



# Solvents

## Application Note

Materials Testing & Research

### Authors

Agilent Technologies, Inc.

### Introduction

Gas chromatography using an Agilent PoraPLOT U column separates six solvents in a test mixture in 12 minutes.



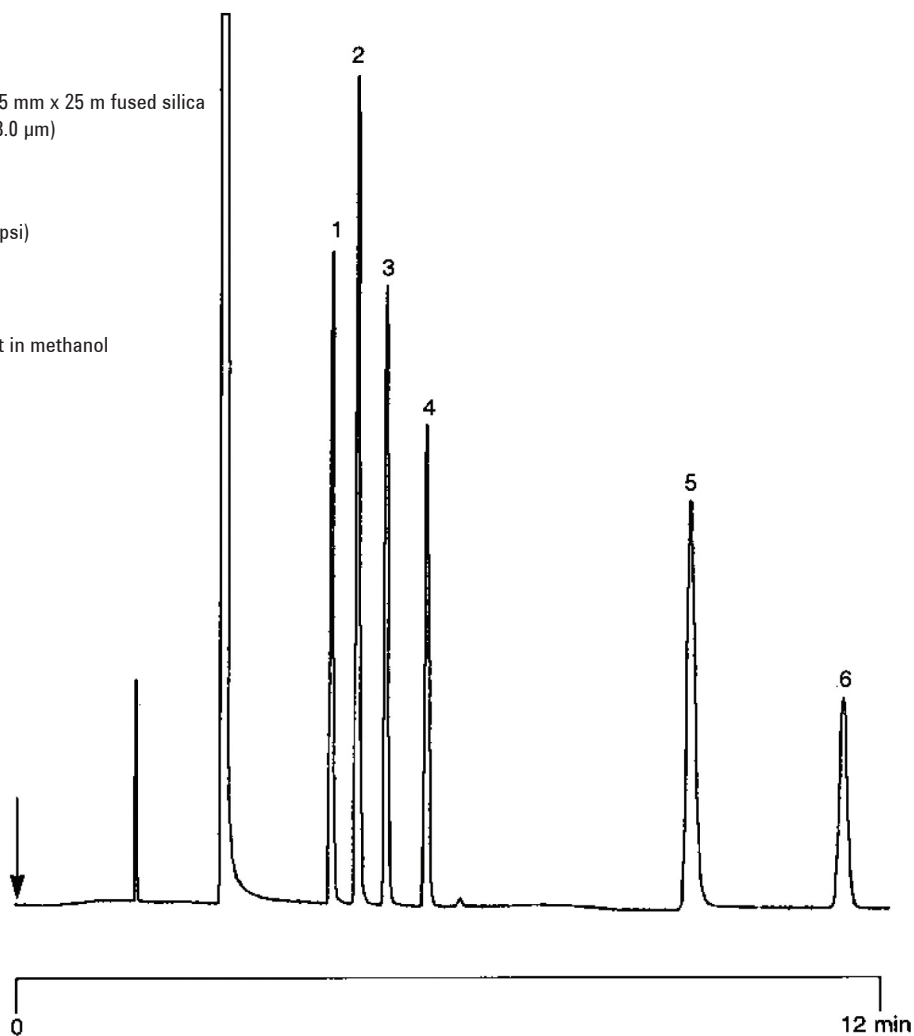
**Agilent Technologies**

## Conditions

Technique : GC-capillary  
Column : Agilent PoraPLOT U, 0.25 mm x 25 m fused silica  
PLOT PoraPLOT U (df= 8.0 µm)  
(Part no. CP7579)  
Temperature : 150 °C  
Carrier Gas : H<sub>2</sub>, 65 kPa (0.65 bar, 9.3 psi)  
Injector : Split, 100 mL/min  
Detector : FID  
Sample : 0.1% of each component in methanol

## Peak identification

1. ethanol
2. n-pentane
3. diethyl ether
4. acetone
5. n-hexane
6. ethyl acetate



[www.agilent.com/chem](http://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

A01127



**Agilent Technologies**