



Enantiomers

Enantiomer separation of some 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 seven underivatized diols in 20 minutes.



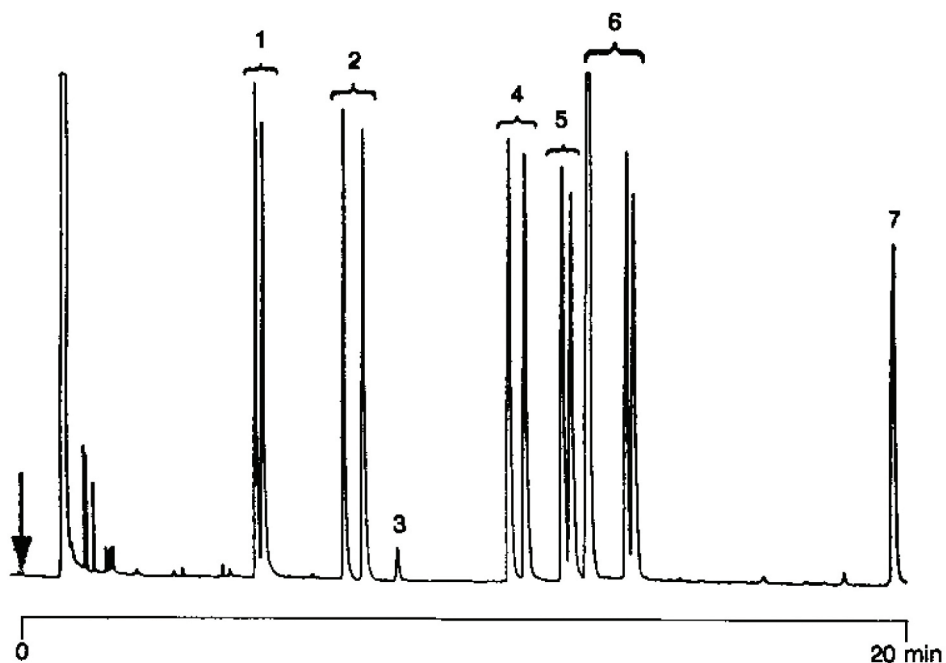
Agilent Technologies

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 : 80 °C (5 min) → 150 °C, 2 °C/min
Carrier Gas : H₂, 70 kPa (0.7 bar, 10 psi)
Injector : Split, 100 mL/min
T = 250 °C
Detector : FID, 4 x 10⁻¹² Afs
T = 300 °C
Concentration Range : 0.01% in CH₃OH

Peak identification

1. 1,2-propanediol
2. 2,3-butanediol
3. meso-2,3-butanediol
4. 1,2-butanediol
5. 1,3-butanediol
6. 2,4-pentanediol
7. 2,5-hexanediol



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

A00968



Agilent Technologies