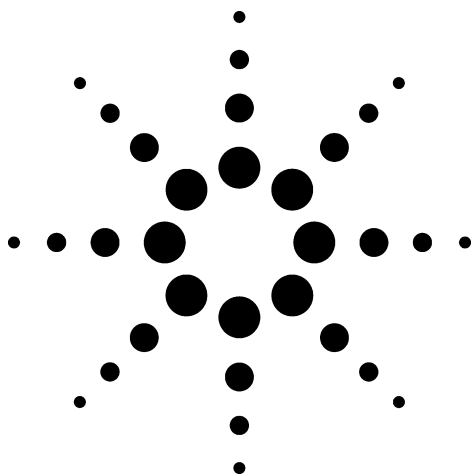




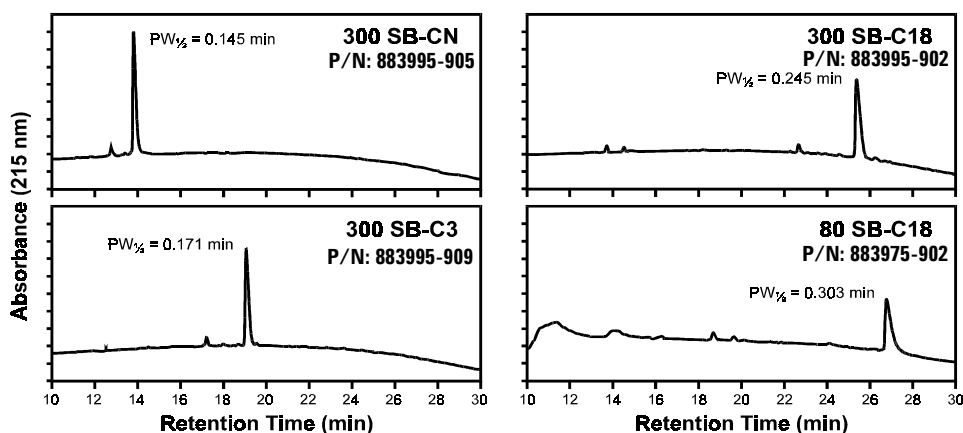
Separation of a Hydrophobic Peptide on Various Bonded Phases

Application
Technical
Robert Ricker

The retention and peak width of the hydrophobic peptide is greater on the more hydrophobic long-chain alkyl bonded-phase packing (300Å, SB-C18), compared to the more polar 300Å, SB-C3 and 300Å, SB-CN column packings. Retention and peak width is greatest on a small pore size, 80Å, (high surface area) SB-C18 column packing.

Highlights

- A range of low pH, high-temperature stable-bonded phases are available; the sterically-protected StableBond column packings.
- The broad range of StableBond column packings permit the choice of bonded-phase functionality that is best for each sample.



Sequence: Ac-Lys Lys Gly (Leu)₁₆ Lys Lys Ala-amide

Courtesy of Drs. R.S. Hodges and C. Mant, Dept. of Biochemistry, University of Alberta, Edmonton, Alberta, Canada

Conditions: LC: Hewlett-Packard 1090 Series II
Columns: 4.6 x 150 mm
Mobile Phase: 5 µg peptide; 40°C; 1.5 ml / min,
25-95% B in 30 min.; A – 0.1% TFA / water, B – 0.09% TFA / ACN
 $PW_{1/2}$ = Peak Width 1/2 peak height



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
5988-6445EN



Agilent Technologies