



Hydrocarbons, $C_1 - C_7$

Analysis of hydrocarbons in natural gas

Application Note

Energy & Fuels

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography using an Agilent CP-Al₂O₃/KCl column separates 23 C_1 to C_7 hydrocarbons in natural gas in 50 minutes.



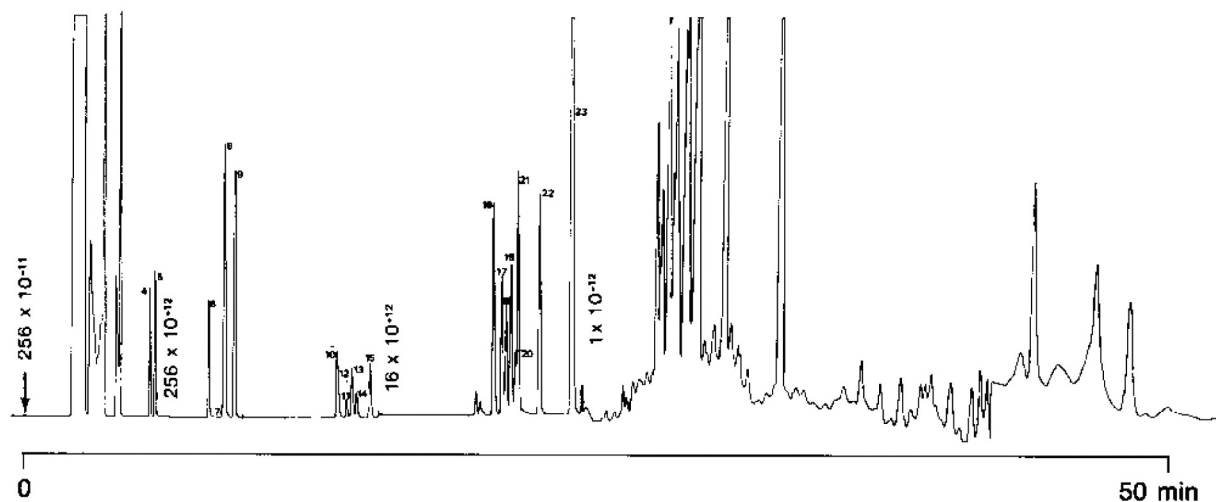
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Al₂O₃/KCl, 0.32 mm x 50 m fused silica
PLOT Al₂O₃/KCl (Part no. CP7515)
Temperature : 100 °C → 200 °C (3 °C/min)
Carrier Gas : N₂, 120 kPa (1.2 bar, 17 psi)
Injector : Splitless (30 s), sample trapped at injector side at
-196 °C
Detector : FID 256 x 10⁻¹¹ Afs
256 x 10⁻¹² Afs after 5.5 min
16 x 10⁻¹² Afs after 14 min
1 x 10⁻¹² Afs after 24 min

Peak identification

1. methane
2. ethane
3. propane
4. isobutane
5. butane
6. 2,2-dimethylpropane
7. cyclopentane
8. 2-methylbutane
9. pentane
10. 2,2-dimethylbutane
11. cyclohexane
12. 2,3-dimethylbutane
13. 2-methylbutane
14. 3-methylpentane
15. hexane
16. 2,2-dimethylpentane and 2,4-dimethylpentane
17. methylcyclohexane
18. 2,2,3-trimethylbutane
19. 3,3-dimethylpentane
20. 2,3-dimethylpentane
21. 2-methylhexane, 3-methylhexane and 3-ethylpentane
22. heptane
23. benzene



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

A00271



Agilent Technologies