



## Hydrocarbons, C<sub>1</sub> – C<sub>6</sub>, other gases

Analysis of hydrocarbons C<sub>1</sub> - C<sub>6</sub>, CO<sub>2</sub>,  
H<sub>2</sub>S and water

### Application Note

Energy & Fuels

#### Authors

Agilent Technologies, Inc.

#### Introduction

Gas chromatography using an Agilent PoraPLOT Q column separates 11 C<sub>1</sub> to C<sub>6</sub> hydrocarbons and other gases in six minutes.



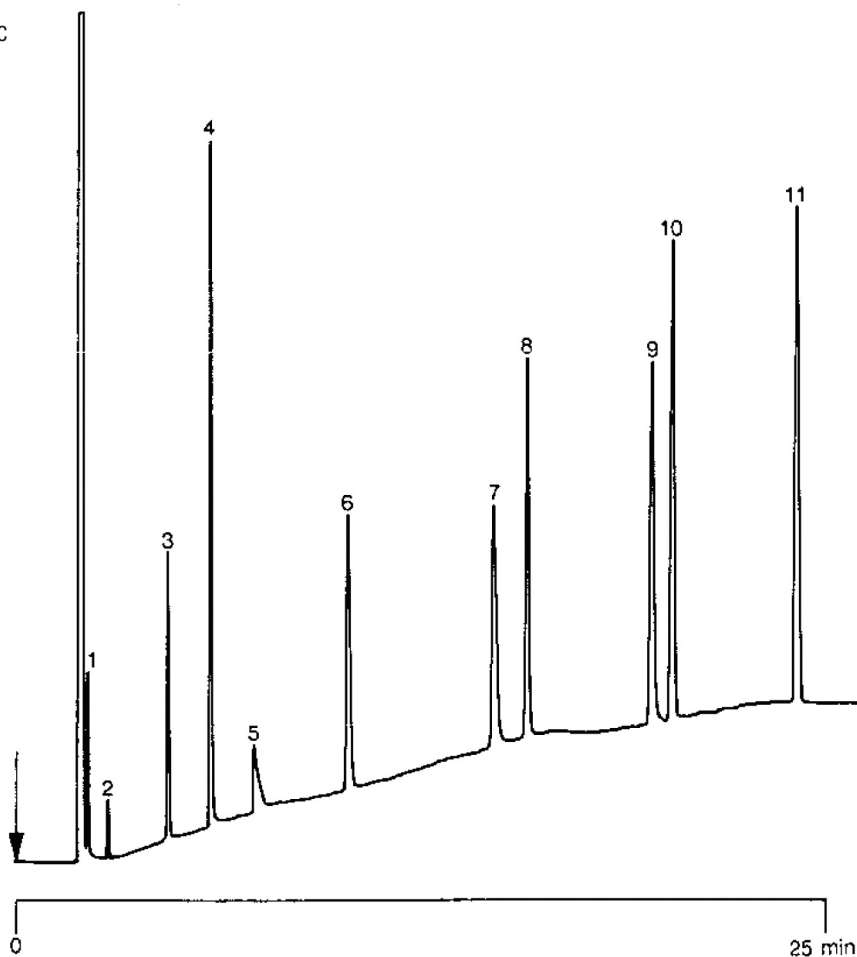
**Agilent Technologies**

## Conditions

Technique : GC-capillary  
Column : Agilent PoraPLOT Q, 0.53 mm x 25 m fused silica  
PLOT PoraPLOT Q (df = 20  $\mu$ m) (Part no. CP7554)  
Temperature : 40 °C  $\rightarrow$  200 °C, 10 °C/min  
Carrier Gas : He, 120 kPa (1.2 bar, 17 psi), 7 mL/min  
Injector : Splitter  
T = 200 °C  
Detector :  $\mu$ -TCD, x4  
T = 220 °C, filaments T = 280 °C  
Sample Size : 0.5 mL  
Courtesy : GL-Sciences, Japan

## Peak identification

1. methane
2. carbon dioxide
3. ethane
4. hydrogen sulfide
5. water
6. propane
7. isobutane
8. n-butane
9. isopentane
10. n-pentane
11. n-hexane



[www.agilent.com/chem](http://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

A00607



**Agilent Technologies**