



Solvents

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography using an Agilent CP-Sil 5 CB column separates 19 solvents in 30 minutes.



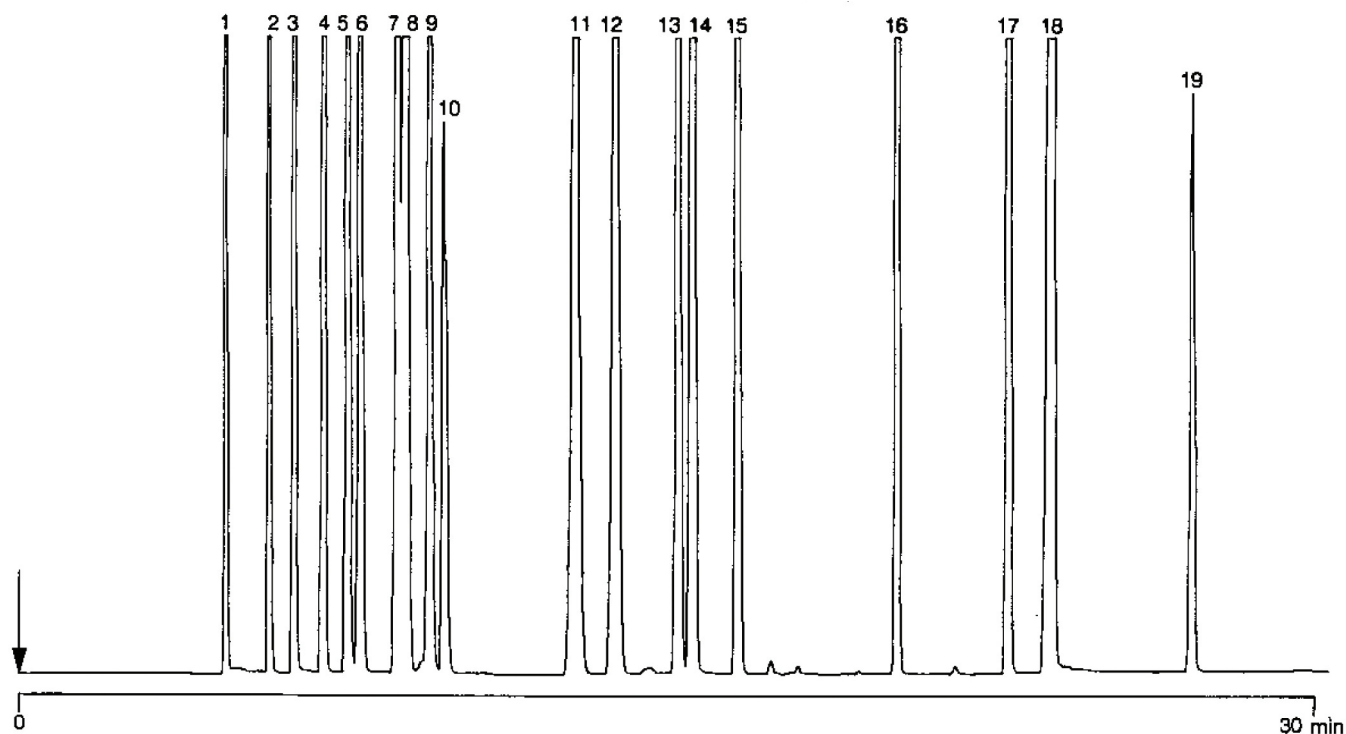
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 5 CB, 0.32 mm x 50 m fused silica
WCOT CP-Sil 5 CB (df = 5.0 µm) (Part no. CP7690)
Temperature : 70 °C (10 min) → 200 °C, 20 °C/min; 200 °C (15 min)
Carrier Gas : He, 80 kPa (0.8 bar, 11.4 psi)
Injector : Splitter, 60 mL/min
T = 250 °C
Detector : FID
T = 250 °C
Sample Size : 1.0 µL
Courtesy : P. de Haan, Analytical Research, Duphar B.V.,
Weesp, The Netherlands

Peak identification

1. methanol
2. methylformate
3. ethanol
4. acetonitrile
5. acetone
6. isopropanol
7. diethylether
8. n-pentane
9. methylacetate
10. dichloromethane
11. methyl-tert-butylether (MTBE)
12. methyl ethylketone (MEK)
13. n-hexane
14. ethyl acetate
15. tetrahydrofuran (THF)
16. n-heptane
17. toluene
18. dimethylsulfoxide (DMSO)
19. 1,4-dichlorobutane



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

A00718



Agilent Technologies