

Aromatic hydrocarbons, C₆-C₈

High resolution separation of xylene isomers

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

The cyclodextrin modification of the Agilent CP-Chirasil DEX CB stationary phase makes high resolution separation of xylene isomers possible. The separation of para- and meta-xylene is very good. The bonded chiral phase will guarantee longest possible lifetime.



Agilent Technologies

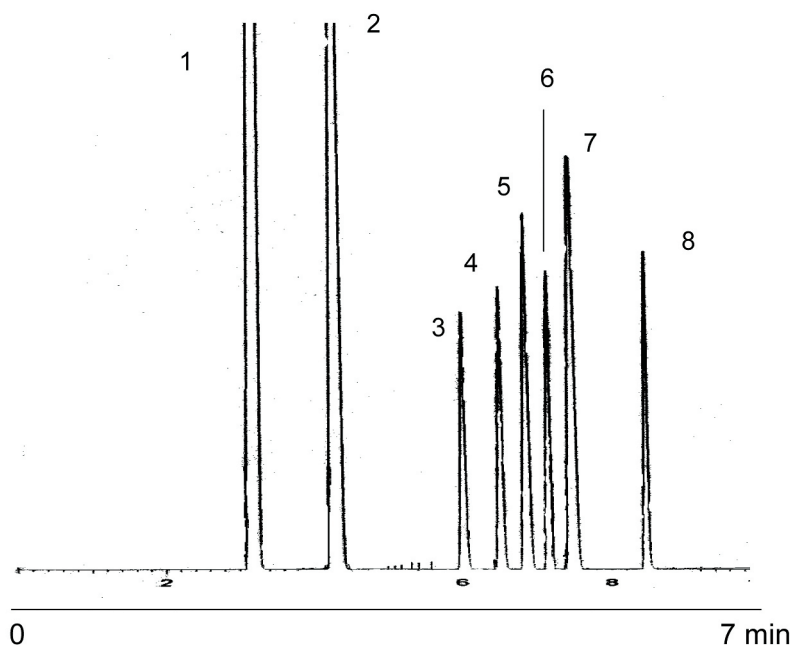
Conditions

Technique : GC
Column : Agilent CP-Chirasil-DEX CB, 0.25 mm x 25 m fused silica (df = 0.25 µm) (Part no. CP7502)
Temperature : 80 °C, (6 min) → 130 °C, 25 °C/min
Carrier Gas : helium, 40kPa, 6 Psi
Injector : Split, T = 210 °C, 1: 20
Detector : FID
T = 230 °C
Sample Size : 0.5 µL
Concentration Range : 10 - 20%

Courtesy : John Valdes, BP Refinery, Texas City Site, Texas City

Peak identification

1. benzene
2. toluene
3. para xylene
4. meta xylene
5. ethyl benzene
6. ortho xylene
7. styrene
8. cumene



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

A01897



Agilent Technologies