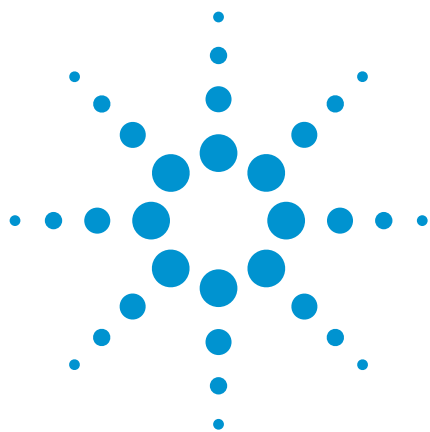




Hydrocarbons, $C_1 - C_6$

Analysis of impurities in propylene

Application Note

Energy & Fuels

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Al₂O₃/Na₂SO₄ column identifies 15 C₁ to C₆ hydrocarbon impurities in propylene in 15 minutes.



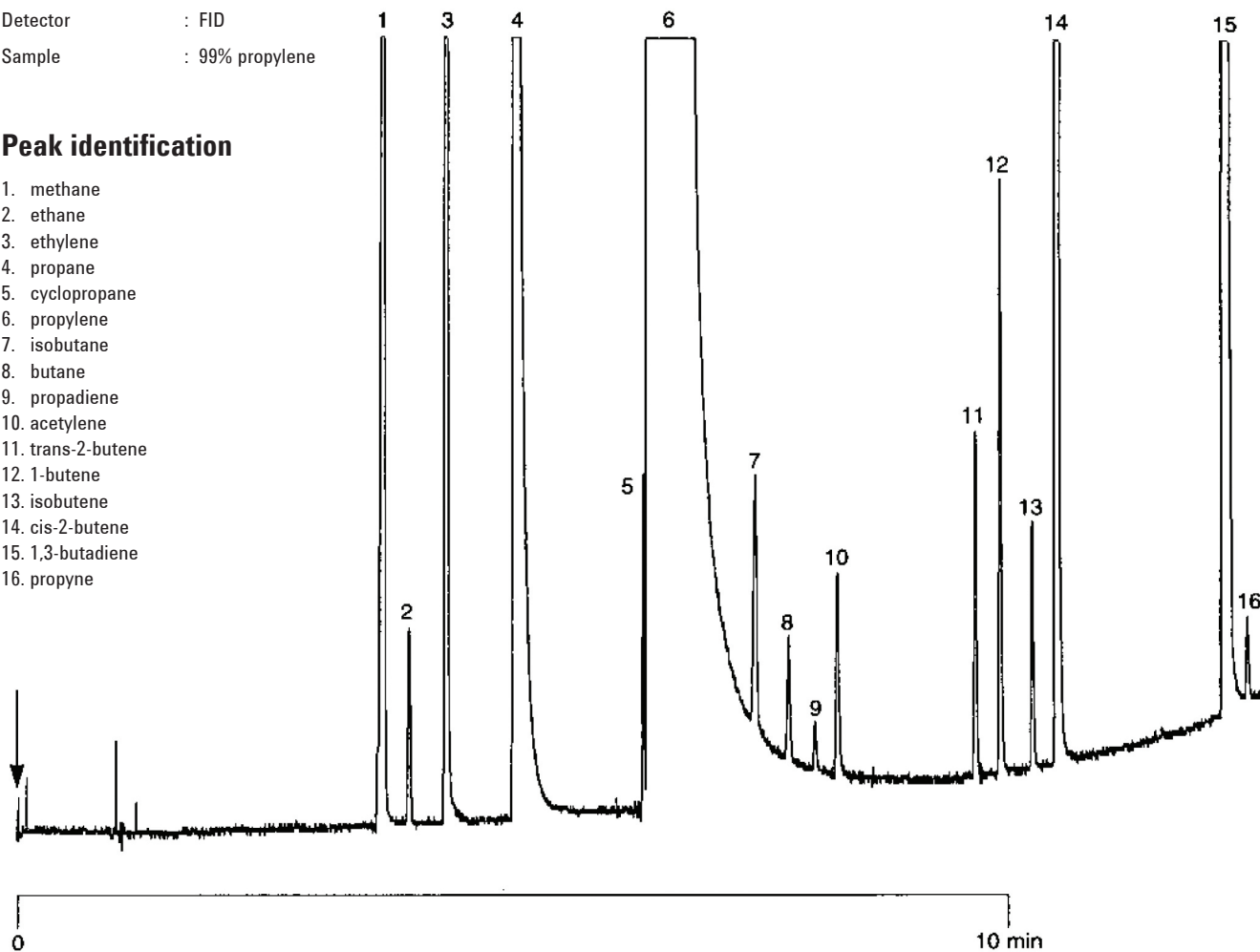
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Al₂O₃/Na₂SO₄, 0.25 mm x 50 m fused silica PLOT CP-Al₂O₃/Na₂SO₄ (df= 4.0 µm) (Part no. CP7587)
Temperature : 50 °C (1 min) → 200 °C, 10 °C/min
Carrier Gas : He, 220 kPa (2.2 bar, 31 psi)
Injector : Split, 30 mL/min
Detector : FID
Sample : 99% propylene

Peak identification

1. methane
2. ethane
3. ethylene
4. propane
5. cyclopropane
6. propylene
7. isobutane
8. butane
9. propadiene
10. acetylene
11. trans-2-butene
12. 1-butene
13. isobutene
14. cis-2-butene
15. 1,3-butadiene
16. 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

A01098



Agilent Technologies