



Diols

Enantiomer separation of underivatized diols

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Enantiomeric separation of seven underivatized diols by gas chromatography with an Agilent CP-Cyclodextrin-B-2,3,6-M-19 column is achieved in 26 minutes.

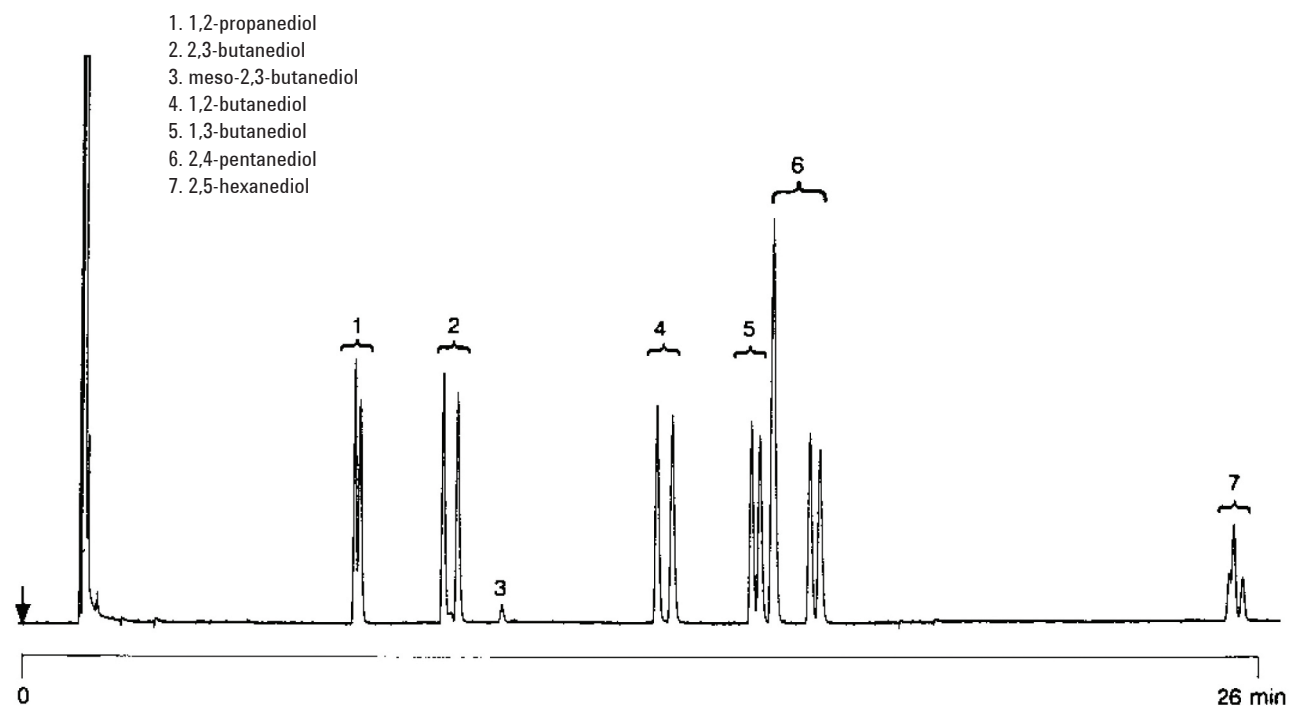


Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Cyclodextrin-B-2,3,6-M-19, 0.25 mm x
25 m fused silica WCOT Cyclodextrin-B-2,3,6-M-19
(df = 25 µm) (Part no. CP7500)
Temperature : 80 °C, 5 min + 2 °C/min → 150 °C
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 = 275 °C
Concentration range : 0.01% in CH₃OH

Peak identification



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

A00498



Agilent Technologies