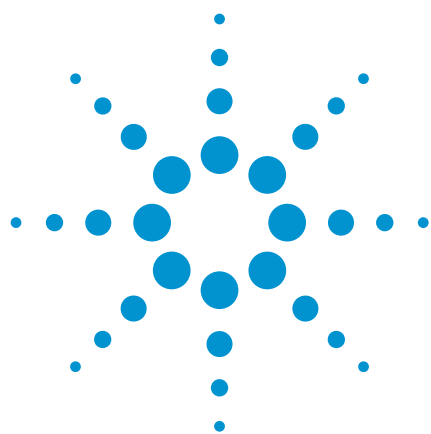




Light hydrocarbons, $C_1 - C_{10}$

Analysis of natural gas on a wide-bore column

Application Note

Energy & Fuels

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography using an Agilent CP-Al₂O₃/KCl column separates ten C_1 to C_{10} light hydrocarbons in natural gas in 20 minutes.



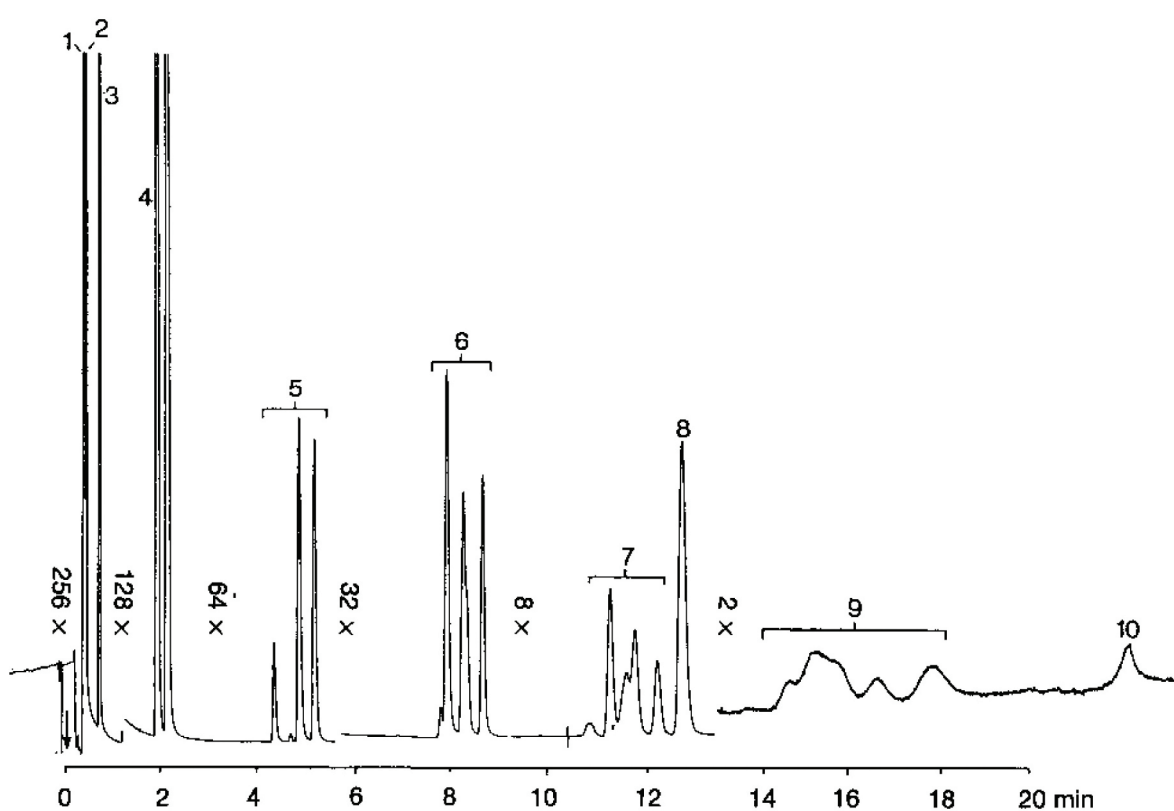
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Al₂O₃/KCl, 0.53 mm x 10 m fused silica
Al₂O₃/KCl PLOT tailor made for the analysis of C₁-C₁₀
hydrocarbons (Part no. CP7516)
Temperature : 30 °C → 200 °C, 10 °C/min
Carrier Gas : H₂, 13 kPa (0.13 bar)
6 mL/min
Injector : Direct Injector
Detector : FID

Peak identification

1. methane
2. ethane
3. propane
4. C₄
5. C₅
6. C₆
7. C₇
8. benzene
9. C₈
10. toluene



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

A00225



Agilent Technologies