



# Hydrocarbons and aromatics, $C_5 - C_{10}$

## Determination of impurities in dicyclopentadiene

### Application Note

Materials Testing & Research

#### Authors

Agilent Technologies, Inc.

#### Introduction

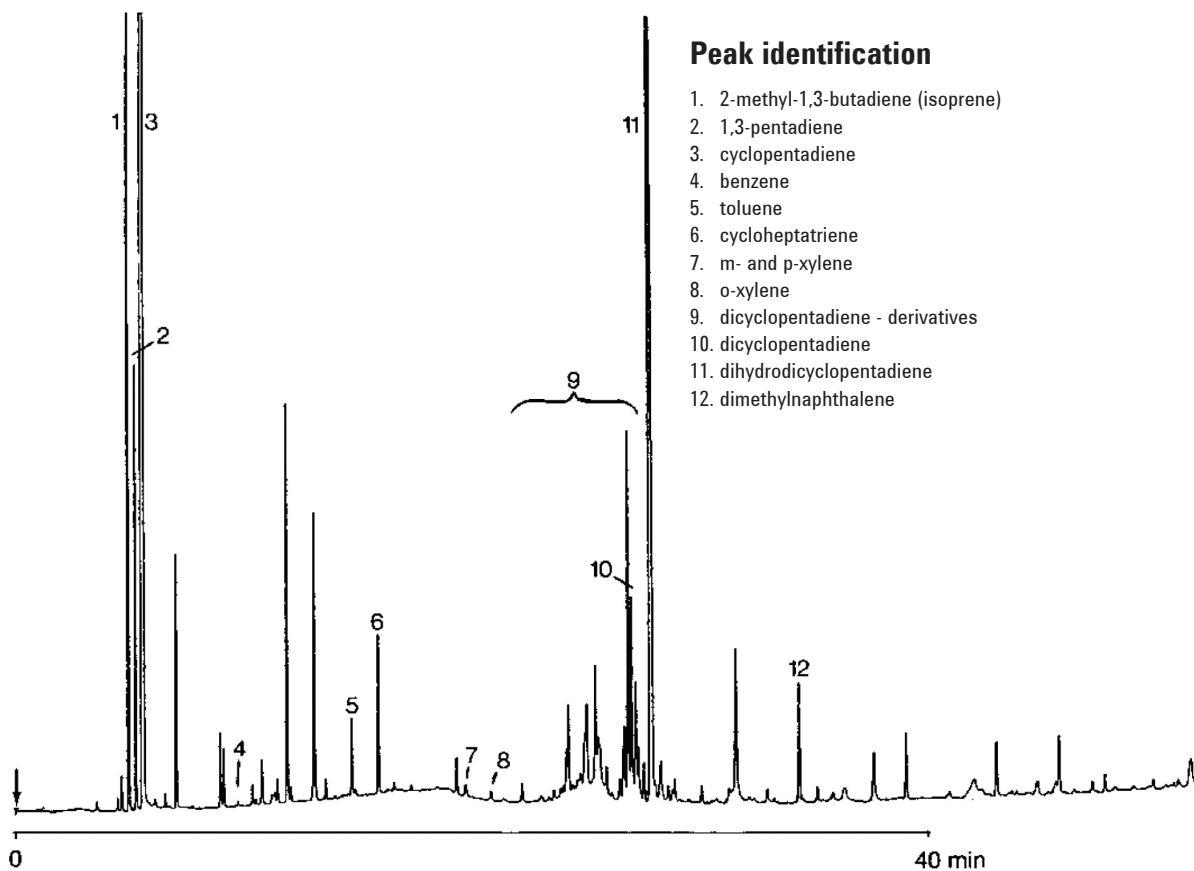
Gas chromatography using an Agilent CP-Sil 5 CB column separates 11  $C_5$  to  $C_{11}$  impurities in dicyclopentadiene in 40 minutes.



**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 $\mu$ m) (Part no. CP7770)
Temperature	: 50 °C (6 min) $\rightarrow$ 275 °C, 5 °C/min
Carrier Gas	: H <sub>2</sub> , 40 kPa (0.4 bar, 5.8 psi), 30 cm/s
Injector	: Splitter, 30 mL/min T = 250 °C
Detector	: FID, 8 x 10 <sup>-12</sup> Afs T = 300 °C
Sample Size	: 0.1 $\mu$ L



## Peak identification

1. 2-methyl-1,3-butadiene (isoprene)
2. 1,3-pentadiene
3. cyclopentadiene
4. benzene
5. toluene
6. cycloheptatriene
7. m- and p-xylene
8. o-xylene
9. dicyclopentadiene - derivatives
10. dicyclopentadiene
11. dihydrodicyclopentadiene
12. dimethylnaphthalene

[www.agilent.com/chem](http://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

A00152



**Agilent Technologies**