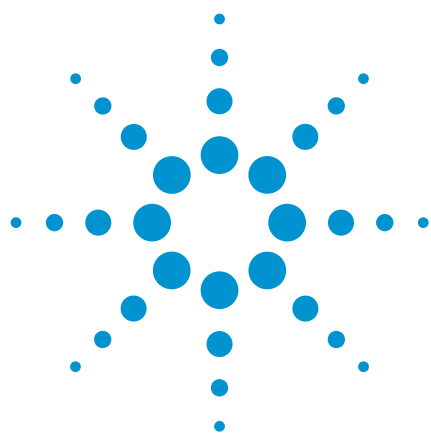




Volatile ketones and alcohols

Separation of volatile ketones and alcohols in water

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 5 CB column separates nine volatile ketones and alcohols in water in 18 minutes.



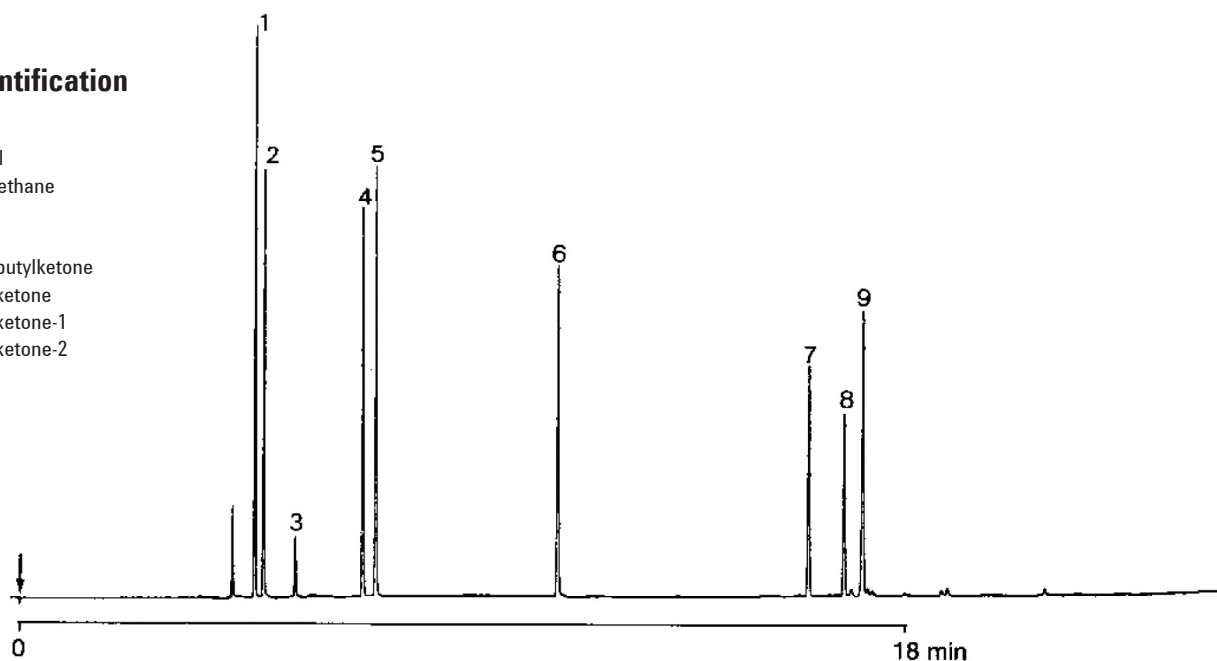
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 5 CB, 0.32 mm x 50 m fused silica
WCOT CP-Sil 5 CB (1.2 μ m) (Part no. CP7770)
Temperature : 30 °C $\rightarrow$ 250 °C, 10 °C/min
Carrier Gas : He, 100 kPa (1.0 bar, 14.5 psi), 23 cm/s
Injector : Splitter, 60 mL/min
T = 250 °C
Detector : FID, 32×10^{-12} Afs
T = 275 °C
Sample Size : 0.1 μ L
Concentration Range : 100 ppm/component
Solvent Sample : water

Peak identification

1. acetone
2. 2-propanol
3. dichloromethane
4. butanone
5. 2-butanol
6. methylisobutylketone
7. ethylamylketone
8. diisobutylketone-1
9. diisobutylketone-2



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

A00120



Agilent Technologies