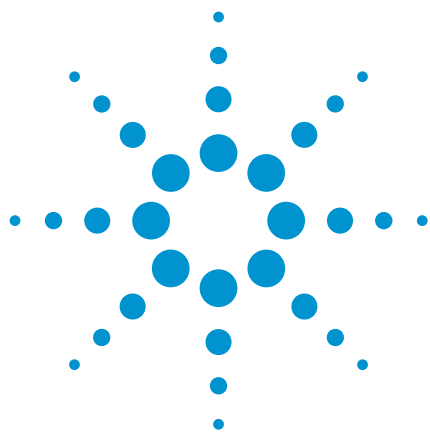




Glycols

Determination of glycerine

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT Q column separates seven glycols and glycerols in eight minutes.



Agilent Technologies

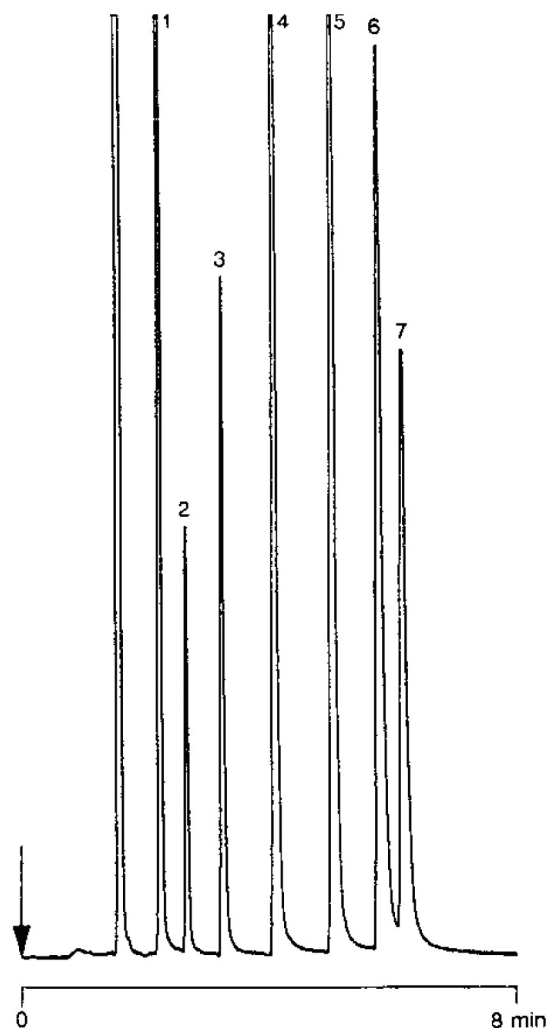
Conditions

Technique : GC-wide-bore
Column : Agilent PoraPLOT Q, 0.53 mm x 25 m fused silica
PLOT PoraPLOT Q (df = 20 µm) (Part no. CP7554)
Temperature : 240 °C
Carrier Gas : H₂, 30 kPa (0.3 bar, 4.3 psi)
Injector : Splitter
T = 220 °C
Detector : FID
T = 250 °C
Sample Size : 1 µL
Concent Range : 1%
Solvent Sample : water/methanol

Courtesy : C. Mariani, Stazione sperimentale per le industrie
degli di e dei grassi, Milano, Italy

Peak identification

1. ethylene glycol
2. 1,2-propylene glycol
3. 1,3-propylene glycol
4. 1,3-butylene glycol
5. 1,4-butylene glycol
6. glycerol (glycerine)
7. diethylene glycol



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

A00770



Agilent Technologies