

Hydrocarbons, $C_1 - C_6$

Analysis of light hydrocarbons

Application Note

Energy & Fuels

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography using an Agilent CP- Al_2O_3/KCl column separates 20 light hydrocarbons in 15 minutes.



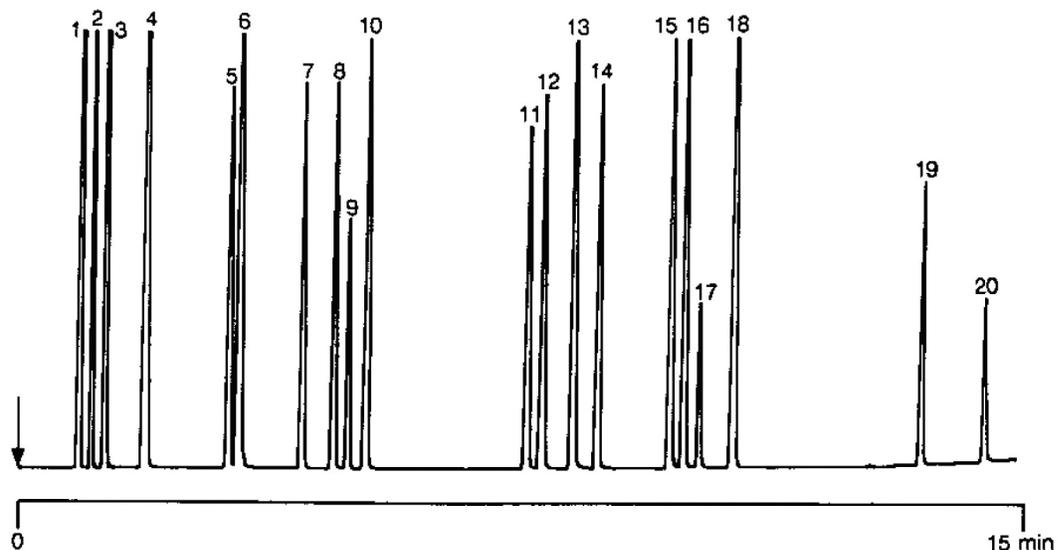
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Al₂O₃/KCl, 0.32 mm x 50 m fused silica
PLOT CP-AL₂O₃/KCl (df = 5 µm) (Part no. CP7515)
Temperature : 70 °C (4 min) → 180 °C, 10 °C/min
Carrier Gas : He, 250 kPa (2.5 bar, 35 psi)
Injector : Splitter, 400 mL/min
Detection : FID
T = 190 °C

Peak identification

1. methane
2. ethane
3. ethylene
4. propane
5. cyclopropane
6. propylene
7. acetylene
8. isobutane
9. propadiene
10. butane
11. trans-2-butene
12. 1-butene
13. isobutene
14. cis-2-butene
15. isopentane
16. propyne
17. pentane
18. 1,3-butadiene
19. ethylacetylene
20. hexane



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

A00677



Agilent Technologies