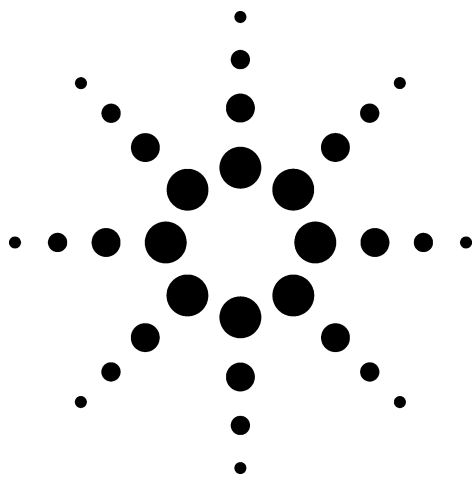




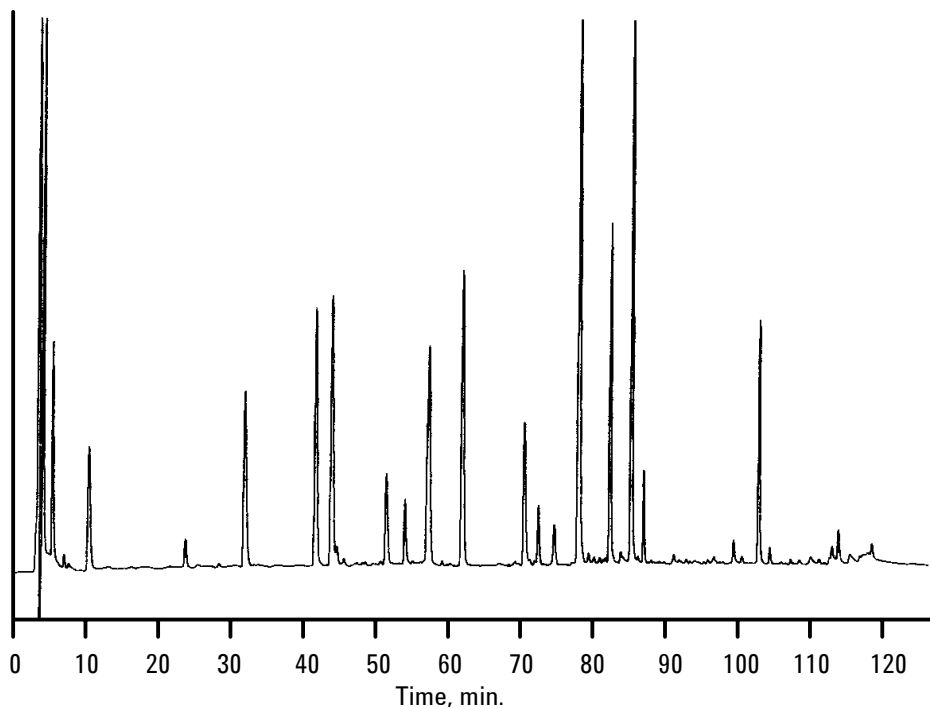
High Resolution Microbore Reversed-Phase HPLC hGH Tryptic Digest

Application
Biochemical
Robert Ricker

The multiple peptide peaks resulting from reversed-phase HPLC of protein digests provide important information about a protein. Additionally, the peaks may be collected and analyzed further (e.g., through mass spectrometry). It is, therefore, important that the peptide sample be chromatographed with high resolution and with reproducibility run-after-run.

Highlights

- *ZORBAX StableBond C18 columns provide excellent resolution of peptides.*
- *Availability of ZORBAX StableBond packings in a 1.0 mm i.d. microbore column allows dependable HPLC results under sample-limited conditions.*
- *ZORBAX StableBond chemistry has excellent stability, especially at low pH.*



Conditions:
ZORBAX 300SB-C18 (1.0 x 250 mm) (MicroTech Scientific)
Mobile Phase: A: H₂O / ACN / TFA 98/2/0.1 B: ACN H₂O / TFA 90/10/0.09
Gradient: 0-60% B at 0.5% B / min.
Injection: 10 µL (70 pmole), 50 µL/min, 50°C, Detect. UV(214 nm)
Sample: Courtesy of Genentech Inc., South San Francisco



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
5988-6318EN



Agilent Technologies