



Hydrocarbons, $C_1 - C_4$

Analysis of light hydrocarbons on a 150 μm column

Application Note

Energy & Fuels

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography using an Agilent CP-Sil 5 CB column separates 13 light hydrocarbons in five minutes.



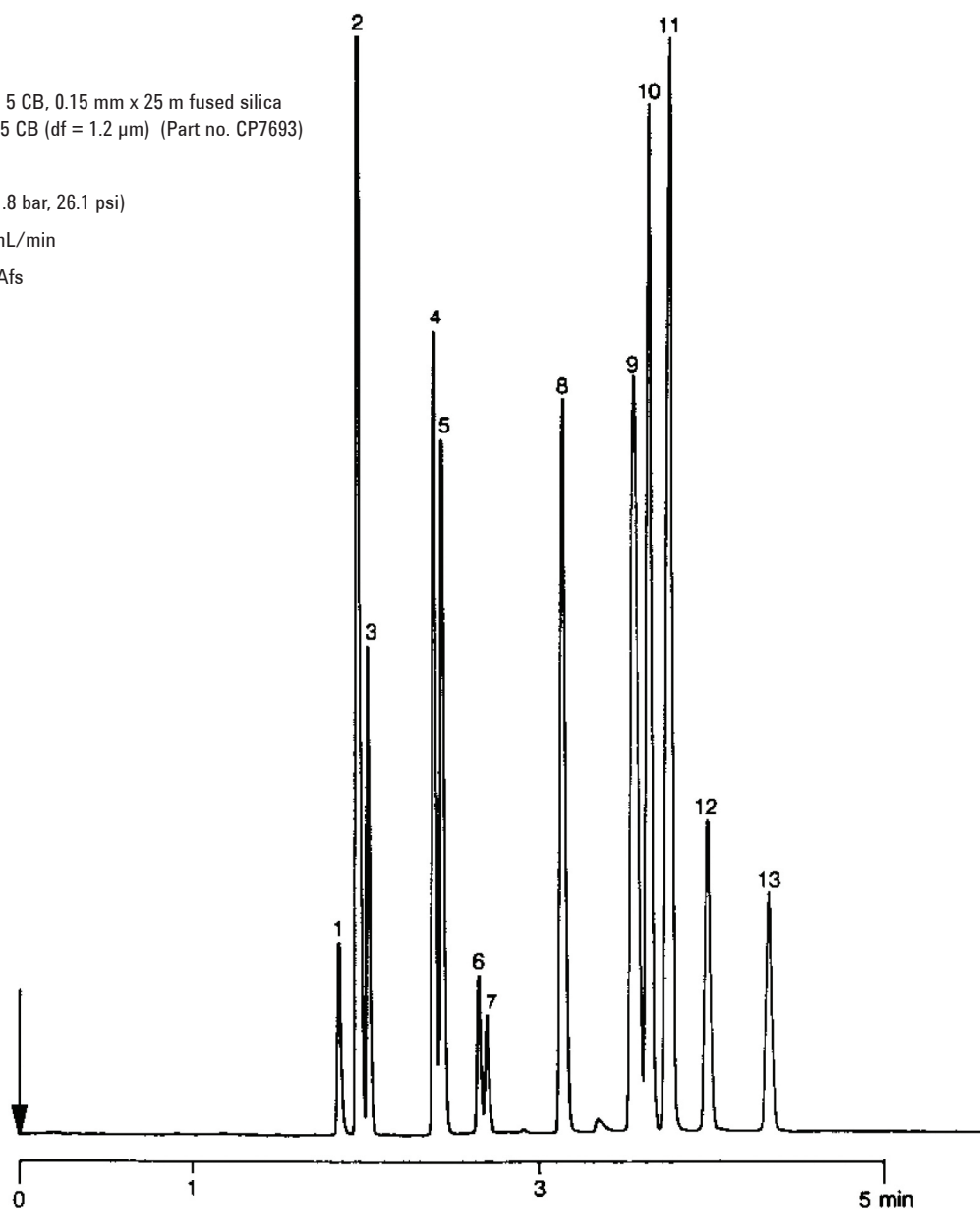
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 5 CB, 0.15 mm x 25 m fused silica
WCOT CP-Sil 5 CB (df = 1.2 μ m) (Part no. CP7693)
Temperature : 25 $^{\circ}$ C
Carrier Gas : H₂, 180 kPa (1.8 bar, 26.1 psi)
Injector : Splitter, 150 mL/min
Detection : FID, 4 X 10⁻¹² Afs

Peak identification

1. methane
2. ethane
3. ethane, acetylene
4. propene
5. propane
6. propadiene
7. propyne
8. isobutane
9. isobutene, 1-butene
10. 1,3-butadiene
11. n-butane
12. trans-2-butene
13. cis-2-butene



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

A00678



Agilent Technologies