



# Hydrocarbons, $C_1 - C_3$

## Natural gas analysis

### Application Note

Energy & Fuels

#### Authors

Agilent Technologies, Inc.

#### Introduction

Gas chromatography with an Agilent PoraPLOT Q column separates six components of natural gas in three minutes.



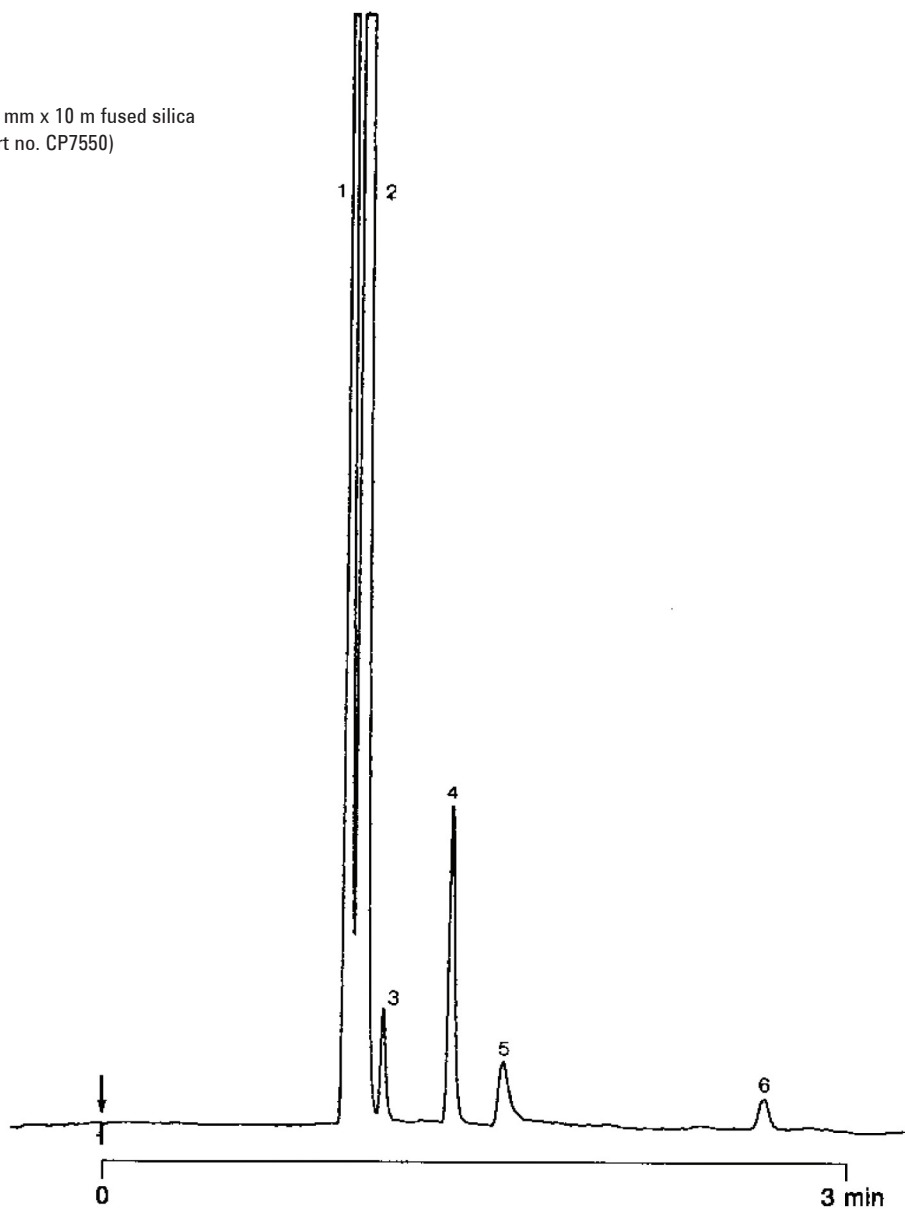
**Agilent Technologies**

## Conditions

Technique : GC-capillary  
Column : Agilent PoraPLOT Q, 0.32 mm x 10 m fused silica  
PoraPLOT Q (10  $\mu$ m) (Part no. CP7550)  
Temperature : 30  $^{\circ}$ C  
Carrier Gas : H<sub>2</sub>, 50 kPa (0.5 bar, 7 psi)  
Injector : Splitter  
Detector :  $\mu$  TCD

## Peak identification

1. nitrogen
2. methane
3. carbon dioxide
4. ethane
5. water
6. propane



[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

A00287



**Agilent Technologies**