



Fatty acids and esters, C₂

Determination of impurities in acetic acid

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Wax 52 CB column separates four C₂ impurities in acetic acid in ten minutes.



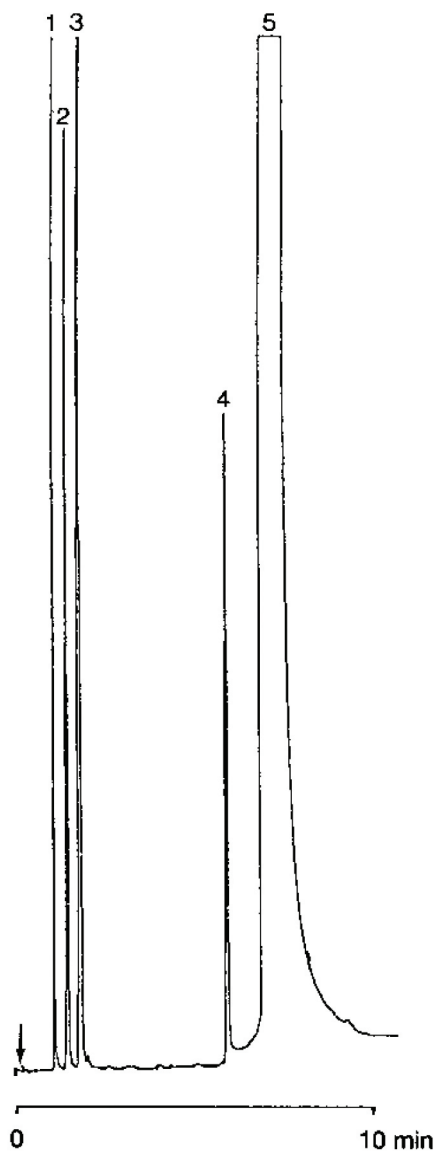
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Wax 52 CB, 0.53 mm x 25 m fused silica
WCOT CP-Wax 52 CB (2.0 μ m) (Part no. CP7658)
Temperature : 50 °C \rightarrow 200 °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 : acetic acid

Peak identification

1. acetaldehyde
2. ethylformate
3. ethylacetate
4. acetic anhydride
5. acetic acid



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

A00184



Agilent Technologies