

## **Solvents**

# Separation of solvent acetate esters

## Application Note

Materials Testing & Research

### **Authors**

Agilent Technologies, Inc.

### **Introduction**

Gas chromatography with an Agilent CP-Select 624 CB column separates 18 solvent acetate esters in 20 minutes



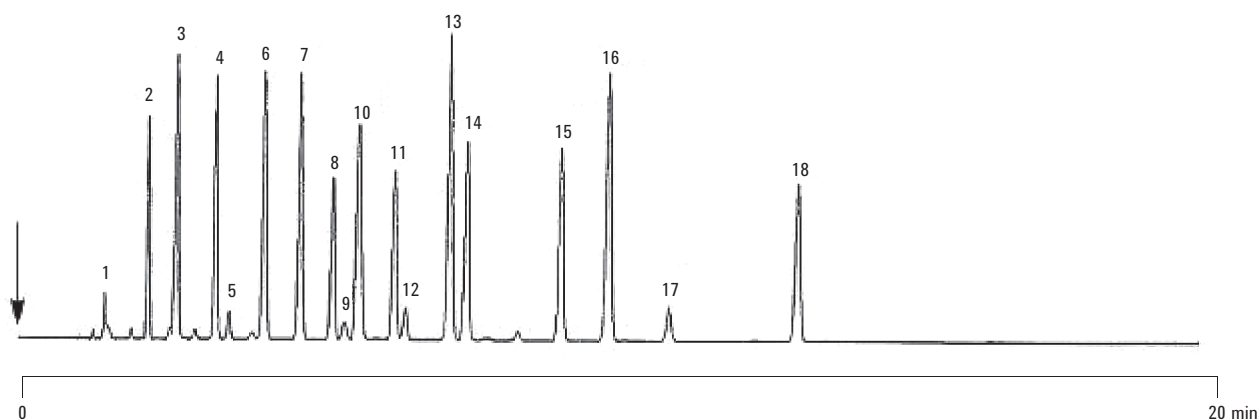
**Agilent Technologies**

## Conditions

Technique : GC-wide-bore  
Column : Agilent CP-Select 624 CB fused silica WCOT  
30 m x 0.53 mm, fused silica WCOT (df = 3.0  $\mu$ m)  
(Part no. CP7416)  
Temperature : 50 °C  $\rightarrow$  200 °C. 10 °C/min  
Carrier Gas : N<sub>2</sub>, 10 mL/min  
Injector : Direct,  
T = 250 °C  
Detector : FID  
T = 250 °C  
Sample Size : 0.02  $\mu$ L  
Solvent Sample : solvents mixture

## Peak identification

1. Methyl acetate
2. Ethyl acetate
3. Isopropyl acetate
4. Propyl acetate
5. Impurity
6. Isobutyl acetate
7. Butyl acetate
8. Methyl glycol acetate
9. Impurity
10. Amyl acetate
11. 2-Ethoxyethyl acetate (ethyl glycol acetate)
12. Impurity
13. 2-Ethylbutyl acetate
14. Methoxybutyl acetate
15. Butyl glycol acetate
16. 2-Ethylhexyl acetate
17. Impurity
18. Butyl diglycol acetate



[www.agilent.com/chem](http://www.agilent.com/chem)

This information is subject to change without notice.

© Agilent Technologies, Inc. 2011, 2017

Printed in the USA

July 18, 2017

First published prior to 11 May, 2010

A01412



**Agilent Technologies**