



Enantiomers

Analysis of aromatic alcohols

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Chirasil-DEX CB column separates enantiomers of phenyl-3-butanol in 20 minutes.



Agilent Technologies

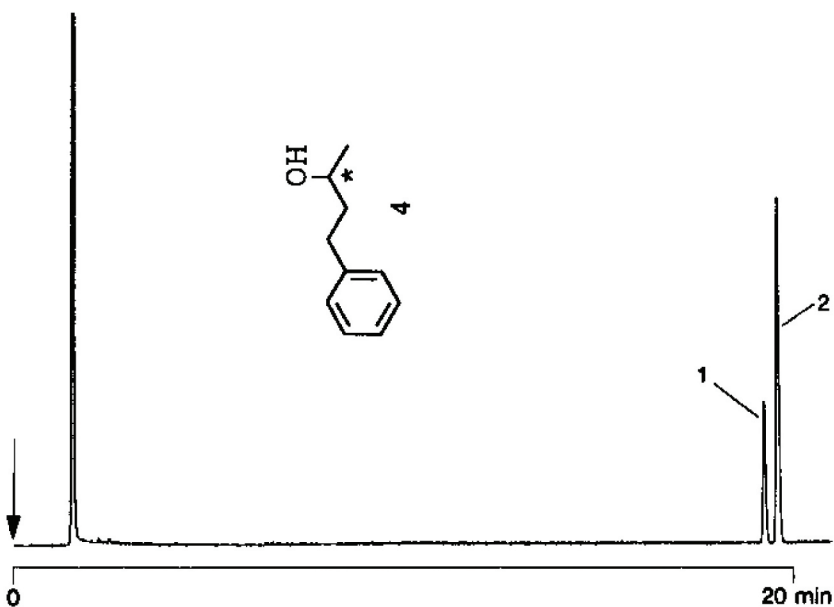
Conditions

Technique : GC-capillary
Column : Agilent CP-Chirasil-DEX CB, 0.32 mm x 25 m fused silica WCOT CP-Chirasil-DEX (df = 0.32 µm)
(Part no. CP7503)
Temperature : 120 °C
Carrier Gas : H₂, 27 kPa (0.27 bar, 4 psi)
Injector : Split, 100 mL/min
T = 300 °C
Detector : FID
T = 270 °C
Sample Size : 2 µL
Concentration Range : 0.2% of each isomer and compound
Solvent Sample : diethyl ether

Courtesy : Inger Cathrine Tveit, Hilde Oppedal and Dr. Anne Fiksdahl, Department of Organic Chemistry, University of Trondheim-NTH, Trondheim, Norway

Peak identification

1. S-1-phenyl-3-butanol
2. R-1-phenyl-3-butanol



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

A00996



Agilent Technologies