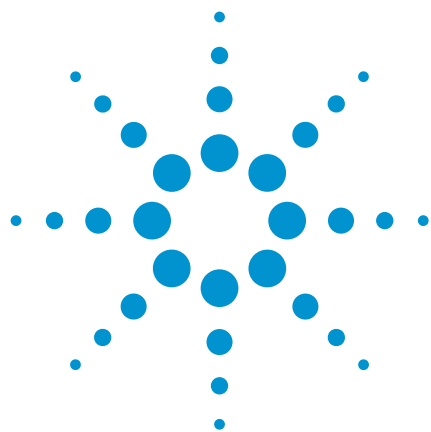




Flavors and aromas

Separation of α -pinene, limonene and linalool

Application Note

Food Testing & Agriculture

Authors

Agilent Technologies, Inc.

Introduction

Agilent CP-Chirasil-Dex CB is one of the few chiral stationary phases that is chemically bonded and combines a high selectivity with a high inertness for optical isomers. The majority of compounds can be analyzed without derivatization. Resolution factors can be improved by using electronic pressure or flow programming.



Agilent Technologies

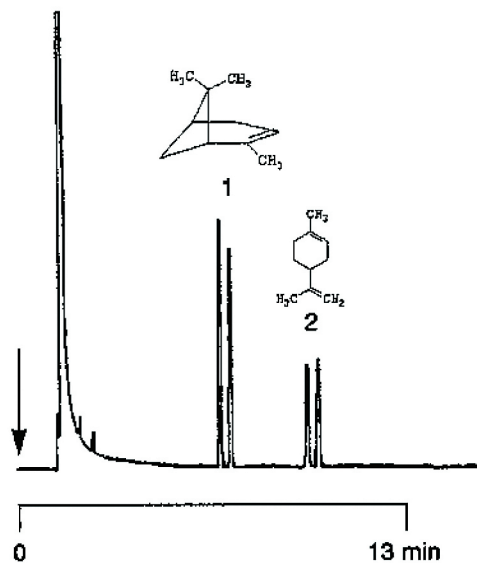
Conditions

Technique	: GC-capillary
Column	: Agilent CP-Chirasil-Dex CB, 0.25 mm x 25 m fused silica WCOT CP-Chirasil-Dex CB (df = 0.25 µm) (Part no. CP7502)
Temperature	: Chromatogram A: 85 °C Chromatogram B: 100 °C
Carrier Gas	: H ₂ , 50 kPa (0.5 bar, 7 psi)
Injector	: Split, 100 mL/min T = 275 °C
Detector	: FID T = 300 °C
Sample Size	: 0.2 µL
Concentration Range	: 0.05 %

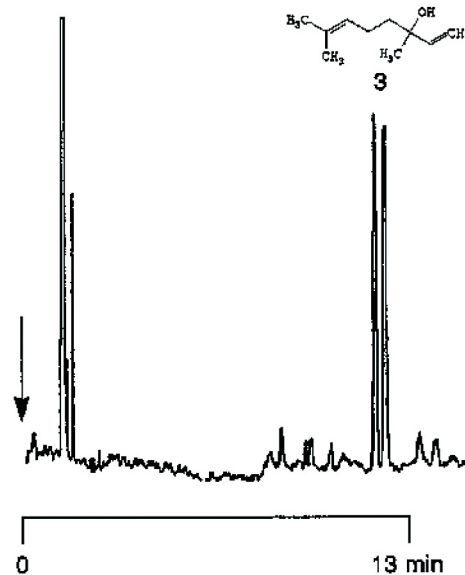
Peak identification

1. α-pinene
2. limonene
3. linalool

Chromatogram A



Chromatogram B



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

A01326



Agilent Technologies