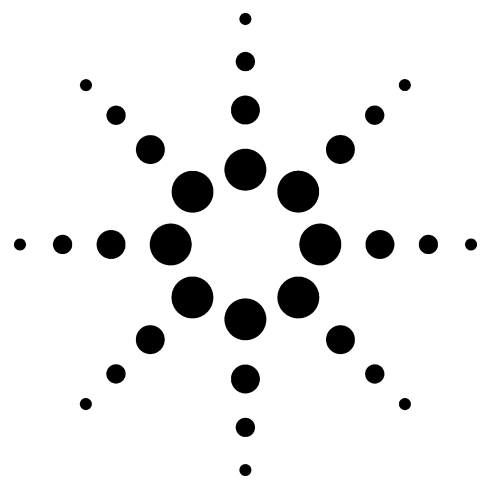
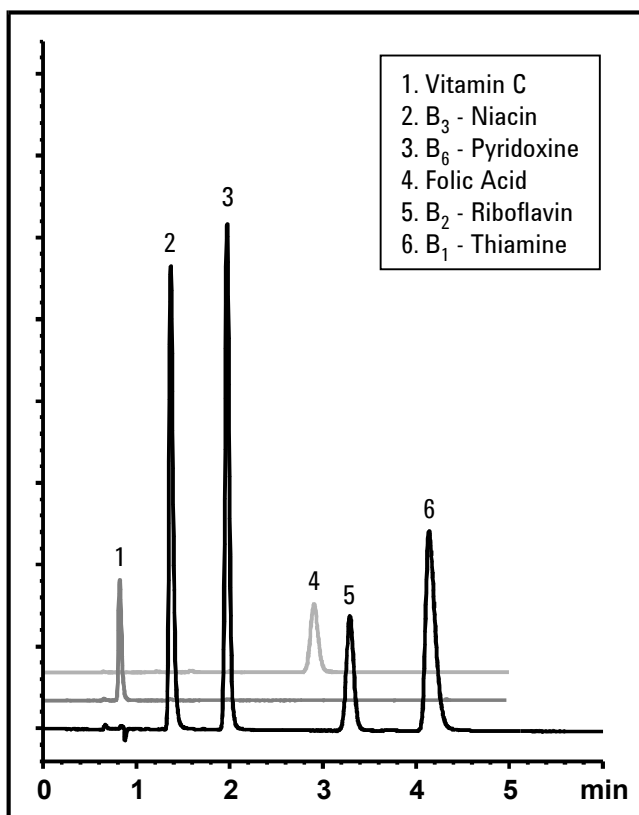




High-Speed Separation of Water-Soluble Vitamins with Ion-Pairing Chromatography

Application
Food Analysis
Robert Ricker

The rapid separation and detection of vitamins is becoming increasingly important to both food and pharmaceutical industries, as well as basic research. Water-soluble vitamins can be easily separated using reverse-phase ion-pairing chromatography. Four B vitamins, folic acid and vitamin C were analyzed in less than 5 minutes using a 75 mm SB-C8 column.



Conditions:
ZORBAX SB-C8 (4.6 x 75 mm, 3.5 μ m) (Agilent P/N: 866953-906)
Mobile Phase: 10mM Hexane Sulfonate with 0.1% Phosphoric Acid:MeOH (74:26)
Injection volume 10 μ L, 1 mL/min, Ambient, Detect: UV(245 nm)

Highlights

- Short run times and excellent peak shape of water-soluble vitamins are achieved using a low-pH mobile phase and a short (75 mm) ZORBAX SB-C8 column with small-particle packing (3.5 μ m).
- ZORBAX StableBond technology offers reproducibility and column stability even at low pH (0.1% phosphoric acid) and with ion-pairing agents.



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
5988-6363EN



Agilent Technologies