



Hydrocarbons, C₁ – C₃, and polar volatiles

Application Note

Energy & Fuels

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT U column separates six C₁ to C₃ hydrocarbons and polar volatiles in five minutes.



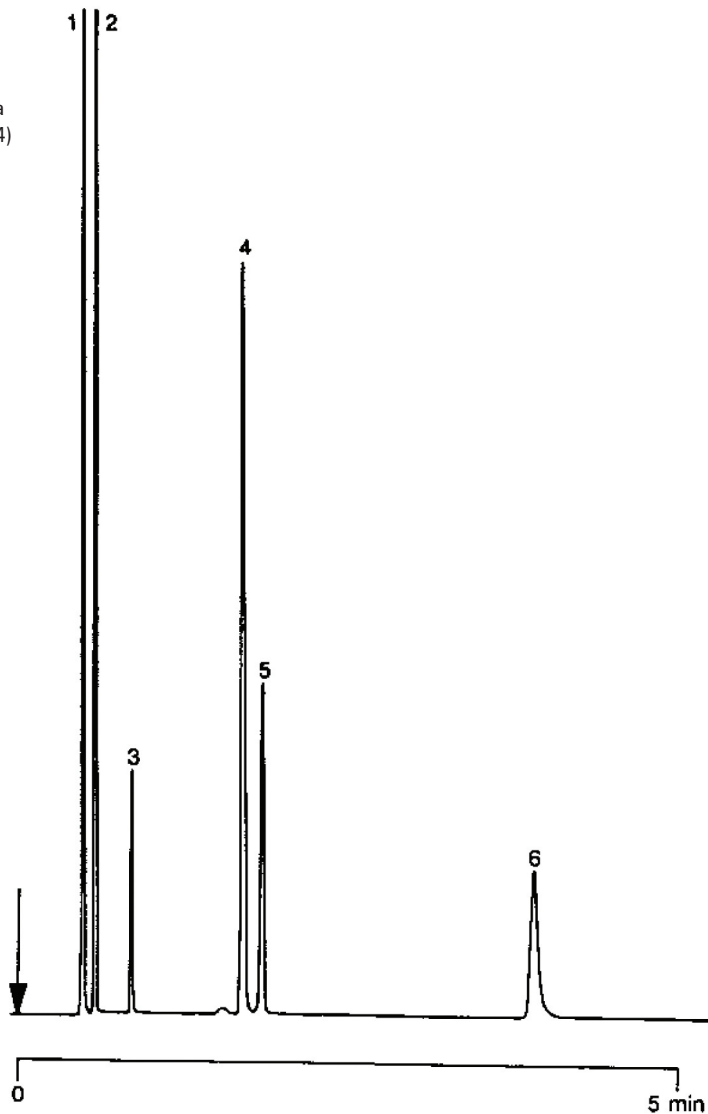
Agilent Technologies

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 : 115 °C
Carrier Gas : H₂, 20 kPa (0.2 bar, 2.9 psi)
Injector : Split, 100 mL/min
Detector : FId, 32 x 10⁻¹² Afs
Sample Size : 100 ppm in H₂

Peak identification

1. methane
2. ethane
3. propane
4. methanol
5. acetaldehyde
6. ethanol



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

A00686



Agilent Technologies