



# Hydrocarbons, $C_1 - C_3$

## Application Note

Energy & Fuels

### Authors

Agilent Technologies, Inc.

### Introduction

Gas chromatography with an Agilent PoraPLOT Q column separates four  $C_1$  to  $C_3$  hydrocarbons in five minutes.



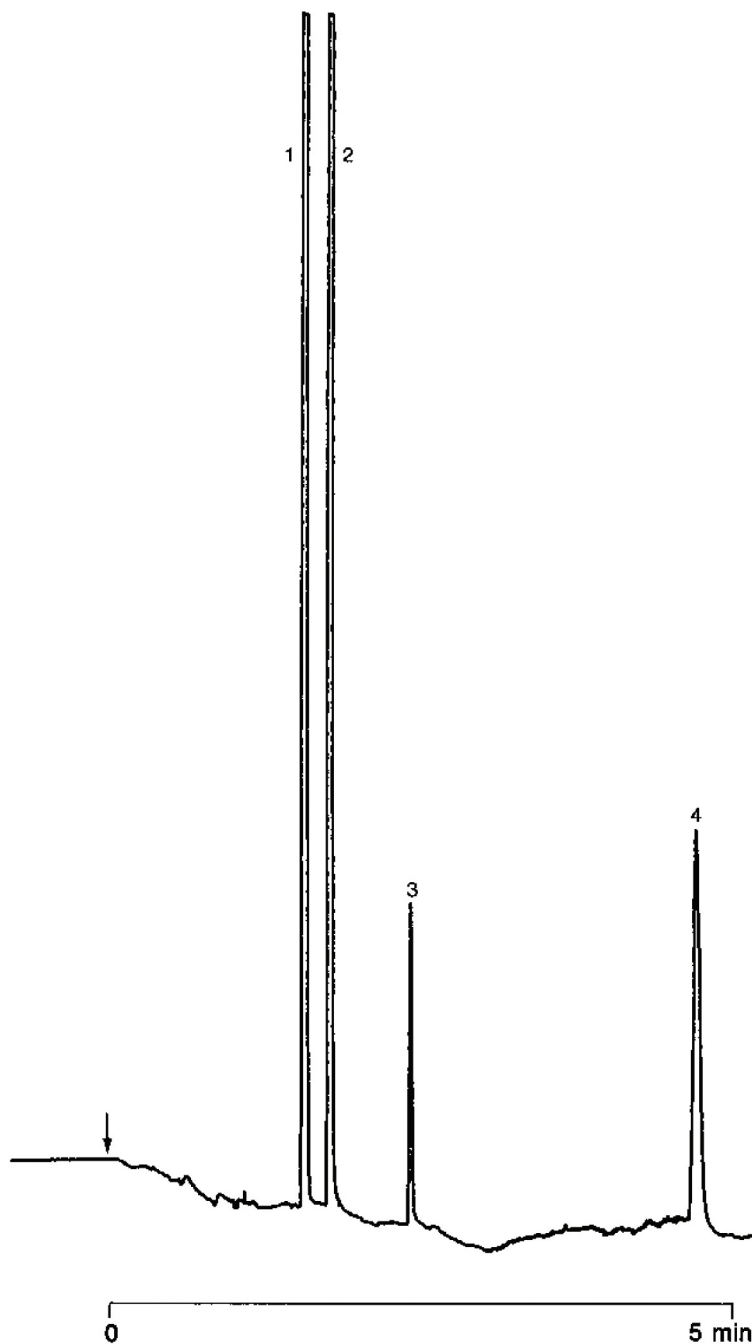
**Agilent Technologies**

## Conditions

Technique : GC-capillary  
Column : Agilent PoraPLOT Q, 0.32 mm x 25 m fused silica PoraPLOT Q (10  $\mu$ m) (Part no. CP7551)  
Temperature : 25  $^{\circ}$ C  
Carrier Gas :  $H_2$ , 100 kPa (1.0 bar, 14.5 psi)  
Injector : Splitter  
Detector :  $\mu$  TCD

## Peak identification

1. nitrogen
2. methane
3. ethane
4. 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

A00290



**Agilent Technologies**