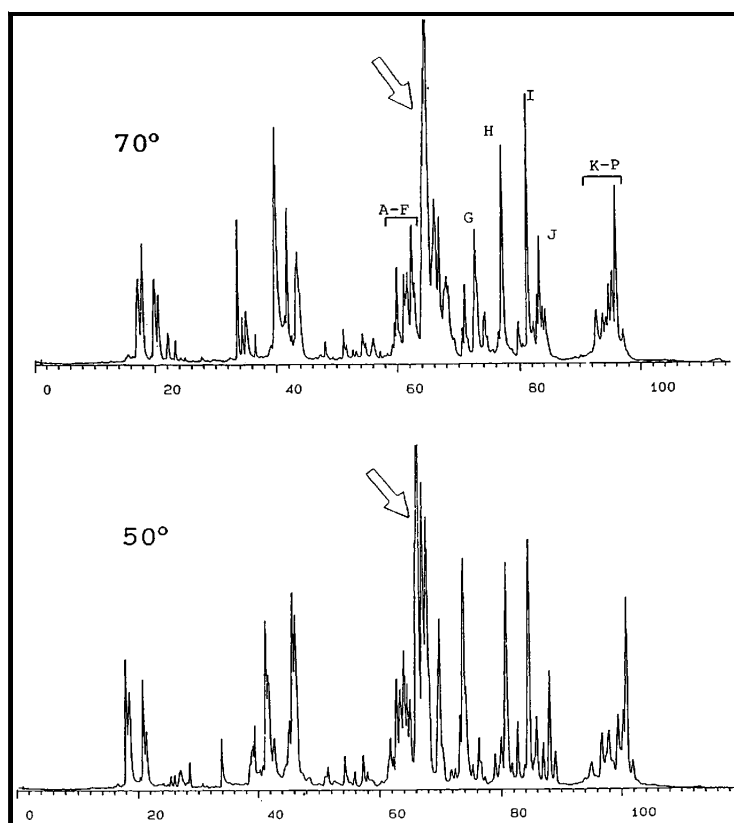


Effect of Temperature on Separation of a Wheat Protein

Application
Food Analysis
Robert Ricker



Conditions:
ZORBAX 300 SB-C8, 4.6 x 150 mm (Agilent P/N: 883995-906)
1 ml/min., 15 µl of Neepawa Wheat Extract
Abs. 210 nm 23-48% B in 120 min.
A=0.1% TFA/water, B=0.1% TFA/ACN

Highlights

- *Wide-pore ZORBAX 300 SB-C8 demonstrates high peak capacity for complex samples of plant protein.*
- *Temperature can be effectively used to increase resolution in reversed-phase separations of complex samples.*



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
5988-6348EN



Agilent Technologies