



## Solvents

# Separation of solvents on a wide bore capillary column

## Application Note

Materials Testing & Research

### Authors

Agilent Technologies, Inc.

### Introduction

Gas chromatography with an Agilent CP-Sil 5 CB column separates 33 solvents and other organics in 13 minutes.



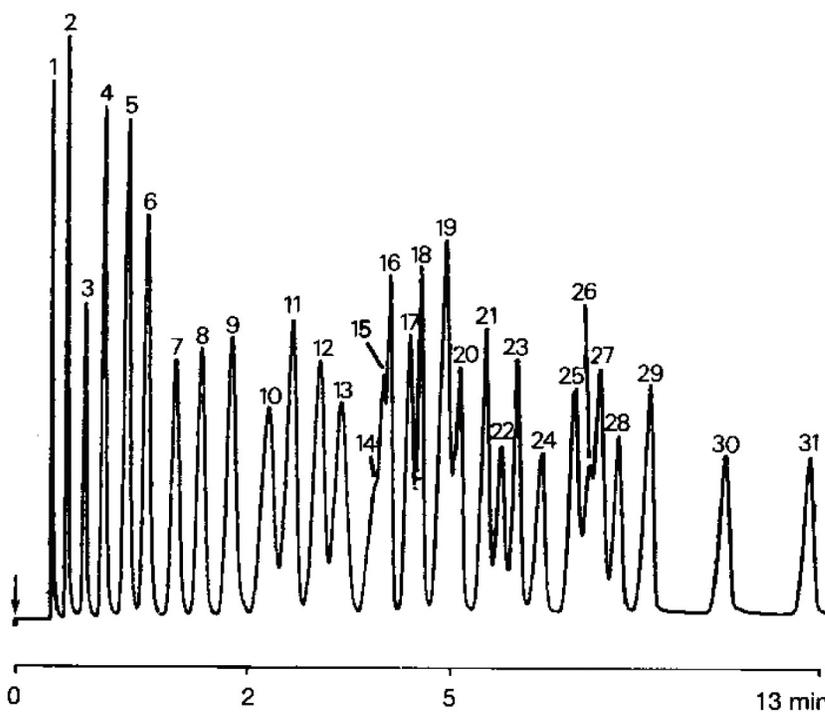
**Agilent Technologies**

## Conditions

Technique : GC-capillary  
Column : Agilent CP-Sil 5 CB, 0.53 mm x 10 m fused silica  
WCOT CP-Sil 5 CB (5.0 µm) (Part no. CP7645)  
Temperature : 50 °C → 250 °C, 5 °C/min  
Carrier Gas : N<sub>2</sub>, 10 kPa (0.1 bar, 1.5 psi), 50 cm/s  
Injector : Direct  
T = 250 °C  
Detector : FID, 100 x 10<sup>12</sup> Afs  
T = 275 °C  
Sample Size : 0.2 µL  
Concentration Range : 3 %/component

## Peak identification

1. methanol
2. acetone
3. nitromethane
4. butanone
5. isobutanol
6. methylisopropylketone
7. isobutylformate
8. propylacetate
9. methylisobutylketone
10. methylisobutylcarbinol
11. cyclopentanone
12. mesityloxiide
13. diisobutylether
14. methoxybutanol
15. 2-methyl-1-pentanol
16. chlorobenzene
17. ethylbenzene
18. p-xylene
19. o-xylene + pentoxone\* + dibutylether
20. 1,2,3-trichloropropane
21. cumene
22. ethylamylketone
23. m-chlorotoluene
24. 1-heptanol
25. tert. butylbenzene
26. 1,4-dichlorobenzene
27. sec. butylbenzene
28. 1,2-dichlorobenzene
29. acetophenone
30. 3,5,5-trimethyl-2-cyclohexen-1-one (isophorone)
31. 1,2,4-trichlorobenzene



[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

A00128



**Agilent Technologies**