



Enantiomers

Enantiomer separation of racemic alkyl bromides

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Chirasil-DEX CB column separates enantiomers of racemic alkyl bromides in 20 minutes.



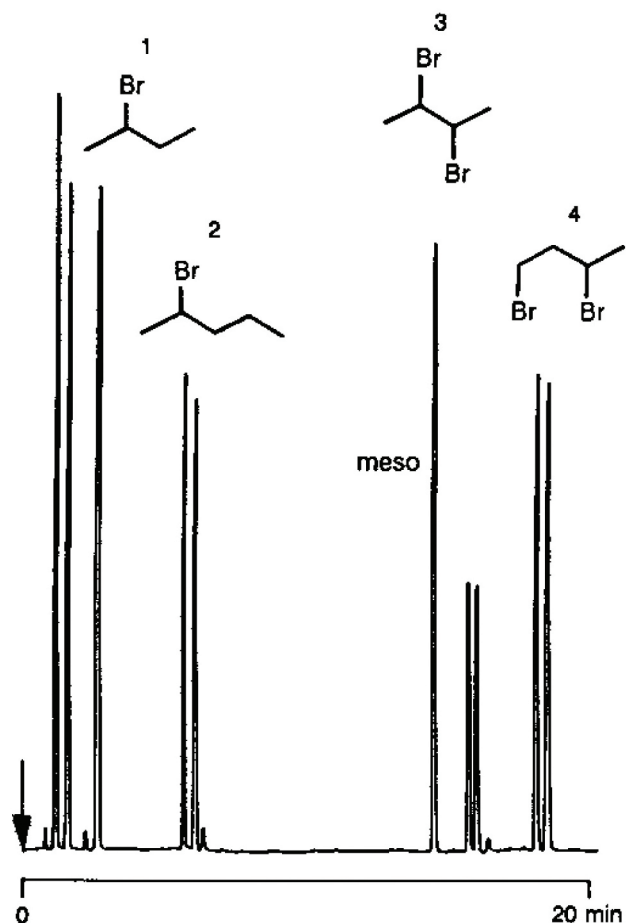
Agilent Technologies

Conditions

Technique	: GC-capillary
Column	: Agilent CP-Chirasil-DEX CB, 0.25 mm x 25 m fused silica WCOT CP-Chirasil-DEX CB (df = 0.25 µm) (Part no. CP7502)
Temperature	: 33 °C → 85 °C, 2.5 °C/min
Carrier Gas	: H ₂ , 100 kPa (1 bar, 14.5 psi)
Injector	: Split
Detector	: FID
Courtesy	: Prof. V. Schurig, Universität Tübingen, Tübingen, Germany

Peak identification

1. (±)-2-bromobutane
2. (±)-2-bromopentane
3. (±)-2,3-dibromobutane
4. (±)-1,3-dibromobutane



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

A00970



Agilent Technologies