



Hydrocarbons, $C_1 - C_6$

Analysis of impurities in cis-2-butene

Application Note

Energy & Fuels

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Al₂O₃/KCl column resolves 21 C₁ to C₆ hydrocarbon impurities in cis-2-butene in 20 minutes.



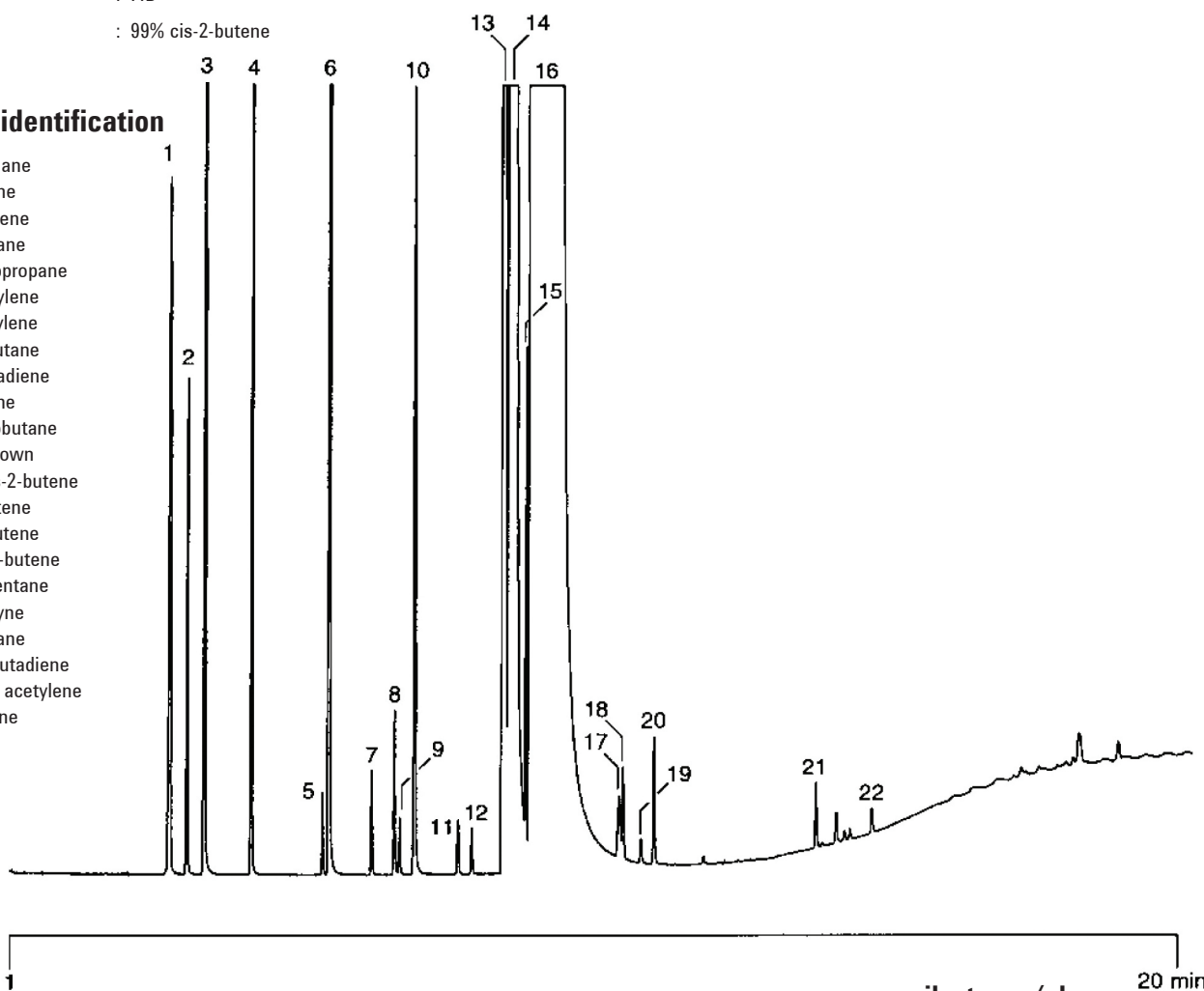
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Al₂O₃/KCl, 0.25 mm x 50 m fused silica
PLOT CP-Al₂O₃/KCl (df= 4.0 μm)
(Part no. CP7577)
Temperature : 50 °C (1 min) → 200 °C, 10 °C/min
Carrier Gas : He, 220 kPa (2.2 bar, 31 psi)
Injector : Split, 20 mL/min
Detector : FID
Sample : 99% cis-2-butene

Peak identification

1. methane
2. ethane
3. ethylene
4. propane
5. cyclopropane
6. propylene
7. acetylene
8. isobutane
9. propadiene
10. butane
11. cyclobutane
12. unknown
13. trans-2-butene
14. 1-butene
15. isobutene
16. cis-2-butene
17. isopentane
18. propyne
19. pentane
20. 1,3-butadiene
21. ethyl acetylene
22. hexane



www.agilent.com/chem

20 min

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

A01099



Agilent Technologies