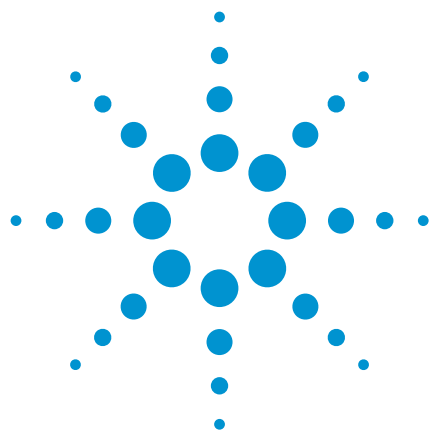




Volatile alcohols and ketones, $C_2 - C_6$

Determination of impurities in 1-propanol

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Wax 57 CB column determines 13 C_2 to C_6 volatile alcohols and ketones impurities in 1-propanol in 25 minutes.



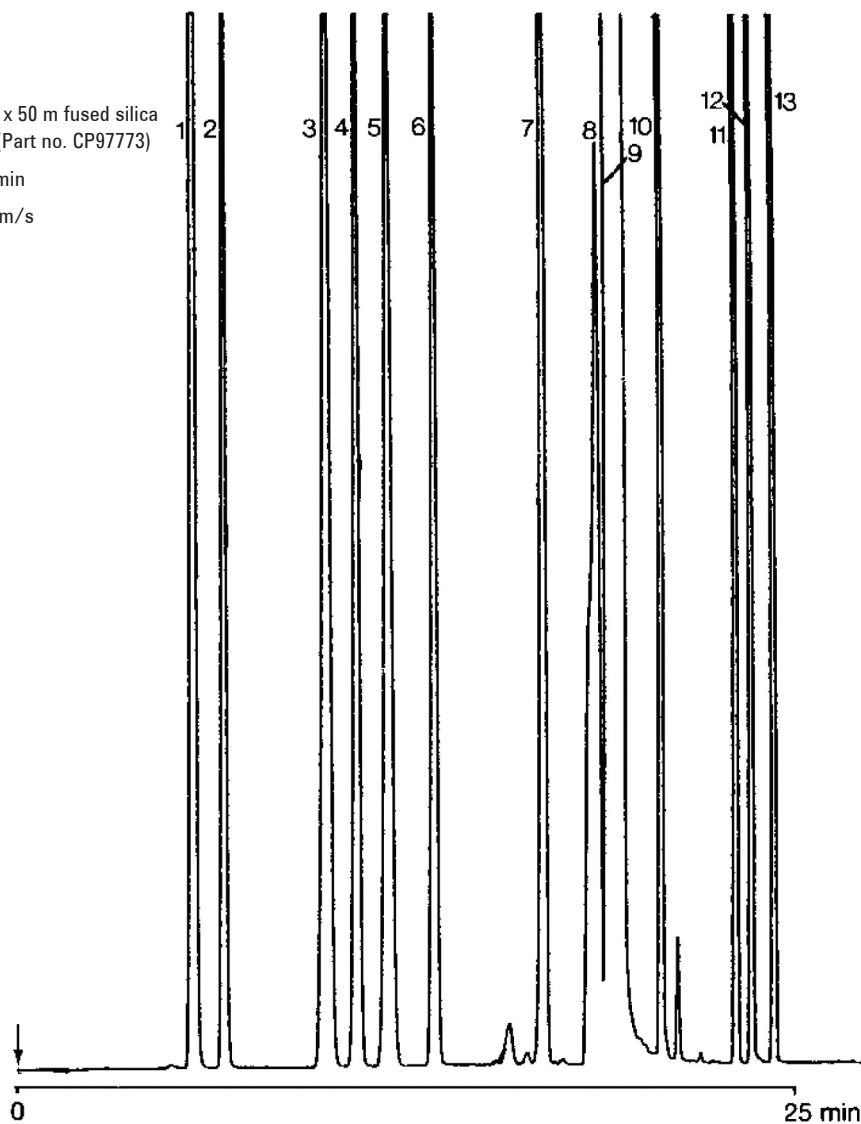
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Wax 57 CB, 0.32 mm x 50 m fused silica
WCOT CP-Wax 57 CB (1.2 μ m) (Part no. CP97773)
Temperature : 40 °C (10 min) $\rightarrow$ 150 °C, 5 °C/min
Carrier Gas : H₂, 40 kPa (0.4 bar, 5.8 psi), 30 cm/s
Injector : Splitter, 50 mL/min
T = 200 °C
Detector : FID, 16 x 10⁻¹² Afs
T = 250 °C
Sample Size : 0.1 μ L
Solvent Sample : 1-propanol

Peak identification

1. n-octane
2. acetone
3. n-nonane
4. butanone
5. tert. butanol
6. 2-propanol
7. methylisobutylketone
8. 2-butanol
9. 1-propanol
10. 2-hexanone
11. mesityloxide
12. 1-butanol
13. methylisobutylcarbinol



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

A00150



Agilent Technologies