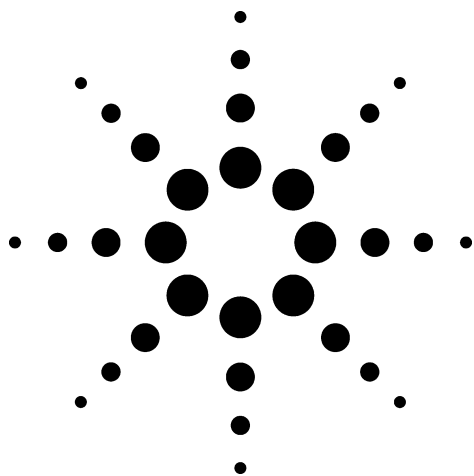
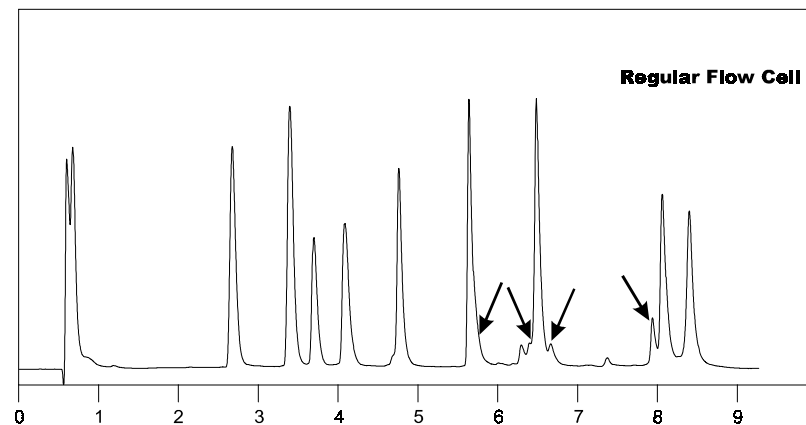
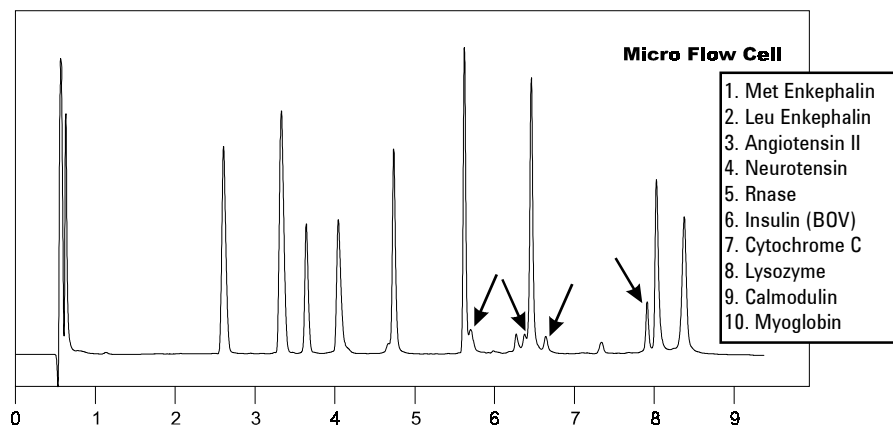




Comparison of Micro and Regular Flow Cell Using ZORBAX 300SB-C8 (3.5 μ m)

Application
Technical
Robert Ricker

The use of a micro flow cell is suggested when using small particle-size column packings. Comparison of flow cells using a 3.5 μ m short ZORBAX 300SB-C8 illustrates reduction of band broadening and improvement of peak resolution for proteins and peptides using a micro flow cell.



Conditions:

ZORBAX 300 SB-C8 (3.5 μ m; 4.6 x 50 mm) (Agilent P/N: 865973-906)

Mobile Phase: A: 95:5 Water:ACN with 0.10% TFA (v/v%) B: 5:95 Water:ACN with 0.085% TFA (v/v%) Gradient: 10-60% B in 10 minutes, posttime 5 minutes

Injection: 10 μ L, 1-6 μ g protein, 1 mL/min, 35°C, Detect. UV(215 nm) micro flow cell

Highlights

- ZORBAX 300SB-C8 (3.5 μ m) with 50-mm column length can reduce time of analysis for peptides and proteins without sacrificing efficiency.
- By reducing the flow cell size, extra-column volume effects are minimized with small particle-size packing.



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
5988-6435EN



Agilent Technologies