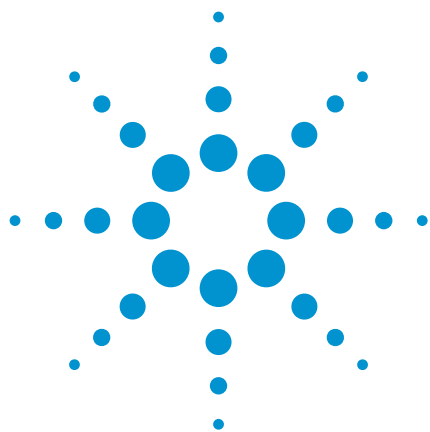




Acetates, $C_1 - C_8$

Separation of acetate esters

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Wax 52 CB column separates 13 C_1 to C_8 acetate esters in 12 minutes.



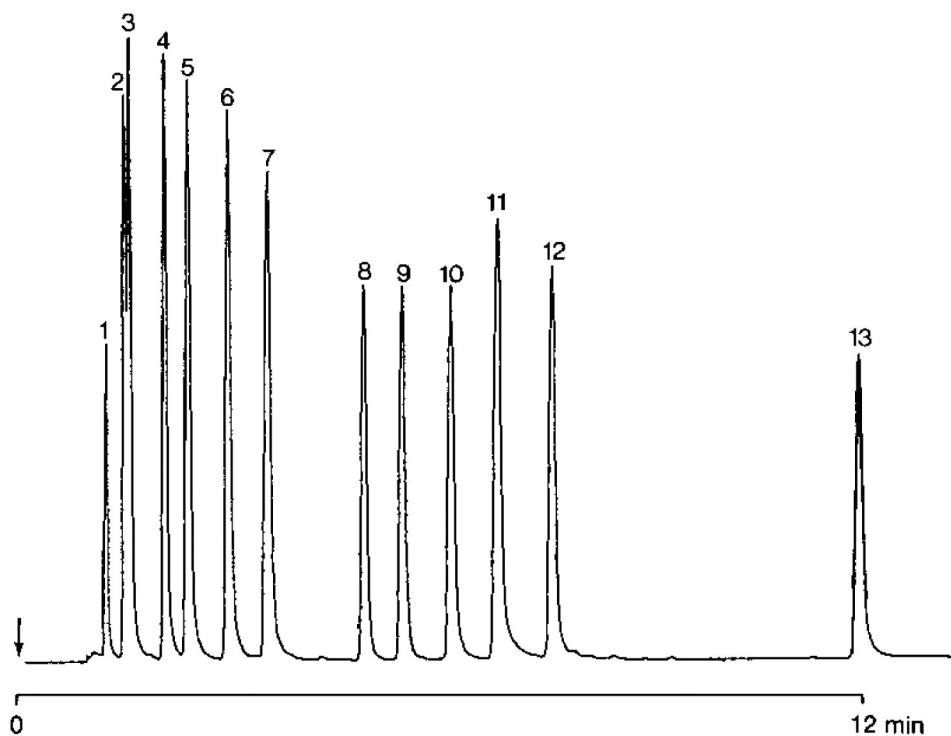
Agilent Technologies

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 : water

Peak identification

1. methyl acetate
2. ethyl acetate
3. isopropyl acetate
4. propyl acetate
5. isobutyl acetate
6. butyl acetate
7. amyl acetate
8. methylglycol acetate
9. ethylglycol acetate
10. methoxybutyl acetate
11. 2-ethylhexyl acetate
12. butylglycol acetate
13. butyldiglycol acetate



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

A00171



Agilent Technologies