



Hydrocarbons, C₁ – C₃

Application Note

Energy & Fuels

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT S column separates eight C₁ to C₃ hydrocarbons in ten minutes.



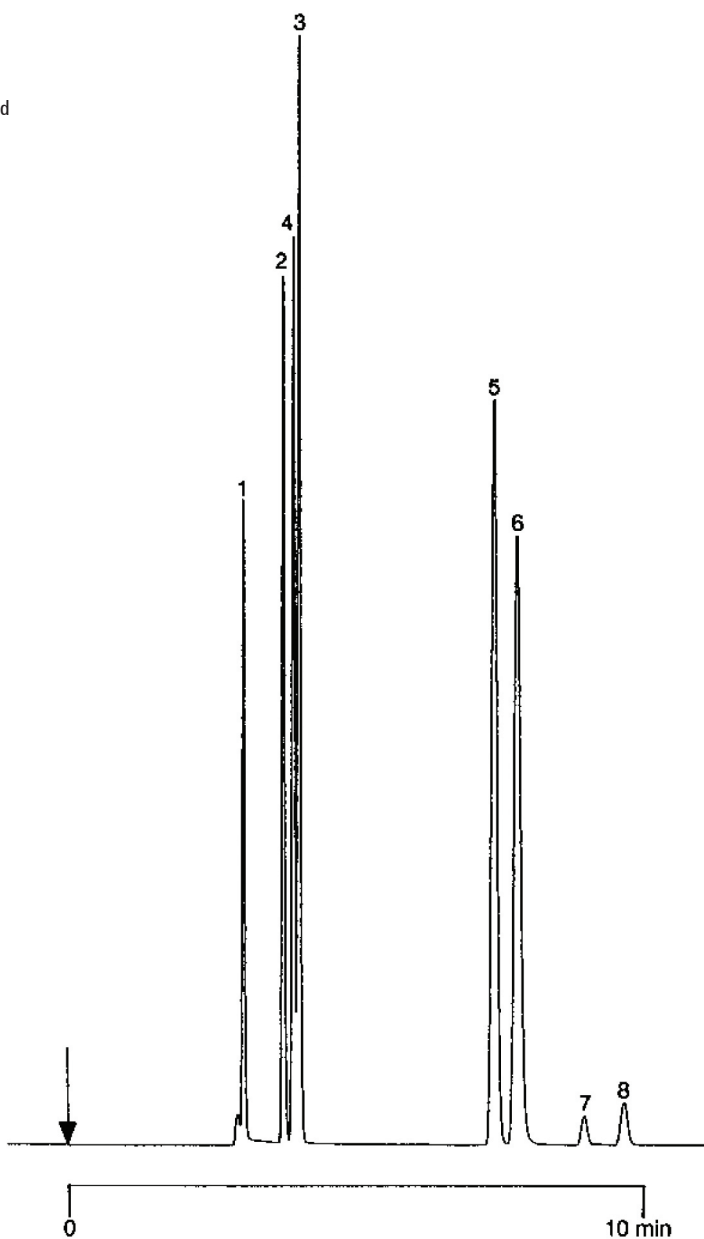
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent PoraPLOT S, 0.53 mm x 25 m (ID x L) fused silica PoraPLOT S (df = 20 μ m) (Part no. CP7574)
Temperature : 80 °C
Carrier Gas : H₂, 30 kPa (0.3 bar, 4.2 psi)
Injector : Split, 100 mL/min
Detector : FID, 64 x 10⁻¹² Afs
Sample : 0.1 - 1% in N₂

Peak identification

1. methane
2. ethylene
3. ethane
4. acetylene
5. propylene
6. propane
7. propadiene
8. propyne



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

A00452



Agilent Technologies