



Solvents

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT Q-HT column separates 18 solvents in 50 minutes.



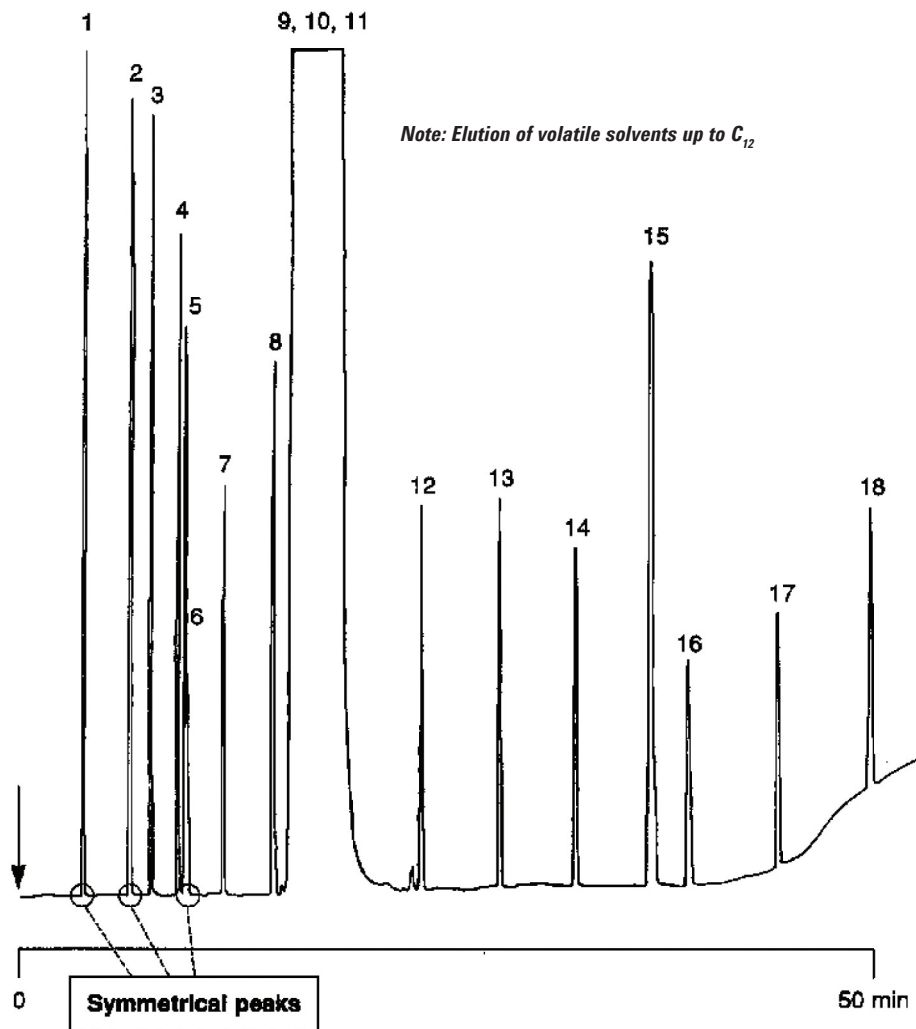
Agilent Technologies

Conditions

Technique : GC-wide-bore
Column : Agilent PoraPLOT Q-HT, 0.53 mm x 25 m fused silica
PLOT PoraPLOT (df = 20 µm) (Part no. CP7559)
Temperature : 100 °C → 250 °C, 5 °C/min
250 °C (10 min) → 290 °C, 5 °C/min
Carrier Gas : H₂, 45 kPa (0.45 bar, 6.4 psi)
Injector : Split, T = 270 °C
Detector : FID, T = 290 °C
Concentration Range : 0.1%
Solvent Sample : cyclohexane

Peak identification

1. methanol
2. ethanol
3. acetonitrile
4. acetone
5. isopropanol
6. dichloromethane
7. pentane
8. ethyl acetate
9. hexane
10. benzene
11. cyclohexane
12. toluene
13. ethylbenzene
14. propylbenzene
15. decane
16. butylbenzene
17. undecane
18. dodecane



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

A01256



Agilent Technologies