



Acetic acid and derivatives

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 8 CB column separates five compounds in a sample of acetic acid in ten minutes.



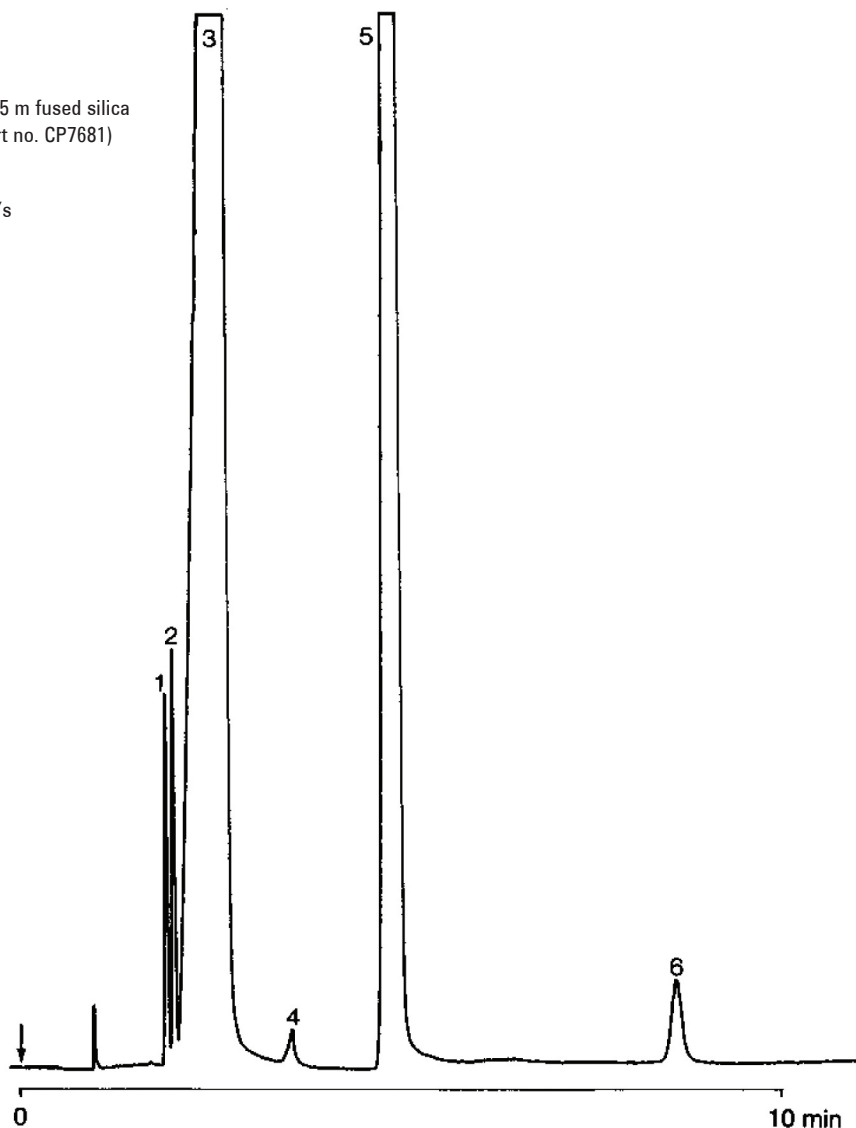
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 8 CB, 0.32 mm x 25 m fused silica
WCOT CP-Sil 8 CB (5.0 μ m) (Part no. CP7681)
Temperature : 95 °C
Carrier Gas : H₂, 40 kPa (0.4 bar, 6 psi) 29 cm/s
Injector : Splitter, 30 mL/min
T = 175 °C
Detector : micro TCD, 16 x 10⁻¹² Afs
T = 180 °C, T filament = 260 °C
Sample Size : 0.12 μ L
Concentration Range : 1% or 48% v/v

Peak identification

- | | | |
|----|------------------|---------|
| 1. | methylacetate | 1% v/v |
| 2. | methyl iodide | 1% v/v |
| 3. | acetic acid | 48% v/v |
| 4. | propionic acid | 1% v/v |
| 5. | acetic anhydride | 48% v/v |
| 6. | unknown | |



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

A00196



Agilent Technologies