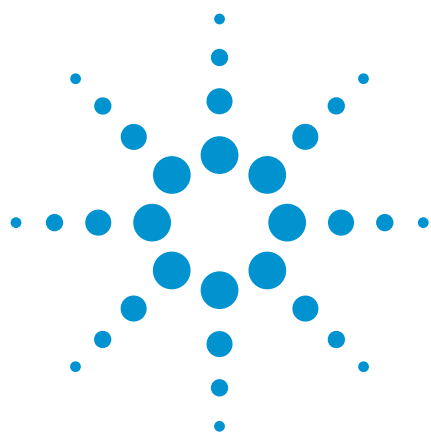




Hydrocarbons, C₁ – C₅

Application Note

Energy & Fuels

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography using an Agilent CP-Al₂O₃/Na₂SO₄ column separates 24 C₁ to C₅ hydrocarbons in 27 minutes.



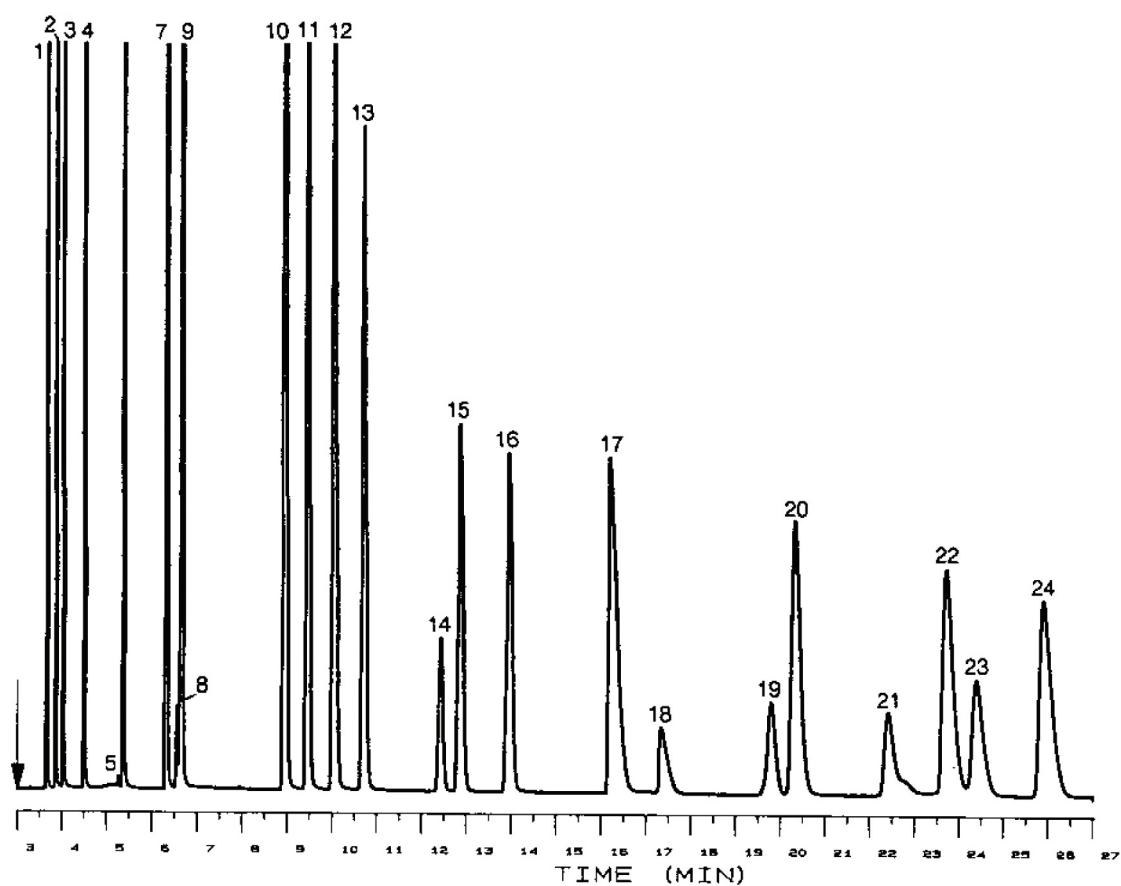
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Al₂O₃/Na₂SO₄, 0.53 mm x 50 m fused silica PLOT CP-Al₂O₃/Na₂SO₄ (df = 10 µm) (Part no. CP7568)
Temperature : 120 °C
Carrier Gas : He, 50 kPa (0.5 bar, 7 psi), 23 cm/s
Injector : Splitter
T = 225 °C
Detector : FID
T = 250 °C

Peak identification

- | | |
|--------------------|-----------------------|
| 1. methane | 13. cis-2-butene |
| 2. ethane | 14. cyclopentane |
| 3. ethylene | 15. isopentane |
| 4. propane | 16. n-pentane |
| 5. cyclopropane | 17. 1,3-butadiene |
| 6. propylene | 18. propyne |
| 7. isobutane | 19. cyclopentene |
| 8. propadiene | 20. trans-2-pentene |
| 9. n-butane | 21. 2-methyl-2-butene |
| 10. trans-2-butene | 22. 1-pentene |
| 11. 1-butene | 23. 2-methyl-1-butene |
| 12. isobutene | 24. cis-2-pentene |



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

A00611



Agilent Technologies