



Vitamins

Separation of vitamin E alcohol/ acetate

Application Note

Food Testing & Agriculture

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography using an Agilent CP-Sil 5 CB column separates vitamin E alcohol and acetate in four minutes.



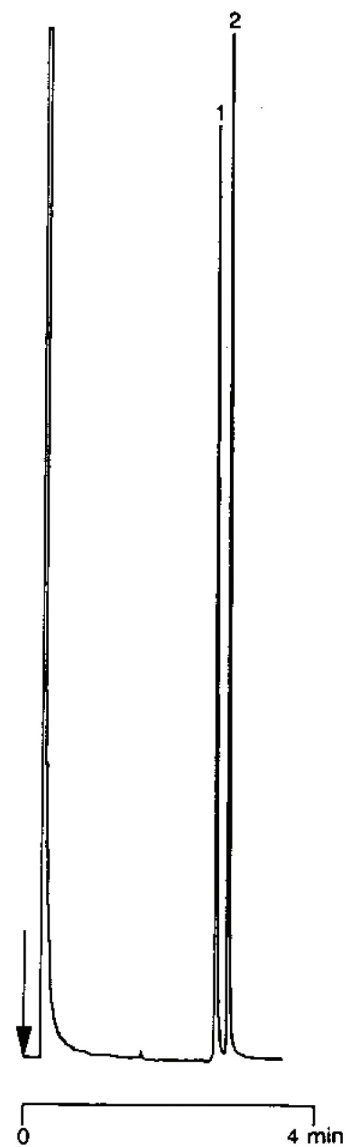
Agilent Technologies

Conditions

Technique	: GC-capillary
Column	: Agilent CP-Sil 5 CB, 0.32 mm x 12.5 m fused silica WCOT CP-Sil 5 CB (df = 0.12 µm) This column is available as a 25m column under Part no. CP7740 (+ retention gap 0.32 mm x 1 m)
Temperature	: 200 °C (0.5 min) → 255 °C, 25 °C/min; 255 °C (1 min)
Carrier Gas	: H ₂ , 25 kPa (0.25 bar, 3.6 psi), 80 cm/s
Injector	: Splitter, 1:15 T = 300 °C
Detector	: FID T = 300 °C
Sample Size	: 1.0 µL
Concentration Range	: 1 mg/mL

Peak identification

1. DL- α -Tocopherol
(Vitamin E alcohol)
2. DL- α -Tocopheryl acetate
(Vitamin E acetate)



www.agilent.com/chem

This information is subject to change without notice.

© Agilent Technologies, Inc. 2011

Printed in the USA

31 October, 2011

First published prior to 11 May, 2010

A00715



Agilent Technologies