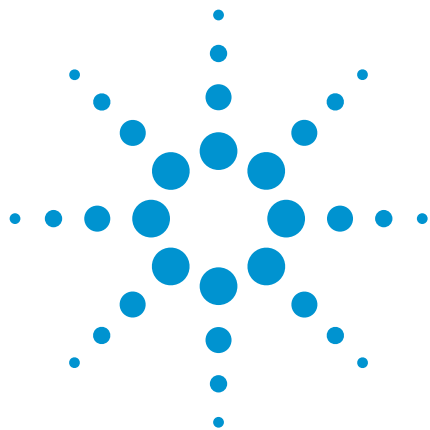




Solvents

Analysis of impurities in acetone

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Wax 52 CB column separates eight impurities in acetone in ten minutes.



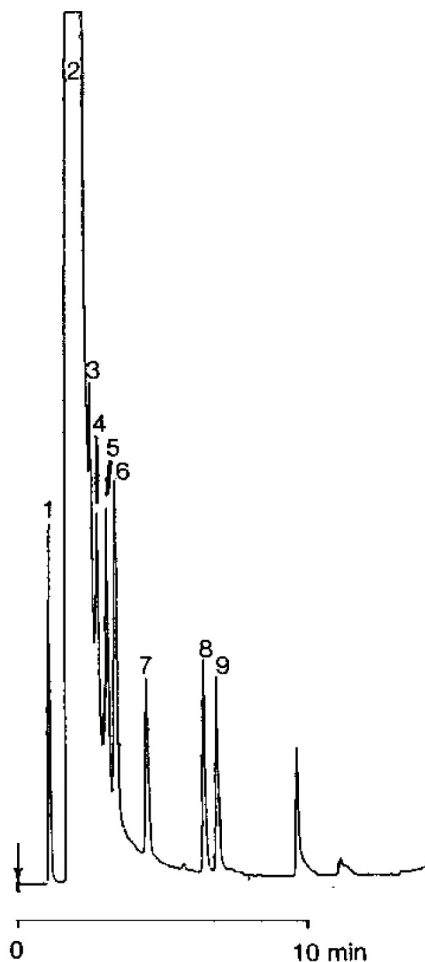
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 52 CB, 0.53 mm x 25 m fused silica
WCOT CP-Sil 52 CB (2.0 μ m) (Part no. CP7658)
Temperature : 35 °C (2 min) $\rightarrow$ 250 °C, 10 °C/min
Carrier Gas : N₂, 47 cm/s (10 mL/min)
Injector : Direct
T = 250 °C
Detector : FID, 100 x 10⁻¹² Afs
T = 275 °C
Sample Size : 0.2 μ L
Solvent Sample : acetone

Peak identification

1. diisopropylether + acetaldehyde
2. acetone
3. methanol
4. 2-butanone
5. ethanol + 2-propanol
6. benzene
7. methylisobutylketone (4-methyl-2-pentanone)
8. mesityloxide
9. methylisobutylcarbinol (4-methyl-2-pentanol)



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

A00170



Agilent Technologies