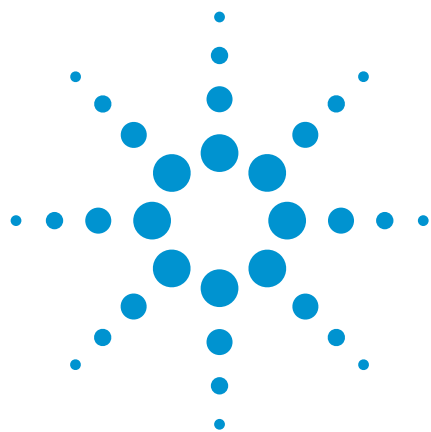




Enantiomers

Enantiomer separation of racemic underivatized diols

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Chirasil-DEX CB column separates enantiomers of racemic underivatized diols in ten minutes.



Agilent Technologies

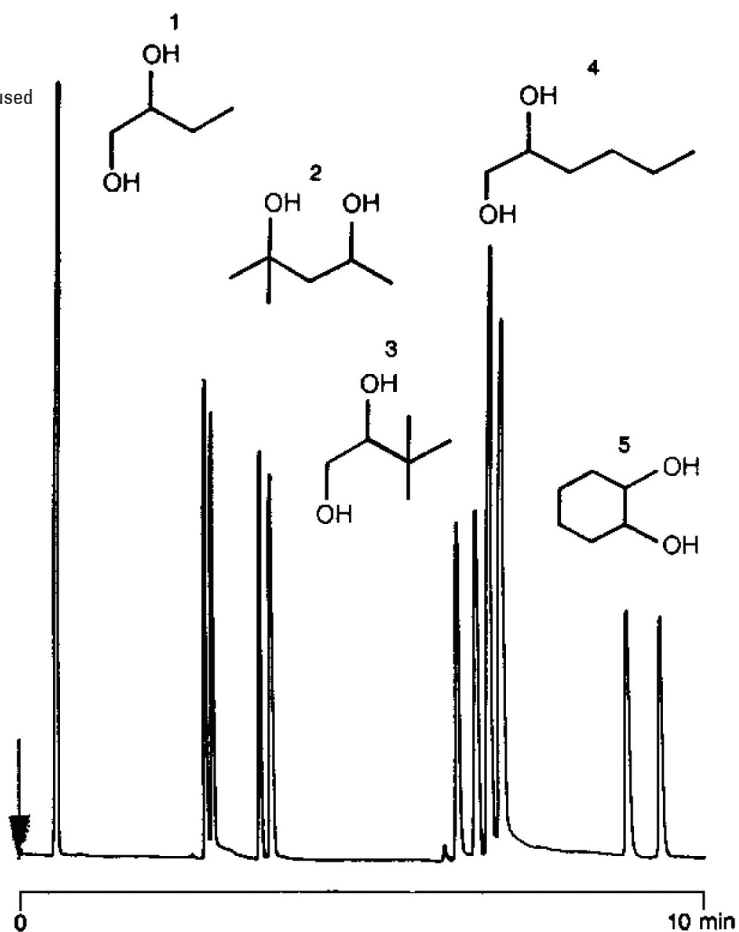
Conditions

Technique : GC-capillary
Column : Agilent CP-Chirasil-DEX CB, 0.25 mm x 25 m fused silica WCOT CP-Chirasil-DEX (df = 0.25 µm) (Part no. CP7502)
Temperature : 90 °C → 120 °C, 2.5 °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. (±)-1,2-butanediol
2. (±)-2-methyl-2,4-pentanediol
3. (±)-3,3-dimethyl-1,2-butanediol
4. (±)-1,2-hexanediol
5. trans-1,2-cyclohexanediol



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

A00971



Agilent Technologies