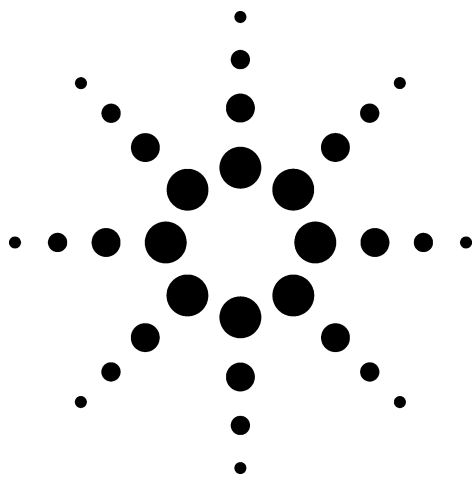




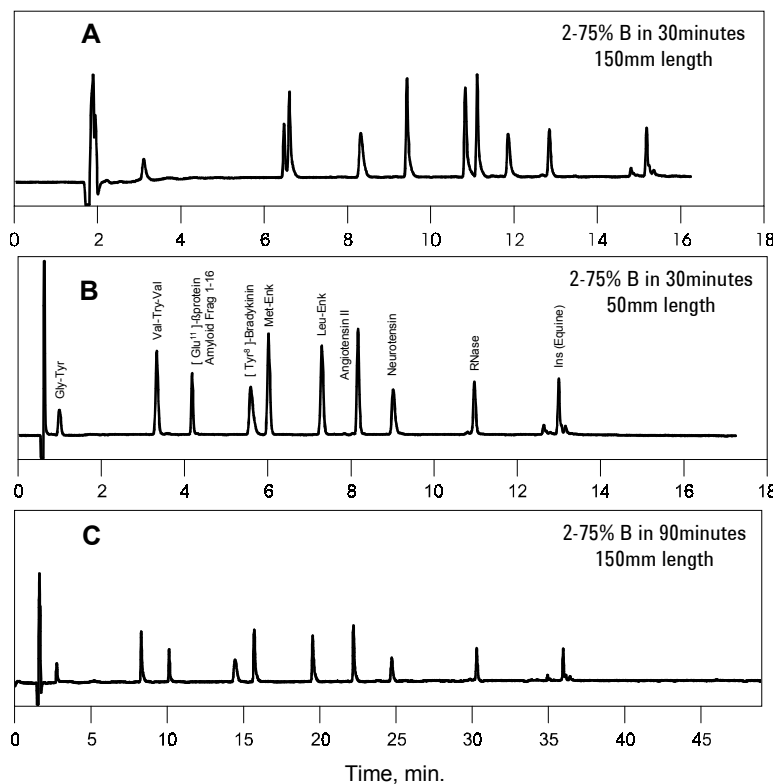
Improving Gradient Resolution by Reducing Column Length: ZORBAX 300SB-C8

Application
Technical
Robert Ricker

Improvement in gradient resolution may be accomplished by decreasing column length (column volume, V_m) without increasing analysis time. A sample containing peptides and proteins was chromatographed on ZORBAX 300SB-C8's of different lengths, keeping all other parameters constant. The results in panel B (50-mm length) show an increase in resolution for most components when compared to Panel A (150-mm column length). Panel C illustrates the increase in run time necessary to achieve the improved separation using a 150-mm column.

Highlights

- Resolution, selectivity, and signal response can be improved by reducing column length instead of changing gradient conditions.
- ZORBAX 300SB-C8 columns are highly stable and reproducible batch-to-batch and for various column configurations.



Conditions:
ZORBAX 300SB-C8, (3.5µm; 4.6 x 50 mm, 150mm) (Agilent P/N: 865973-906, 863973-906)
Mobile Phase: A: 95:5 Water:ACN, 0.10% TFA (v/v%) B: 5:95 Water:ACN, 0.085% TFA (v/v%)
Gradient: as indicated in panels
Injection: 10µL, 1-6 µg protein, 1 mL/min, 35°C, Detect. UV(215 nm) micro flow cell



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
5988-6436EN



Agilent Technologies