



Solvents

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT Q column separates 11 solvents in four minutes.



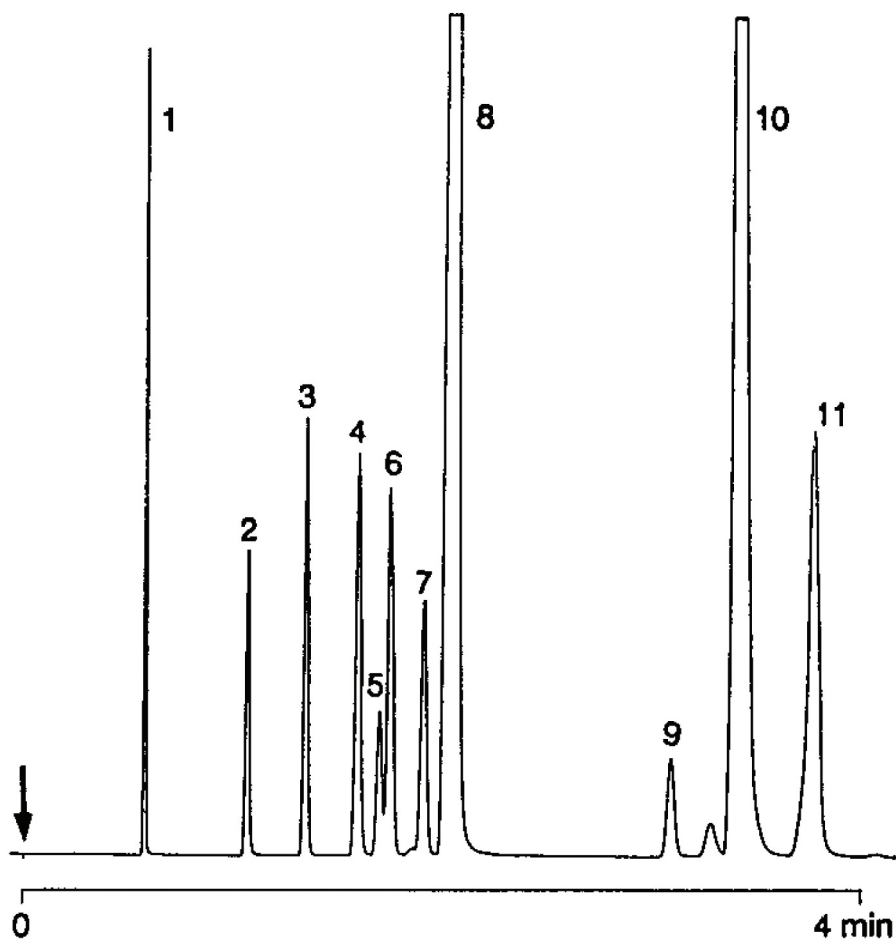
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent PoraPLOT Q, 0.53 mm x 25 m fused silica
PoraPLOT Q (df = 20 μ m) (Part no. CP7554)
Temperature : 100 °C \rightarrow 200 °C, 10 °C/min
Carrier Gas : H₂, 40 kPa (0.4 bar, 6psi); 15 mL/min
Injector : Direct
Detector : FID, 128 x 10⁻¹² Afs
Sample Size : 10 μ L headspace
Concentration range : 1 - 3%

Peak identification

1. methanol
2. ethanol
3. acetonitrile
4. acetone
5. isopropanol
6. dichloromethane
7. methyl acetate
8. n-pentane
9. ethyl acetate
10. n-hexane
11. benzene



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

A00438



Agilent Technologies