



Enantiomers

Analysis of racemic aromatic alcohols (phenyl substituted alcohols)

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Chirasil-DEX CB column separates enantiomers of six racemic aromatic alcohols (phenyl substituted alcohols) in 15 minutes.



Agilent Technologies

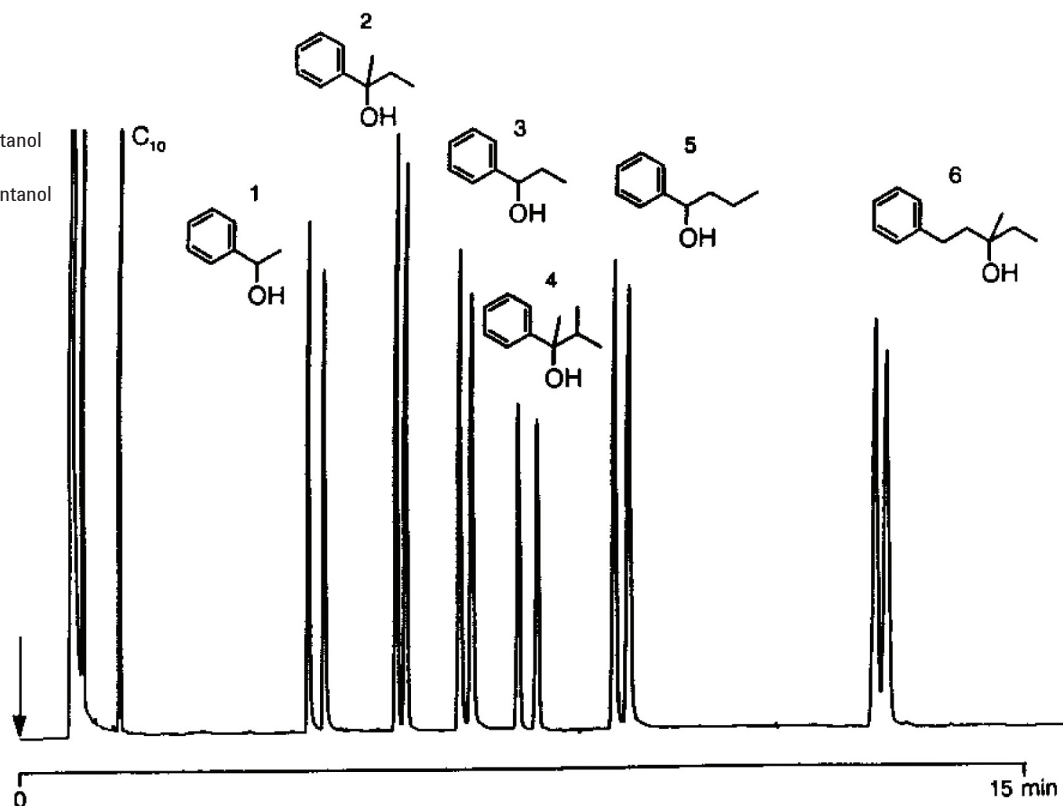
Conditions

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. (\pm)- α -methylbenzyl alcohol
2. (\pm)-2-phenyl-2-butanol
3. (\pm)-1-phenyl-1-propanol
4. (\pm)-2-phenyl-3-methyl-2-butanol
5. (\pm)-1-phenyl-1-butanol
6. (\pm)-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

A00967