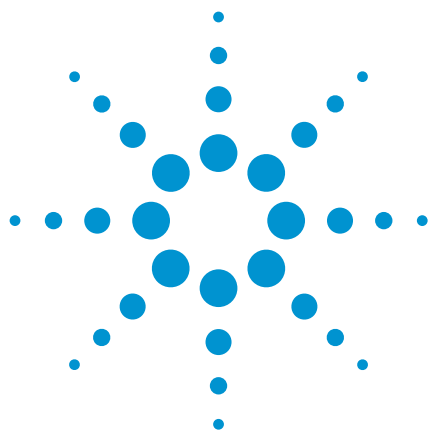




Enantiomers

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Chirasil-DEX CB column separates enantiomers of chlorophenol in 12 minutes.



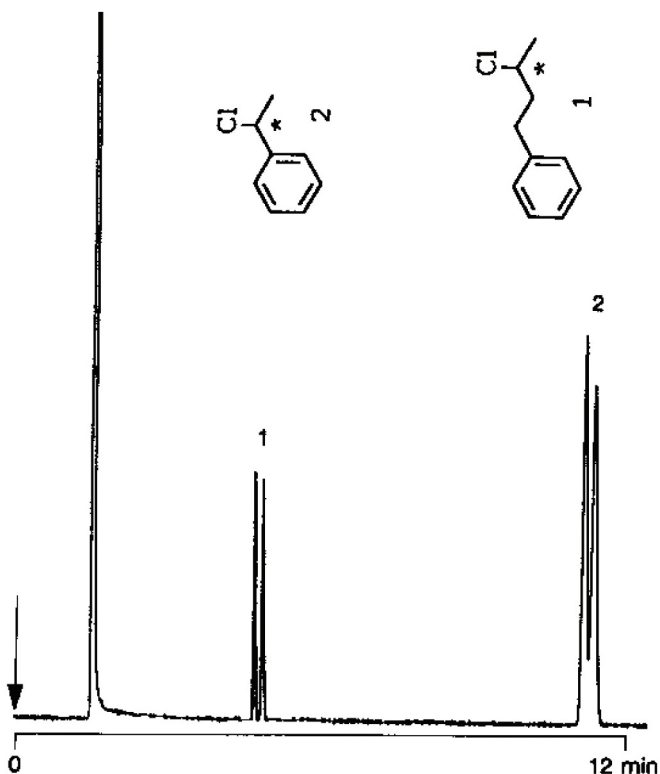
Agilent Technologies

Conditions

Technique	: GC-capillary
Column	: Agilent CP-Chirasil-DEX CB, 0.32 mm x 25 m fused silica WCOT CP-Chirasil-DEX (df = 0.32 µm) (Part no. CP7503)
Temperature	: 120 °C
Carrier Gas	: H ₂ , 27 kPa (0.27 bar, 4 psi)
Injector	: Split, 100 mL/min T = 300 °C
Detector	: FID T = 270 °C
Sample Size	: 2 µL
Concentration Range	: 0.2% of each isomer and compound
Solvent Sample	: diethyl ether
Courtesy	: Inger Cathrine Tveit, Hilde Oppedal and Dr. Anne Fiksdahl, Department of Organic Chemistry, University of Trondheim-NTH, Trondheim, Norway

Peak identification

1. R- and S-(3-chlorobutyl)benzene
2. R- and S-(1-chloroethyl)benzene



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

A00994



Agilent Technologies