



Aromatic hydrocarbons

Analysis of impurities in p-, m- and o-xylenes

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Xylenes column identifies four impurities in p-, m-, and o-xylene in 15 minutes.



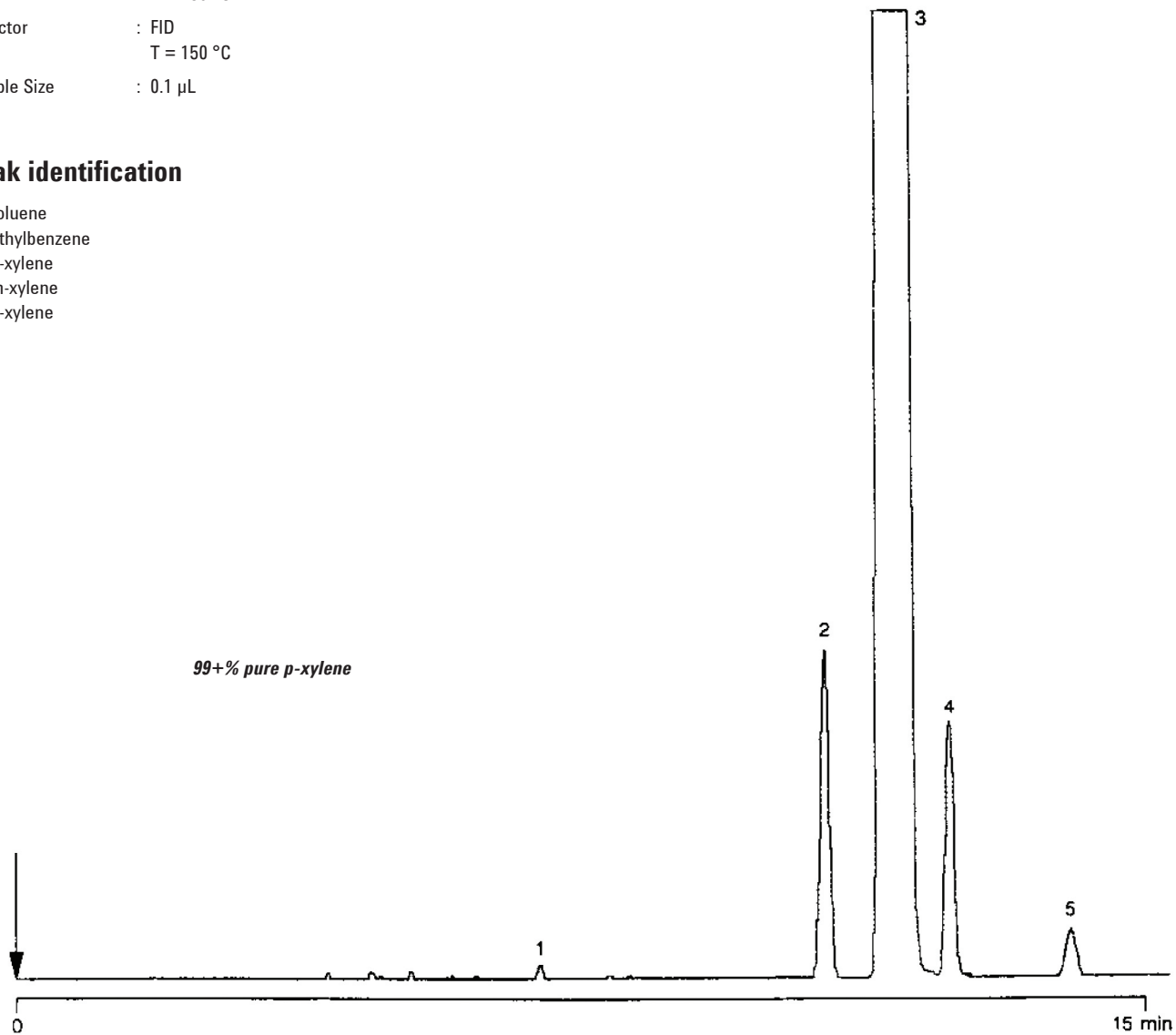
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Xylenes, 0.32 mm x 50 m fused silica
PLOT CP-Xylenes (Part no. CP7426)
Temperature : 35 °C
Carrier Gas : He, 65 kPa (0.65 bar, 9.3 psi)
Injector : Splitter, 60 mL/min
T = 150 °C
Detector : FID
T = 150 °C
Sample Size : 0.1 µL

