



Alcohols, C₁ – C₃

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT Q column separates four C₁ to C₃ alcohols in six minutes.



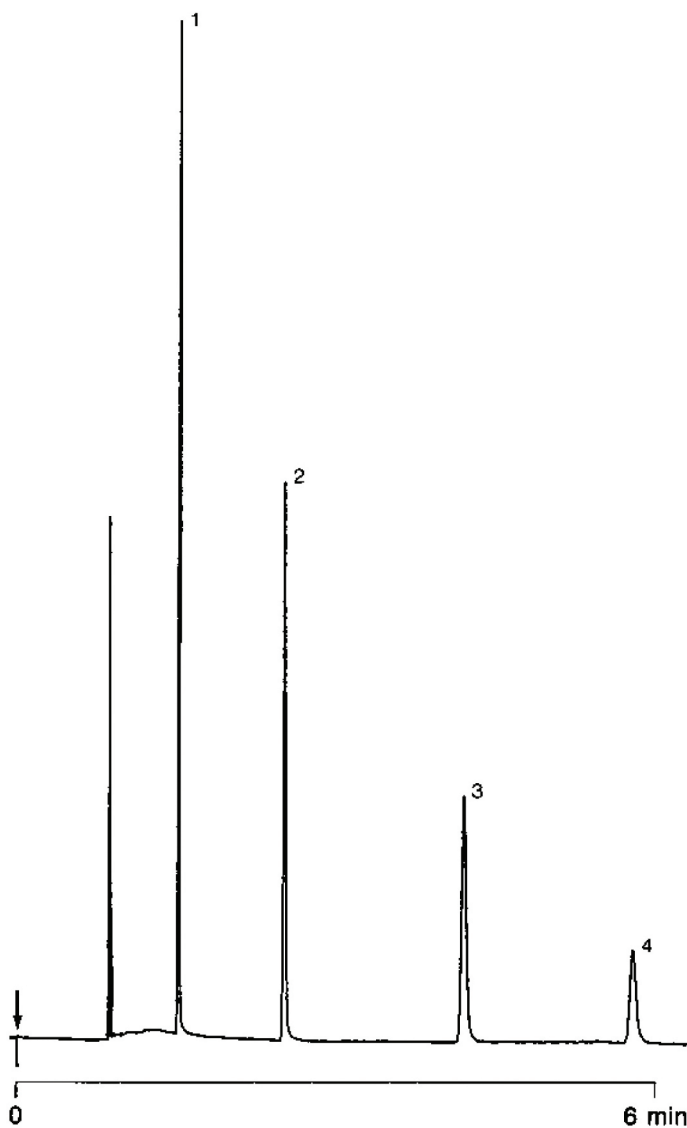
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent PoraPLOT Q, 0.32 mm x 25 m fused silica
PoraPLOT Q (10 μ m) (Part no. CP7551)
Temperature : 100 °C
Carrier Gas : H₂, 100 kPa (1.0 bar, 14.5 psi)
Injector : Splitter
Detector : FID

Peak identification

1. methanol
2. ethanol
3. isopropanol
4. propanol



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

A00289



Agilent Technologies