



Alcohols, $C_1 - C_5$

Analysis of alcohols in water

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT Q column separates 14 C_1 to C_5 alcohols in a sample of water in seven minutes.



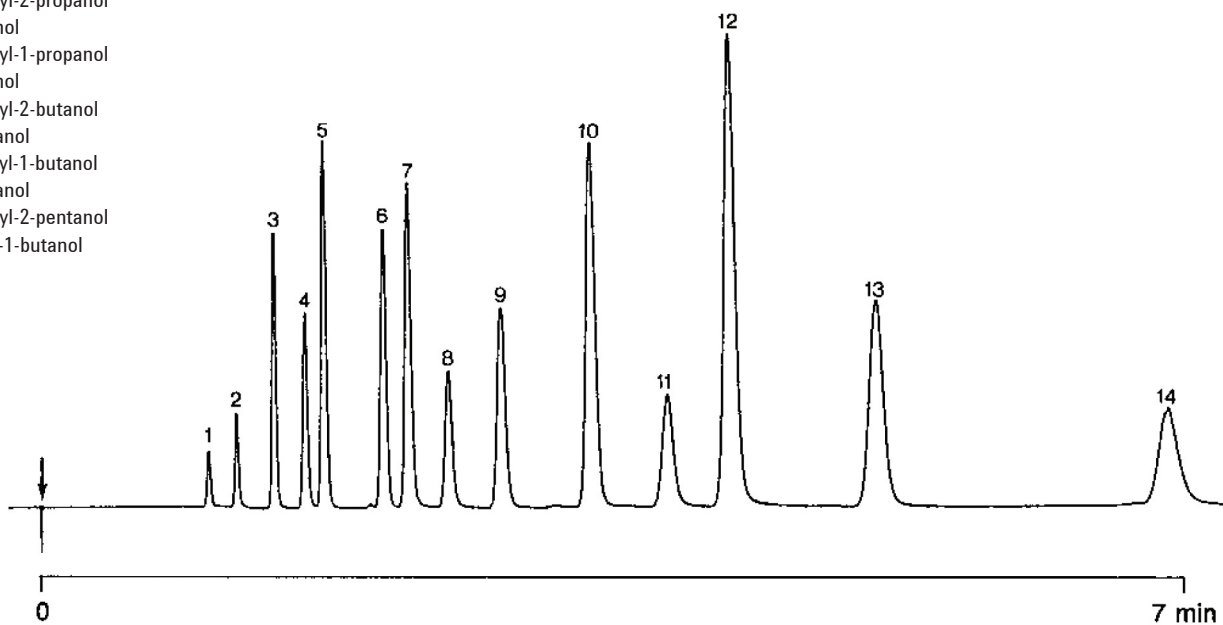
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent PoraPLOT Q, 0.32 mm x 10 m fused silica
PoraPLOT Q (10 μ m) (Part no. CP7550)
Temperature : 200 °C
Carrier Gas : H₂, 40 kPa (0.4 bar, 5.8 psi)
Injector : Splitter
Detector : FID
Concentration Range : 1%

Peak identification

1. methanol
2. ethanol
3. 2-propanol
4. 1-propanol
5. 2-methyl-2-propanol
6. 2-butanol
7. 2-methyl-1-propanol
8. 1-butanol
9. 2-methyl-2-butanol
10. 2-pentanol
11. 3-methyl-1-butanol
12. 1-pentanol
13. 4-methyl-2-pentanol
14. 2-ethyl-1-butanol



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

A00280



Agilent Technologies