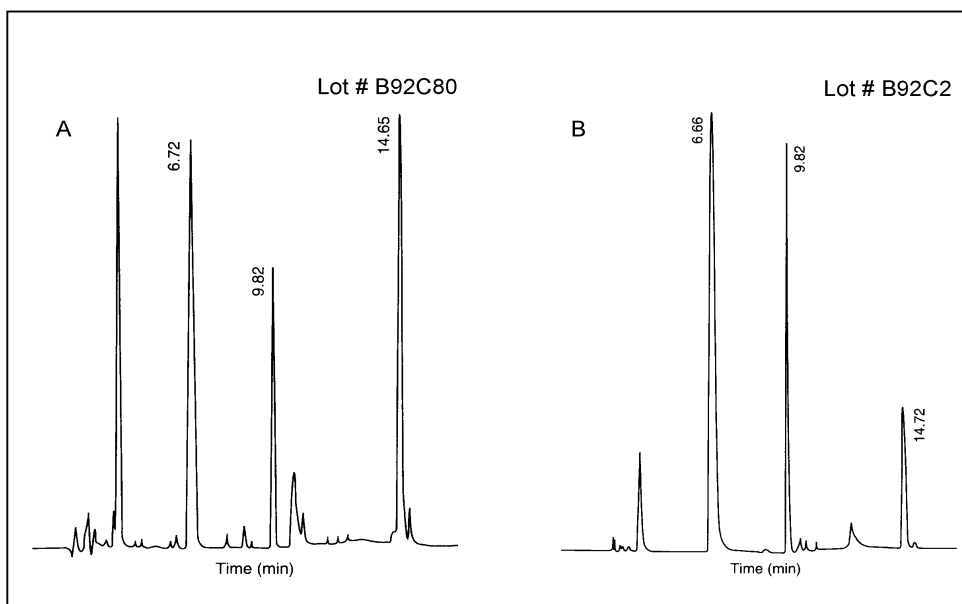




Lot-to-Lot Reproducibility Neuro EndoPeptidase (NEP) Inhibitor

Application
Technical
Robert Ricker



Conditions:
ZORBAX Rx-C18 (4.6 X 150 mm) (Agilent P/N: 883967-902)
1 mL/min; 40°C; Detect. UV(254 nm)
Gradient: 38-70% B in 10 min.
Mobile Phase: A = 0.1 M NaClO₄, pH 2.3
B = 26% A + 74% ACN

Highlights

- In this gradient-elution separation on different lots of ZORBAX Rx-C18 silica, there was less than 1% difference in elution times.
- High reproducibility in elution patterns of basic drugs demonstrates the careful lot-to-lot manufacture of ZORBAX Rx and StableBond columns.
- Consistent elution and good peak shape of these 'silanol-sensitive' basic drugs results from the extremely homogenous silanol surface and low acidity of ZORBAX Rx-SIL.



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
5988-6472EN



Agilent Technologies