



Solvents

Analysis of solvents in butanone

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Wax 57 CB column separates ten alcohol, ketone, alkane and ether solvents in butanone in 30 minutes.



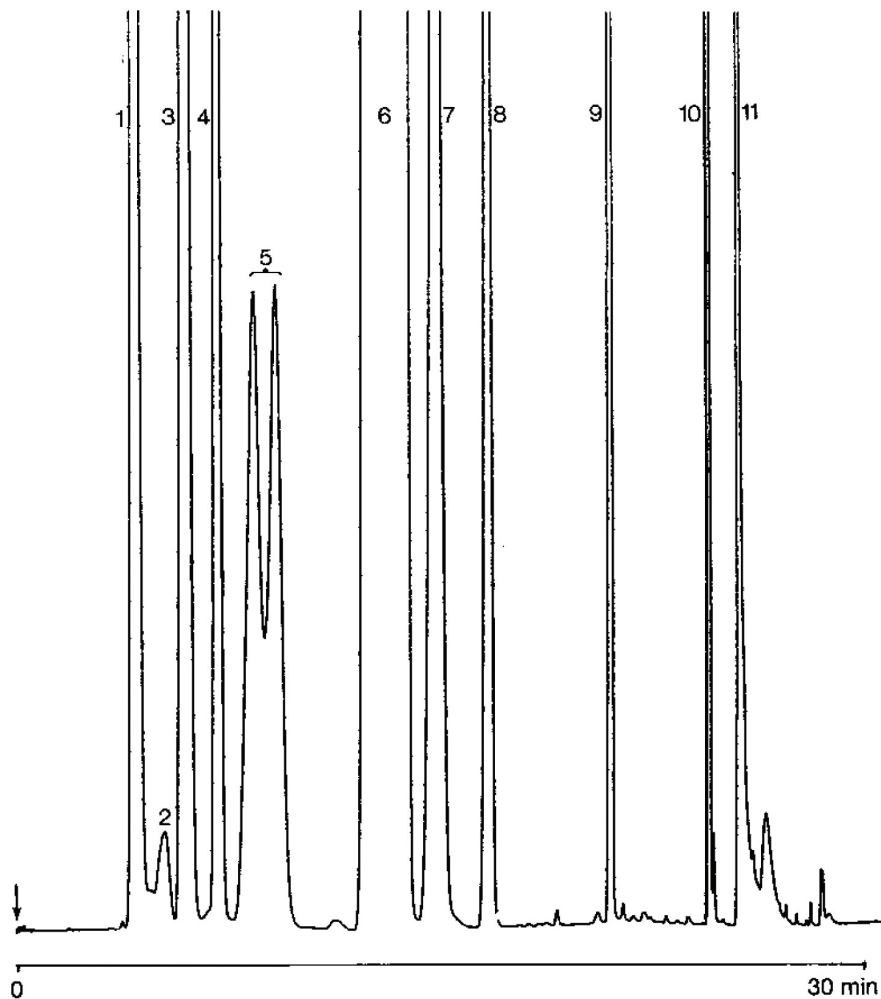
Agilent Technologies

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 : 35 °C (15 min) $\rightarrow$ 175 °C, 10 °C/min
Carrier Gas : H₂, 55 kPa (0.55 bar, 8 psi), 30 cm/s
Injector : Splitter, 40 mL/min
T = 250 °C
Detector : FID, 10 x 10⁻¹² Afs
T = 250 °C
Sample Size : 1 μ L
Concentration Range : appr. 0.5 %
Solvent Sample : butanone

Peak identification

1. cyclohexane
2. C₈-hydrocarbons
3. n-octane
4. acetone
5. sec-butylether
6. butanone
7. tert. butanol
8. 2-propanol
9. 2-butanol
10. 1-butanol
11. 5-methyl-3-heptanone



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

A00167



Agilent Technologies