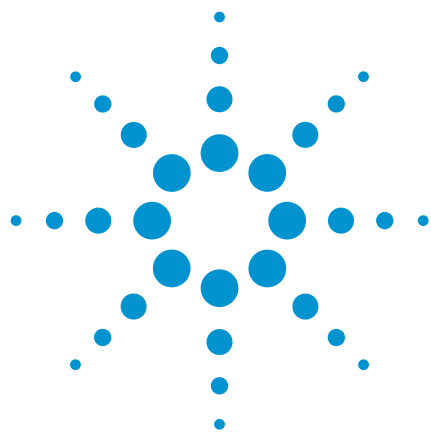




Aromatics, C₆ – C₉

Determination of impurities in styrene

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography using an Agilent CP-Sil 5 CB column separates nine C₆ to C₉ aromatic impurities in styrene in 25 minutes.



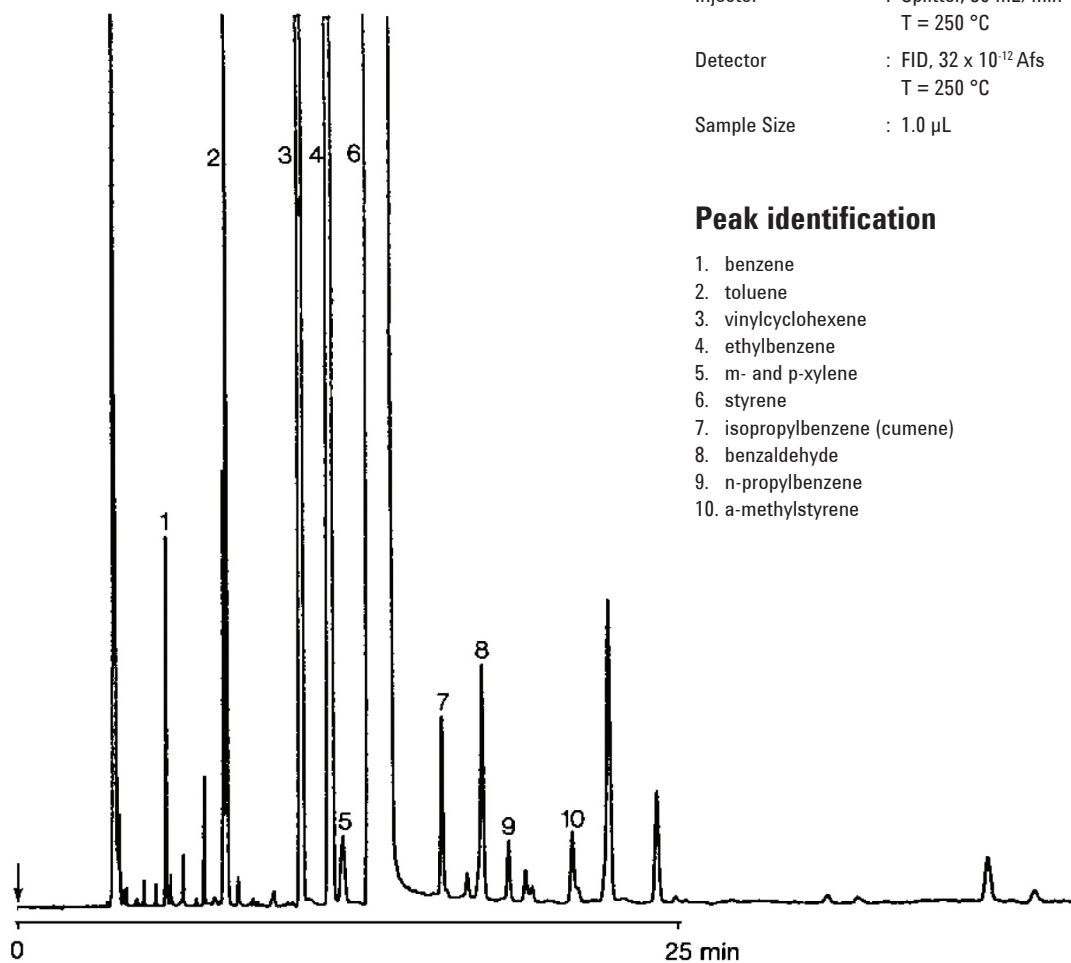
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 5 CB, 0.32 mm x 50 m fused silica
WCOT CP-Sil 5 CB (1.2 μ m) (Part no. CP7770)
Temperature : 100 °C
Carrier Gas : H₂, 100 kPa (1.0 bar, 14.5 psi), 30 cm/s
Injector : Splitter, 30 mL/min
T = 250 °C
Detector : FID, 32 x 10⁻¹² Afs
T = 250 °C
Sample Size : 1.0 μ L

Peak identification

1. benzene
2. toluene
3. vinylcyclohexene
4. ethylbenzene
5. m- and p-xylene
6. styrene
7. isopropylbenzene (cumene)
8. benzaldehyde
9. n-propylbenzene
10. a-methylstyrene



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

A00151



Agilent Technologies