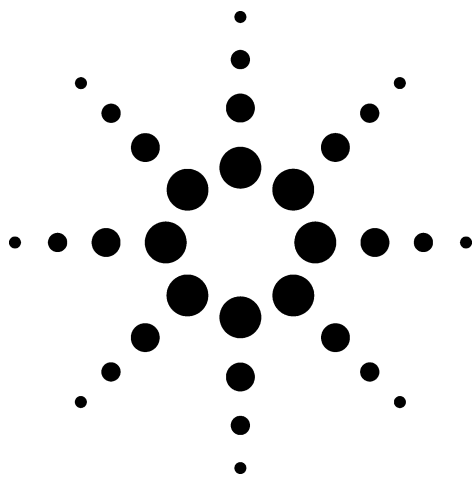




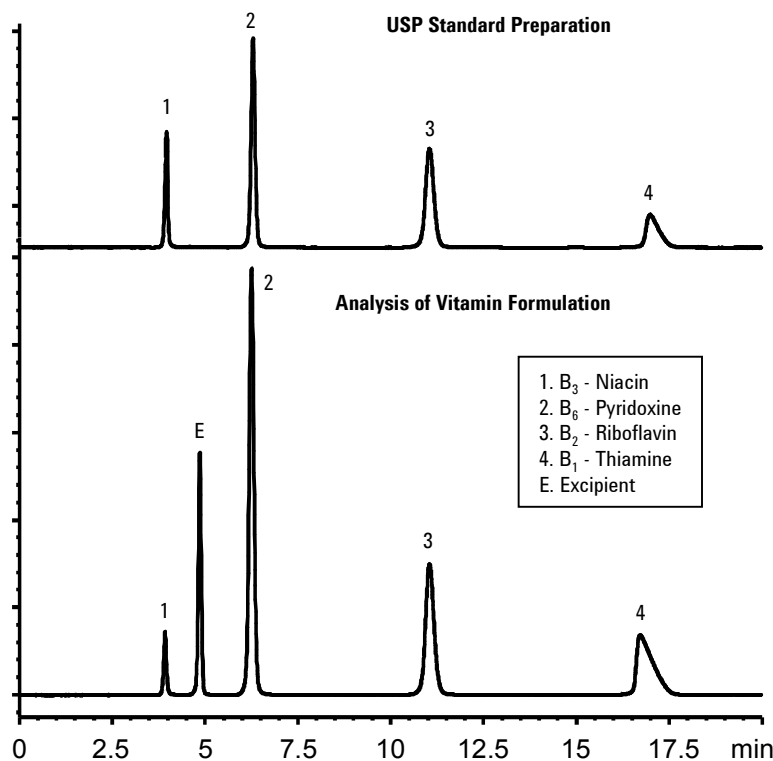
Separation of Water-Soluble Vitamins Using the USP 23 Method

Application
Food Analysis
Robert Ricker

Recently, USP-recommended methods for vitamin analysis have become increasingly important. Analysis and quantitation of water-soluble vitamins using USP-recommended chromatographic methods can become difficult when column performance is questionable. Routine analyses performed by analysts in the pharmaceutical industry require the use of reproducible and stable columns.

Highlights

- *StableBond columns provide analysts with long-term stability and reliability for the demands of USP-generated methods.*
- *Lot-to-lot reproducibility of StableBond column packings assures the chromatographer of consistency in USP-generated methods.*
- *Details of the USP Method appear on page 2168 of USP 23.*



Conditions:
ZORBAX SB-C18 (L1 packing) (4.6 x 250 mm) (Agilent P/N: 880975-902)
Mobile Phase: 7.2 mM Hexane Sulfonate:MeOH:Acetic Acid (73:27:1) (ratio to 101%)
Injection: 20 µL, 1 mL/min, 30°C, Detect. UV(280 nm)



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
5988-6364EN



Agilent Technologies