

# Hydrocarbons, $C_3 - C_{11}$

## Application Note

Energy & Fuels

### Authors

Agilent Technologies, Inc.

### Introduction

Gas chromatography with an Agilent CP-Sil PONA CB column separates 31 components in unleaded gasoline according to ASTM D5769 in 58 minutes.



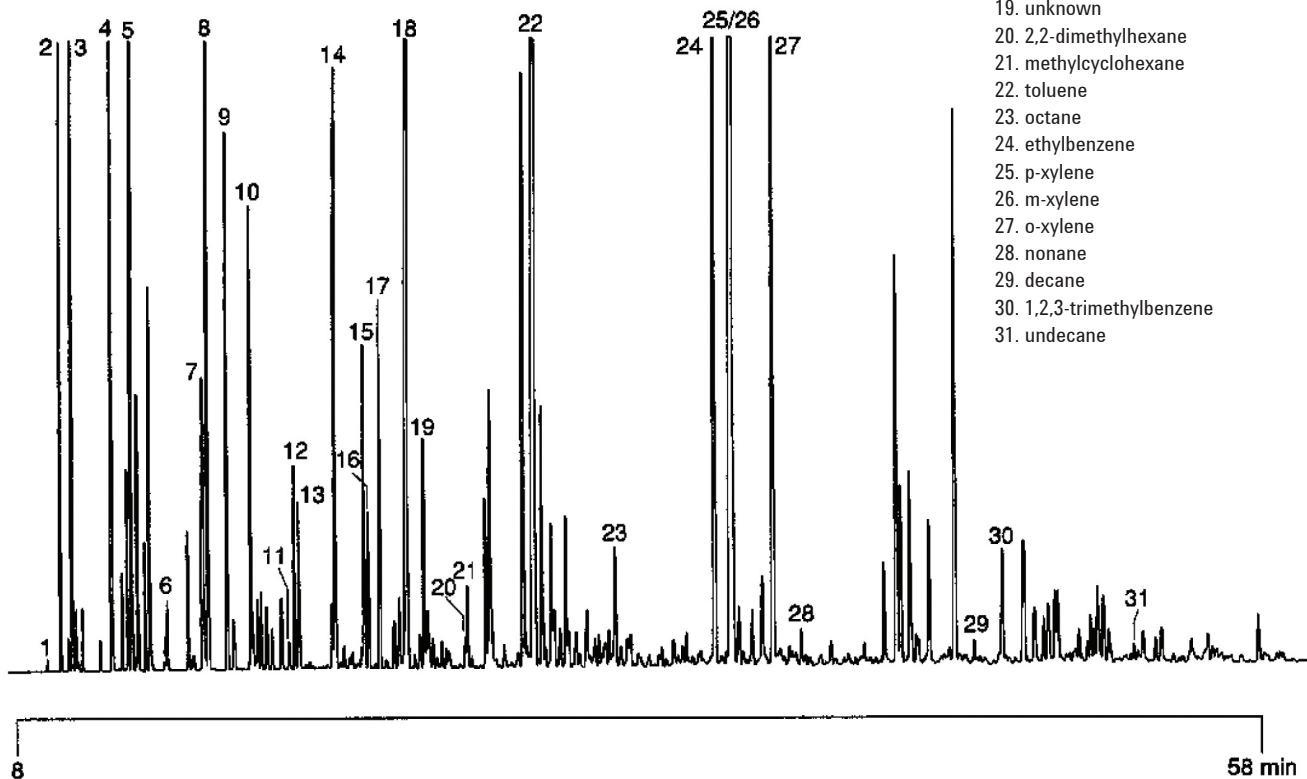
**Agilent Technologies**

## Conditions

Technique : GC-capillary  
Column : Agilent CP-Sil PONA CB, 0.25 mm x 100 m fused silica WCOT (df = 0.5 µm) (Part no. CP7530)  
Temperature : 35 °C (7 min) → 250 °C, 3 °C/min  
Carrier Gas : He, 240 kPa (2.4 bar, 34 psi)  
Injector : Split, 80 mL/min  
T = 275 °C  
Detector : FID  
T = 275 °C  
Sample Size : 0.1 µL

## Peak identification

1. propane
2. isobutane
3. butane
4. 2-methylbutane (isopentane)
5. pentane
6. 2,2-dimethylbutane
7. 2,3-dimethylbutane
8. 2-methylpentane
9. 3-methylpentane
10. hexane
11. 2,2-dimethylpentane
12. methylcyclopentane
13. 2,4-dimethylpentane
14. benzene
15. 2-methylhexane
16. 2,3-dimethylpentane
17. 3-methylhexane
18. tert. amyl methyl ether (TAME)
19. unknown
20. 2,2-dimethylhexane
21. methylcyclohexane
22. toluene
23. octane
24. ethylbenzene
25. p-xylene
26. m-xylene
27. o-xylene
28. nonane
29. decane
30. 1,2,3-trimethylbenzene
31. undecane



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

A01274



**Agilent Technologies**