



Aromatic hydrocarbons, C₇ – C₉

Analysis of impurities in o-xylene

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Wax 52 CB column separates nine C₇ to C₉ aromatic hydrocarbon impurities in o-xylene in 16 minutes.



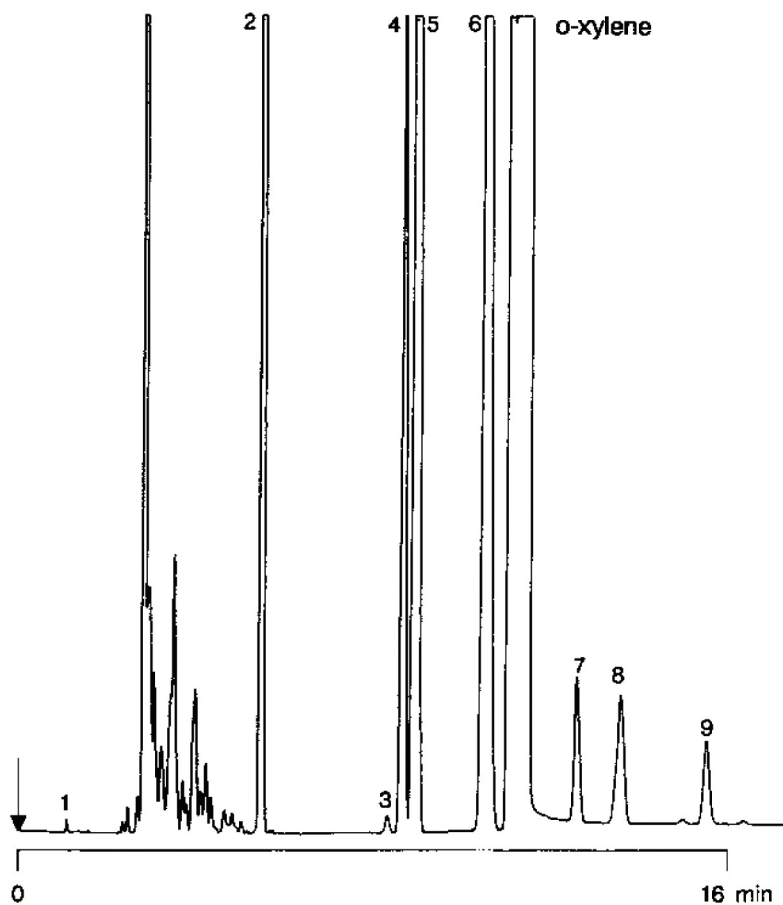
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Wax 52 CB, 0.32 mm x 25 m fused silica
WCOT CP-Wax 52 (df = 1.2 µm) (Part no. CP7763)
Temperature : 100 °C (isothermal)
Carrier Gas : nitrogen
Injector : Splitter 1:20
T = 250 °C
Detector : FID
T = 250 °C

Peak identification

1. C₁ - C₄ paraffins
2. toluene
3. ethylbenzene
4. p-xylene
5. m-xylene
6. iso-propylbenzene
7. n-propylbenzene
8. ethyltoluene
9. styrene



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

A00418



Agilent Technologies