



Alcohols, C₁ – C₆

Separation of volatile alcohols in 1-hexanol

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography using an Agilent CP-Sil 5 CB column separates ten C₁ to C₆ volatile alcohols in 1-hexanol in 15 minutes.



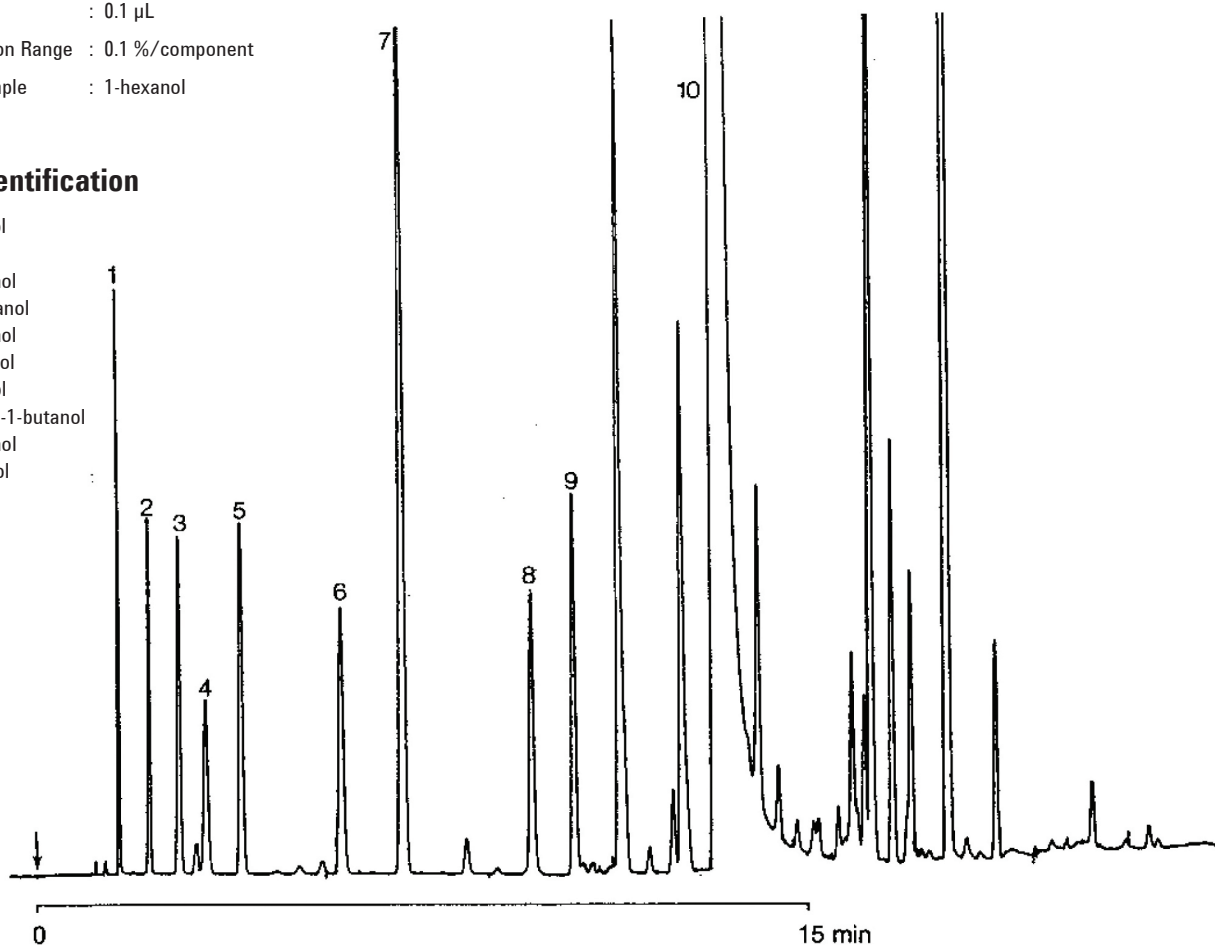
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 5 CB, 0.53 mm x 50 m fused silica
WCOT CP-Sil 5 CB (5.0 μ m) (Part no. CP7685)
Temperature : 200 °C
Carrier Gas : H₂, 50 kPa (0.5 bar, 7.25 psi), 25 cm/s
Injector : Splitter, 55 mL/min
T = 240 °C
Detector : FID, 16 x 10⁻¹² Afs
T = 250 °C
Sample Size : 0.1 μ L
Concentration Range : 0.1 %/component
Solvent Sample : 1-hexanol

Peak identification

1. methanol
2. ethanol
3. 2-propanol
4. tert. butanol
5. 1-propanol
6. isobutanol
7. 1-butanol
8. 2-methyl-1-butanol
9. 1-pentanol
10. 1-hexanol



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

A00162



Agilent Technologies