



Enantiomers

Analysis of epoxides

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Chirasil-DEX CB column separates enantiomers of four epoxides in 14 minutes.



Agilent Technologies

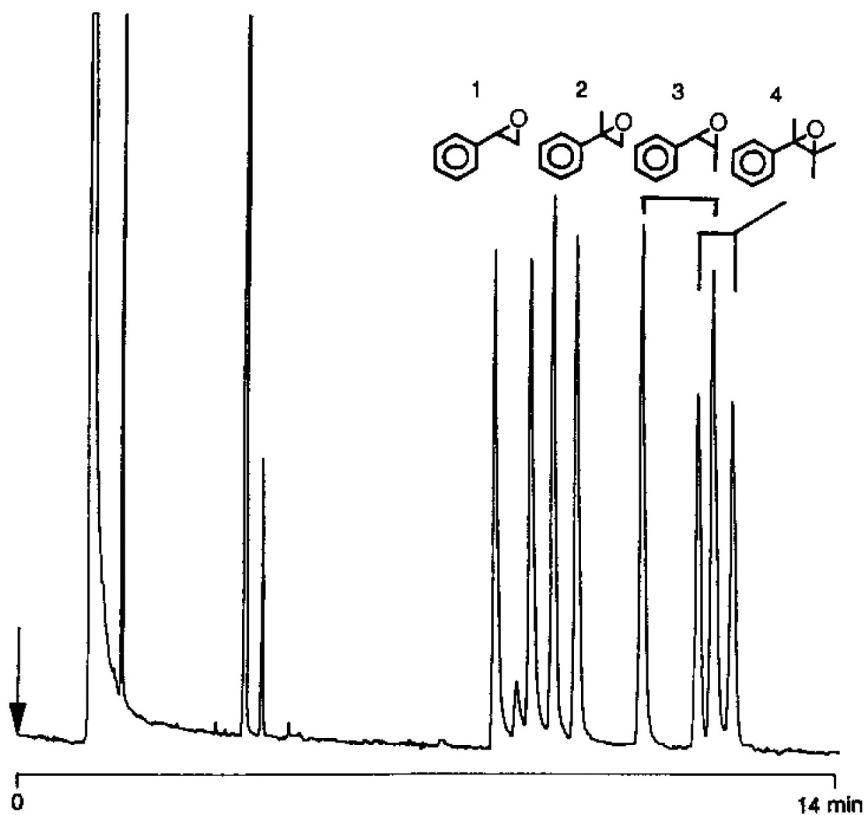
Conditions

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
Carrier Gas : H₂, 60 kPa (0.6 bar, 8.6 psi)
Injector : Split
Detector : FID

Courtesy : Prof. V. Schurig,
Universität Tübingen,
Tübingen, Germany

Peak identification

1. (±)-styrene oxide
2. (±)-2-phenylpropylene oxide
3. (±)-1-phenylpropylene oxide
4. (±)-1,1,2-trimethyl-2-phenyl oxirane



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

A00963



Agilent Technologies