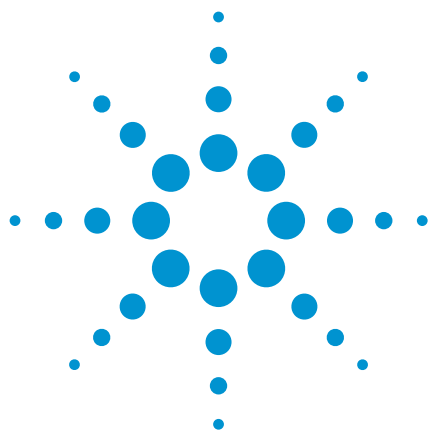




Enantiomers

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Chirasil-DEX CB column separates racemic aromatic alcohols in 14 minutes.



Agilent Technologies

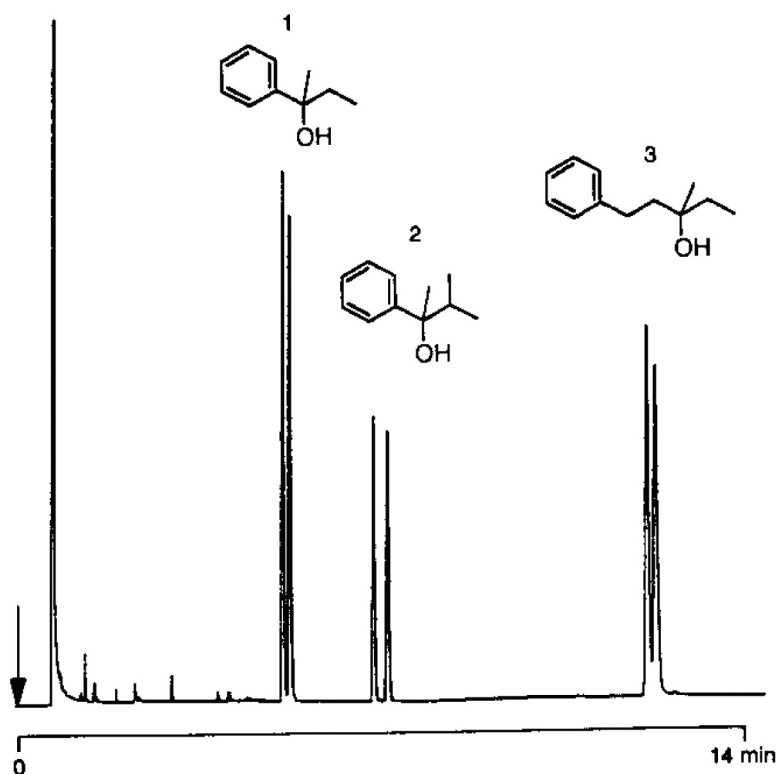
Conditions

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 μ m) (Part no. CP7502)
Temperature : 100 °C $\rightarrow$ 130 °C, 2 °C/min
Carrier Gas : H₂, 100 kPa (1 bar, 14.5 psi)
Injector : Split
Detector : FID

Courtesy : Prof. V. Schurig, Universität Tübingen, Tübingen, Germany

Peak identification

1. (±)-2-phenyl-2-butanol
2. (±)-2-phenyl-3-methyl-2-butanol
3. (±)-3-methyl-5-phenyl-3-pentanol



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

A00975



Agilent Technologies