



Solvents

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT Q column separates eight solvents in 25 minutes.



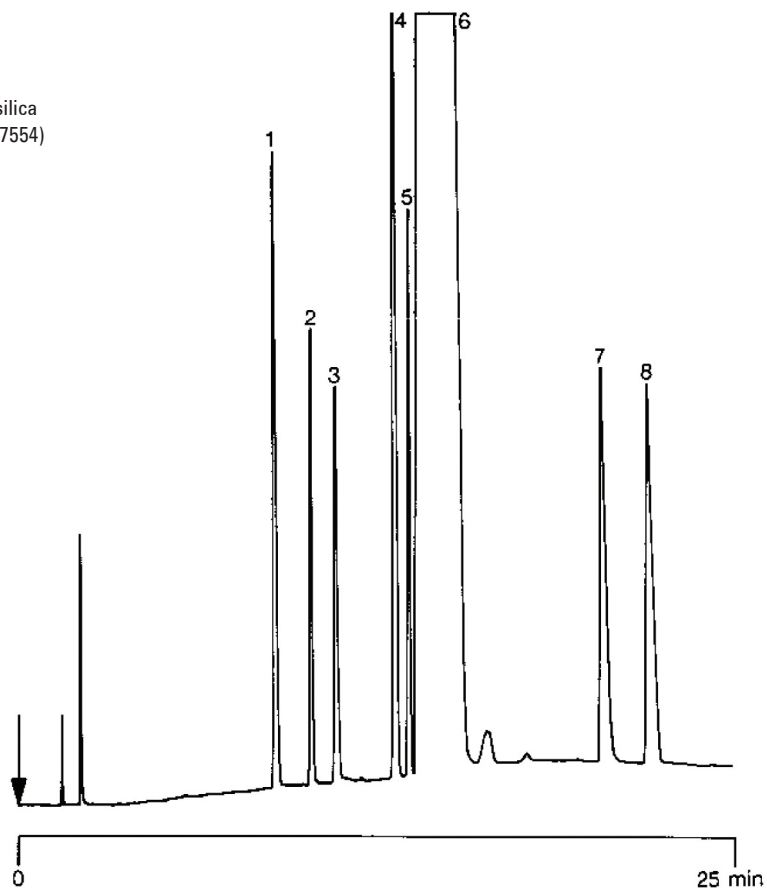
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent PoraPLOT Q, 0.53 mm x 25 m fused silica
PLOT PoraPLOT Q (df = 20 μ m) (Part no. CP7554)
Temperature : 100 °C (3 min) \rightarrow 200 °C, 7 °C/min
Carrier Gas : He, 50 kPa (0.5 bar, 7 psi), 4.5 mL/min
Injector : Splitter, 1:5
T = 200 °C
Detector : μ -TCD, x4
T = 200 °C, filaments T = 280 °C
Sample Size : 1 μ L

Peak identification

1. acetone
2. n-pentane
3. methyl acetate
4. MEK (methyl ethylketone)
5. ethyl acetate
6. n-hexane
7. MIBK (methyl isobutylketone)
8. MBK (methyl butylketone)



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

A00608



Agilent Technologies