



# Hydrocarbons, C<sub>1</sub> – C<sub>6</sub>, and solvents

## Application Note

Energy & Fuels

### Authors

Agilent Technologies, Inc.

### Introduction

Gas chromatography with an Agilent PoraPLOT U column separates 11 C<sub>1</sub> to C<sub>6</sub> hydrocarbons and solvents in less than five minutes.



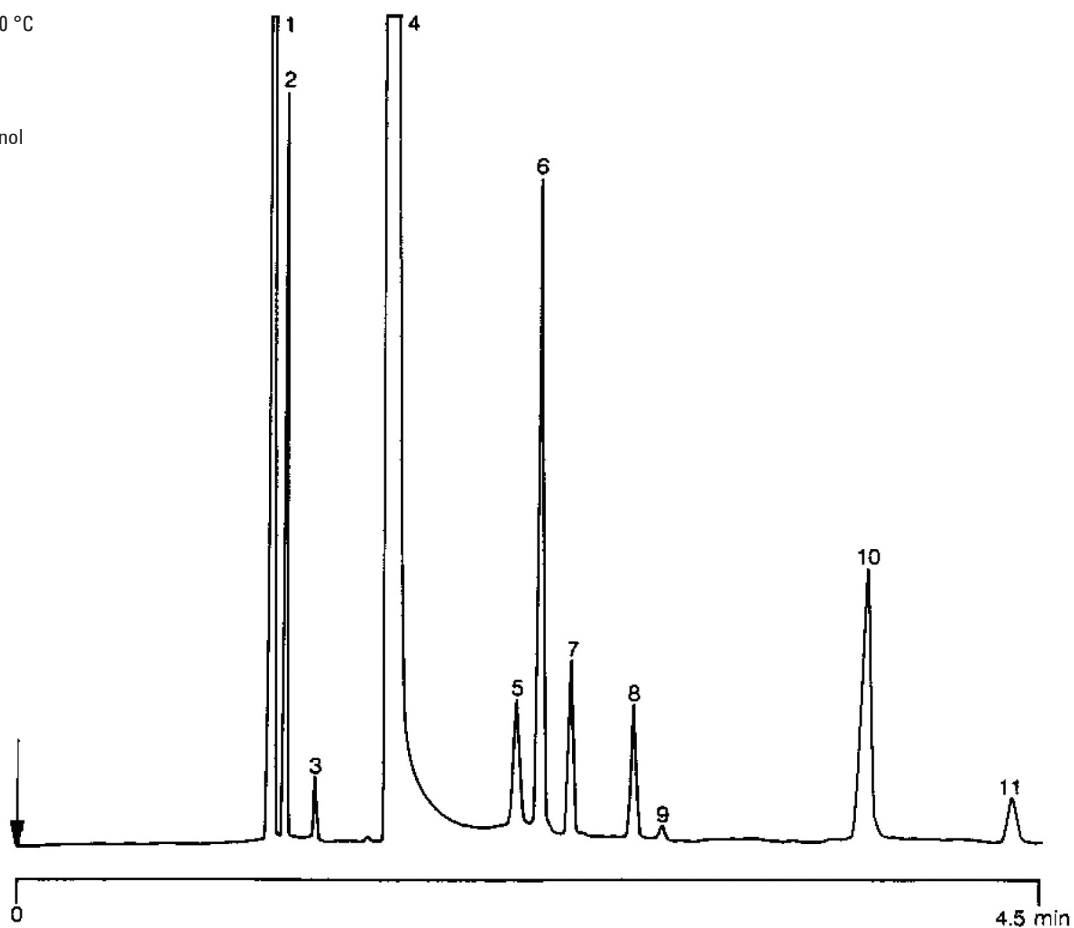
**Agilent Technologies**

## Conditions

Technique : GC-capillary  
Column : Agilent PoraPLOT U, 0.32 mm x 25 m fused silica  
PLOT PoraPLOT U (df = 10 µm) (Part no. CP7581)  
Temperature : 175 °C  
Carrier Gas : H<sub>2</sub>, 45 kPa (0.45 bar, 6.4 psi)  
Injector : Splitter, 100 mL/min;  
T = 200 °C  
Detector : FID, 8;  
T = 250 °C  
Sample Size : 0.1 µL  
Concentration Range : 0.1%  
Solvent Sample : methanol

## Peak identification

1. methane
2. ethane
3. propane
4. methanol + butane
5. ethanol
6. acetone
7. diethyl ether
8. pentane
9. unknown
10. ethyl acetate
11. 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

A00670



**Agilent Technologies**