



Solvents

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Wax 52 CB column separates 35 solvents in 12 minutes.



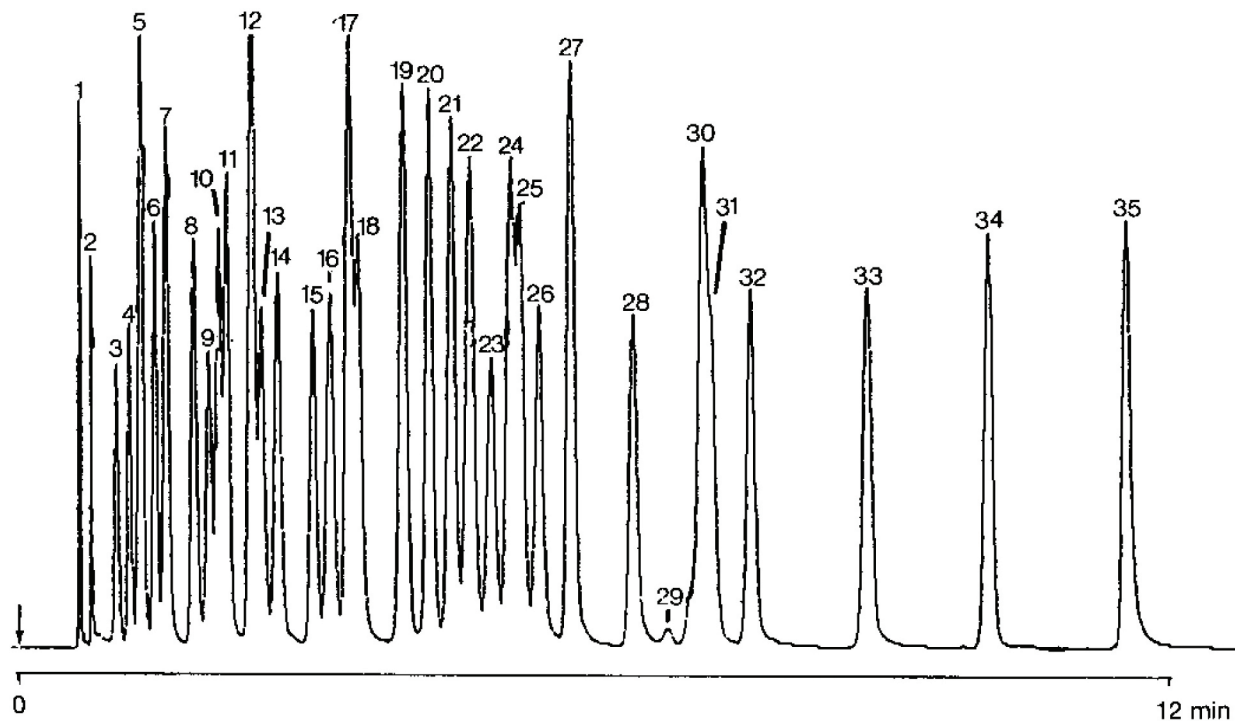
Agilent Technologies

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
Concentration Range : equal conc.

Peak identification

- | | |
|-----------------------------|--|
| 1. diethylether | 20. chlorobenzene |
| 2. 1,1-dichloroethene | 21. 4-ethyltoluene |
| 3. methylacetate | 22. tert. butylbenzene |
| 4. trans-1,2-dichloroethene | 23. pentoxone
(= 4-methoxy-4-methylpentanone-2) |
| 5. tert. butanol | 24. cumene (isopropylbenzene) |
| 6. ethanol | 25. cyclohexanone |
| 7. methylisopropylketone | 26. 2-heptanol |
| 8. propylacetate | 27. α -methylstyrene |
| 9. trichloroethene | 28. cyclohexanol |
| 10. 2-butanol | 29. 1,3 dichlorobenzene |
| 11. 1-propanol | 30. 1,2,3,5-tetramethylbenzene |
| 12. toluene | 31. butylglycolacetate |
| 13. 1,2 dichloroethane | 32. 1,2-dichlorobenzene |
| 14. butylacetate | 33. isophorone
(= 3,5,5-trimethyl-2-cyclohexen-1-one) |
| 15. amylacetate | 34. α,α -dichlorotoluene |
| 16. 2-nitropropane | 35. benzylalcohol |
| 17. p-xylene | |
| 18. 4-methyl-2-pentanol | |
| 19. o-xylene | |



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

A00185



Agilent Technologies