



Hydrocarbons, $C_1 - C_2$

Application Note

Energy & Fuels

Authors

Agilent Technologies, Inc.

Introduction

Fast gas chromatography with an Agilent PoraPLOT U column separates four C_1 to C_4 hydrocarbons in two minutes at 30 °C and 1 minute at 60 °C.



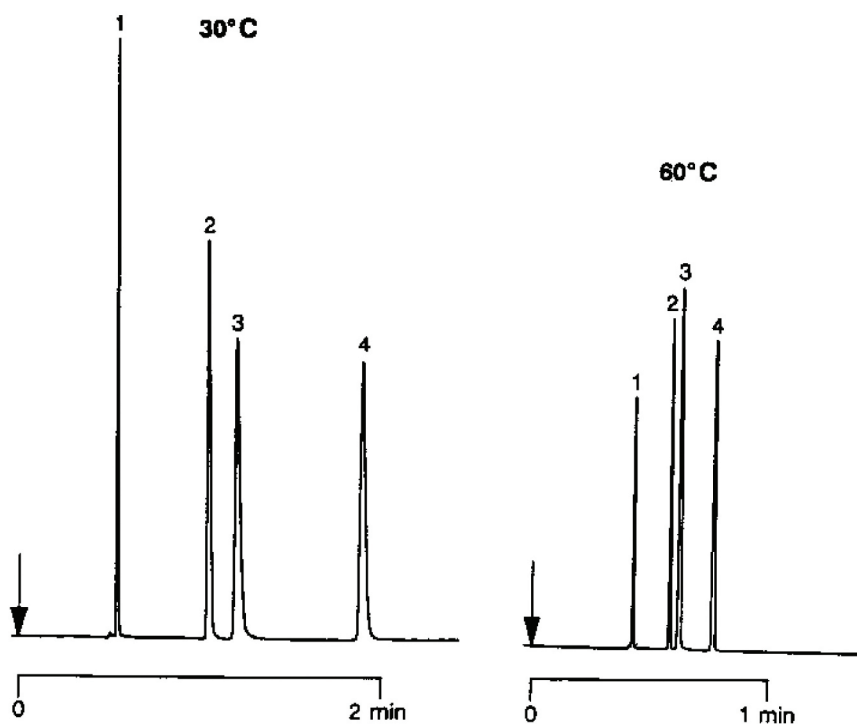
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent PoraPLOT U, 0.53 mm x 25 m fused silica
PLOT PoraPLOT U (df = 20 μ m) (Part no. CP7584)
Temperature : A 30 °C; B 60 °C
Carrier Gas : H₂
Injector : Split, 100 mL/min
Detector : FID, 64 X 10⁻¹² Afs
Sample Size : 1% in N₂

Peak identification

1. methane
2. ethylene
3. ethane
4. acetylene



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

A00685



Agilent Technologies