



# Solvents

## Application Note

Materials Testing & Research

### Authors

Agilent Technologies, Inc.

### Introduction

Gas chromatography with an Agilent CP-Select 624 CB column separates 25 constituents in a toluene sample in 17 minutes.



**Agilent Technologies**

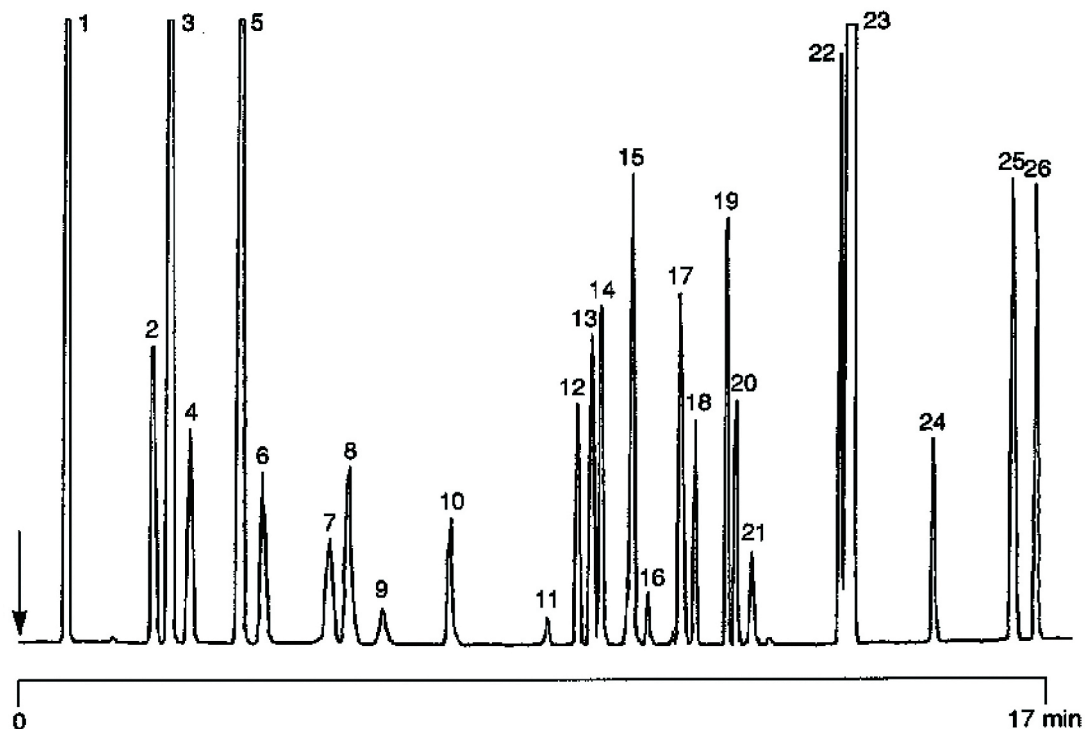
## Conditions

Technique : GC-capillary  
Column : Agilent CP-Select 624 CB 0.32 mm x 30m CB fused silica WCOT CP-Select 624 CB (df = 1.8 µm)  
(Part no. CP7414)  
Temperature : 40 °C (7 min) → 245 °C , 30 °C /min  
Carrier Gas : He, 50 kPa (0.5 bar, 7.2 psi)  
Injector : Splitter, 50 mL/min  
T = 250 °C  
Detector : FID  
T = 250 °C  
Sample Size : 0.1 µL  
Concentration Range : 1 %, except methanol, ethanol and acetone 5%  
Solvent Sample : toluene

Courtesy : Mr. B.M.M. van Genugten Diosynth B.V.,  
The Netherlands

## Peak identification

1. methanol
2. pentane
3. ethanol
4. diethyl ether
5. acetone
6. 2-propanol
7. dichloromethane
8. 2-methyl-2-propanol
9. C<sub>6</sub>-isomer
10. C<sub>6</sub>-isomer
11. C<sub>6</sub>-isomer
12. ethyl acetate
13. 2-butanol
14. tetrahydrofuran
15. cyclohexane
16. C<sub>7</sub>-isomer
17. triethylamine
18. C<sub>7</sub>-isomer
19. 1-butanol
20. trichloroethylene
21. C<sub>7</sub>-isomer
22. pyridine
23. toluene
24. N,N-dimethylformamide
25. cyclohexanol
26. cyclohexanone



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

A01281



**Agilent Technologies**