



Solvents

The effect of column length

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with Agilent PoraPLOT Q columns of different dimensions analyzes ten solvents in 20 minutes to demonstrate the effect of column length on separation.



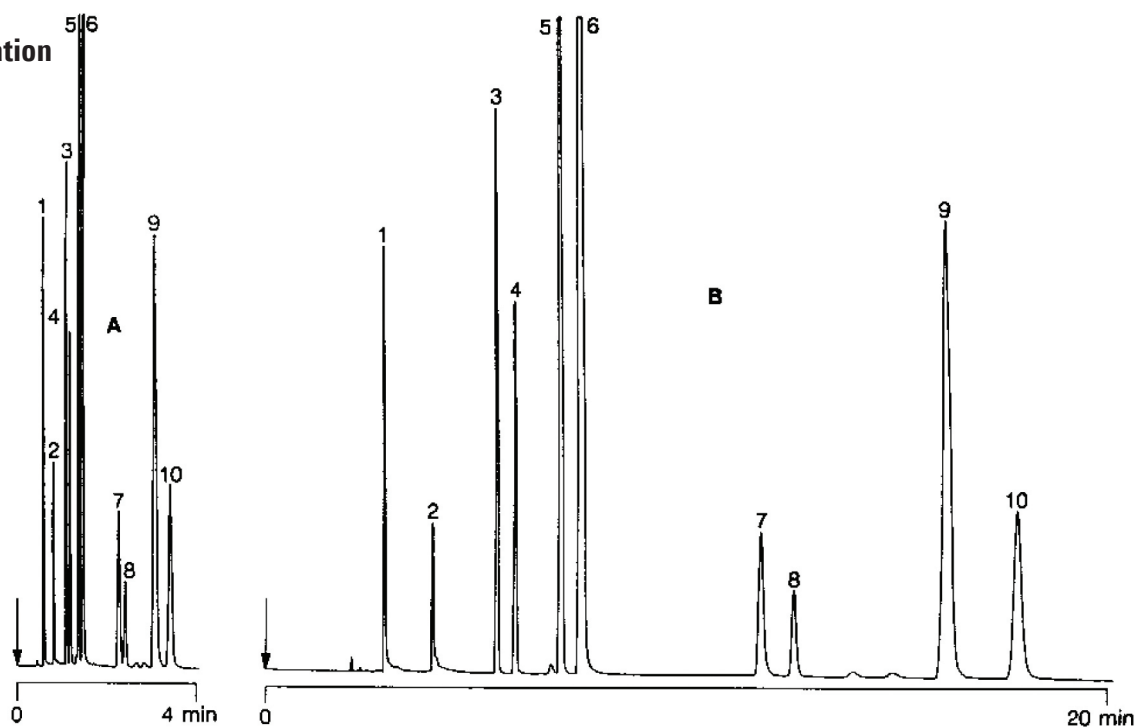
Agilent Technologies

Conditions

Technique : GC-capillary
Column : A: Agilent PoraPLOT Q, 0.53 mm x 10 m fused silica
PLOT PoraPLOT Q (df = 20 μ m) (Part no. CP7553)
B: Agilent PoraPLOT Q, 0.53 mm x 50 m fused silica
PLOT PoraPLOT Q (df = 20 μ m) (Part no. CP7555)
Temperature : 150 °C
Carrier Gas : H₂, 5 mL/min
Injector : Splitter, 100 mL/min;
T = 250 °C
Detector : FID
T = 300 °C
Sample Size : 1.5 μ L headspace
Concentration range : 1 - 10%

Peak identification

1. methanol
2. ethanol
3. acetone
4. dichloromethane
5. diethyl ether
6. pentane
7. THF
8. ethyl acetate
9. hexane
10. 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

A00667



Agilent Technologies