Peak identification

1. toluene
2. ethylbenzene
3. p-xylene
4. m-xylene
5. o-xylene



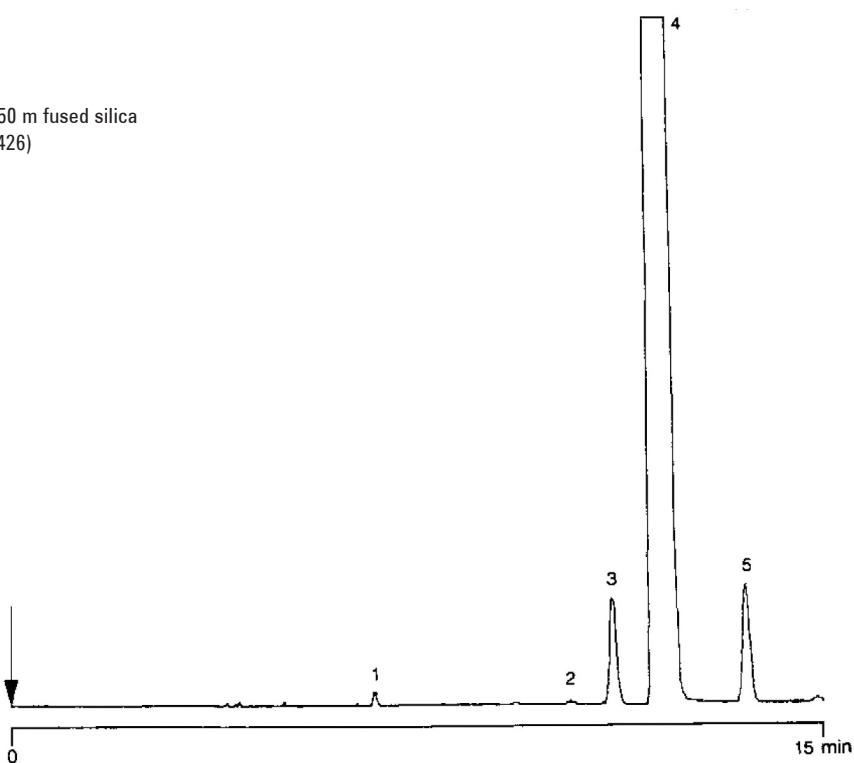
Conditions

Technique : GC-capillary
Column : Agilent CP-Xylenes, 0.32 mm x 50 m fused silica
PLOT CP-Xylenes (Part no. CP7426)
Temperature : 35 °C
Carrier Gas : He, 65 kPa (0.65 bar, 9.3 psi)
Injector : Splitter, 60 mL/min
T = 150 °C
Detector : FID
T = 150 °C
Sample Size : 0.1 µL

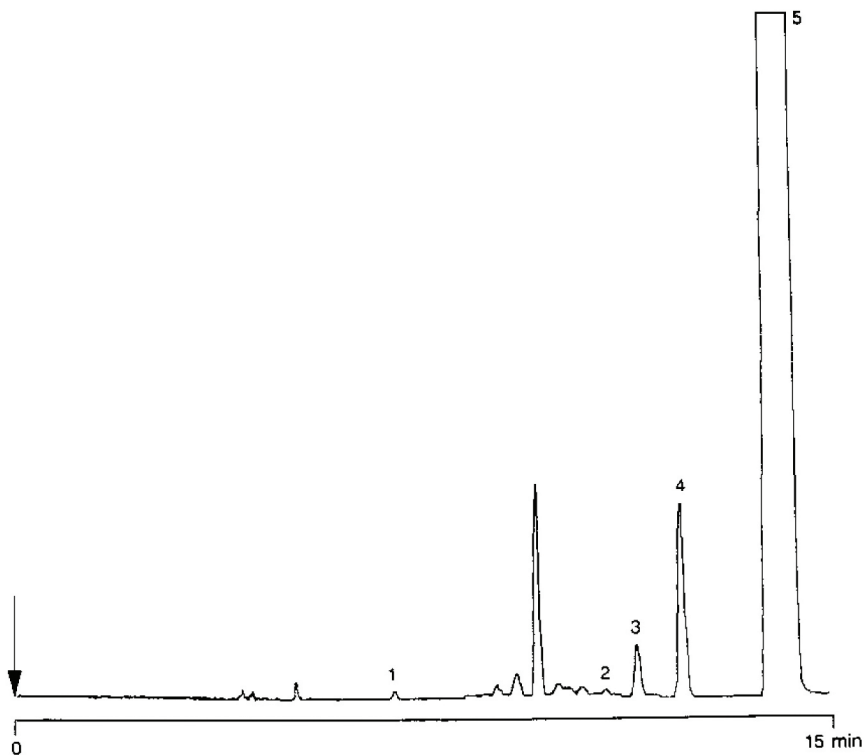
Peak identification

1. toluene
2. ethylbenzene
3. p-xylene
4. m-xylene
5. o-xylene

99+% pure m-xylene



99+% pure o-xylene



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

A00768



Agilent Technologies