



Alcohols C₁ - C₂

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT Q column separates two C₁ to C₂ alcohols in water in three minutes.



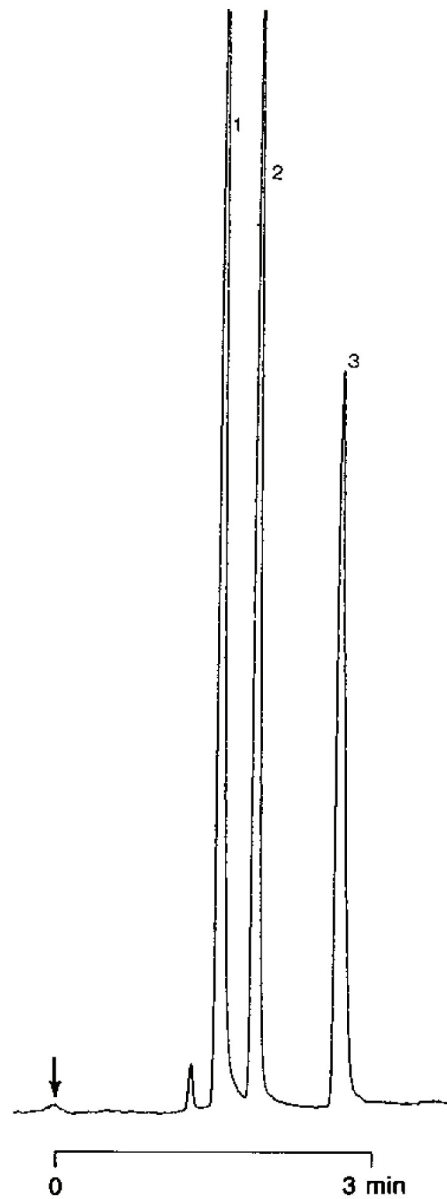
Agilent Technologies

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 : 175 $^{\circ}$ C
Carrier Gas : H₂, 50 kPa (0.5 bar, 7 psi)
Injector : Splitter
Detector : μ TCD

Peak identification

1. water
2. methanol
3. ethanol



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

A00285



Agilent Technologies