



Alcohols and acetates

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Wax for Volatile Amines and Diamines column separates 17 alcohols and acetates in ten minutes.



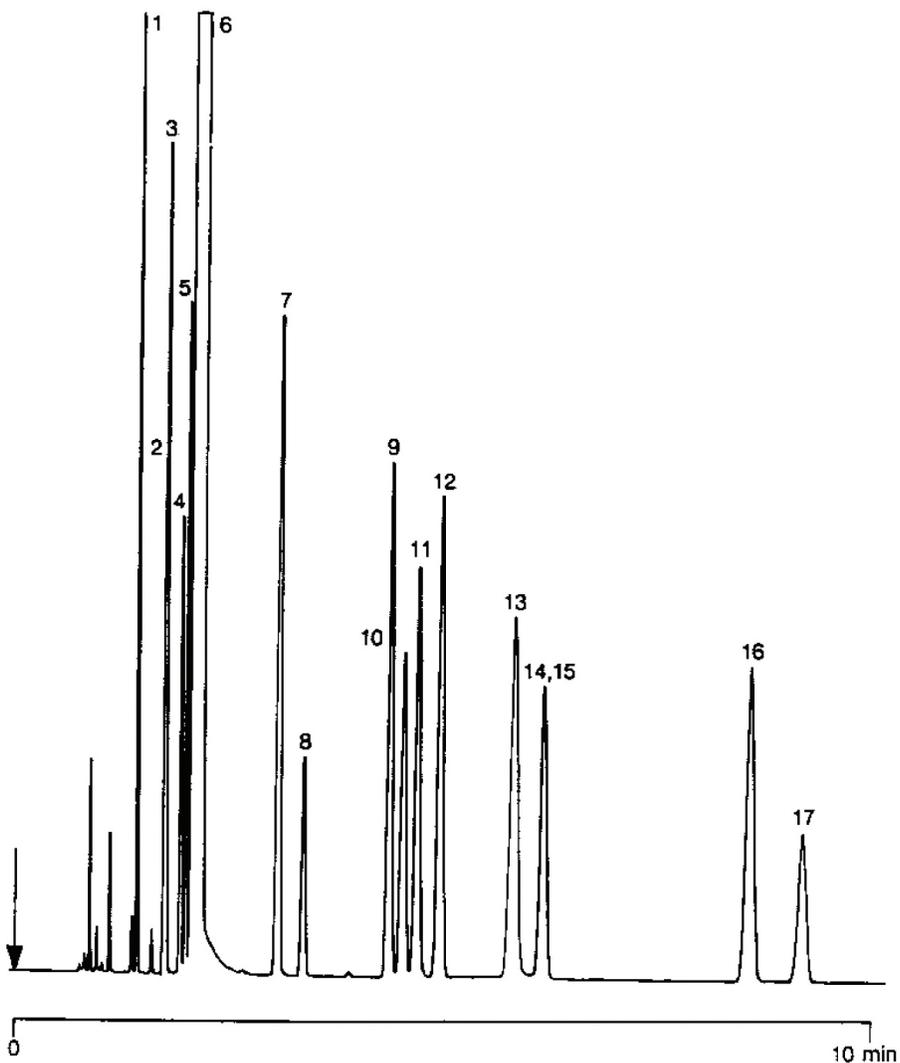
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Wax for volatile amines and diamines,
0.32 mm x 25 m fused silica WCOT CP-Wax
(df = 1.2 μ m) (Part no. CP7422)
Temperature : 120 °C
Carrier Gas : N₂, 40 kPa (0.4 bar, 5.7 psi)
Injector : Split, 120 mL/min
Detector : FID, 16 x 10⁻¹² Afs
Sample Size : 0.05 μ L
Concentration range : 0.01%
Solvent sample : CH₂CL₂

Peak identification

1. methyl acetate
2. methanol
3. ethyl acetate
4. 2-propanol
5. ethanol
6. dichloromethane
7. methyl iso-butylketone
8. 1-propanol
9. butyl acetate
10. isobutanol
11. 2-methyl-2-pentanol
12. 3-pentanol
13. iso-amyl acetate
14. 1-butanol
15. mesityloxiide
16. 3-hexanol
17. 2-methyl-1-butanol +
3-methyl-1-butanol



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

A00696



Agilent Technologies