



Glycols

Mono, di-,tri-and tetraethylene glycol analysis

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Wax 57 CB for Glycols and Alcohols column separates mono-, di-, tri-, and tetraethylene glycols in 15 minutes.



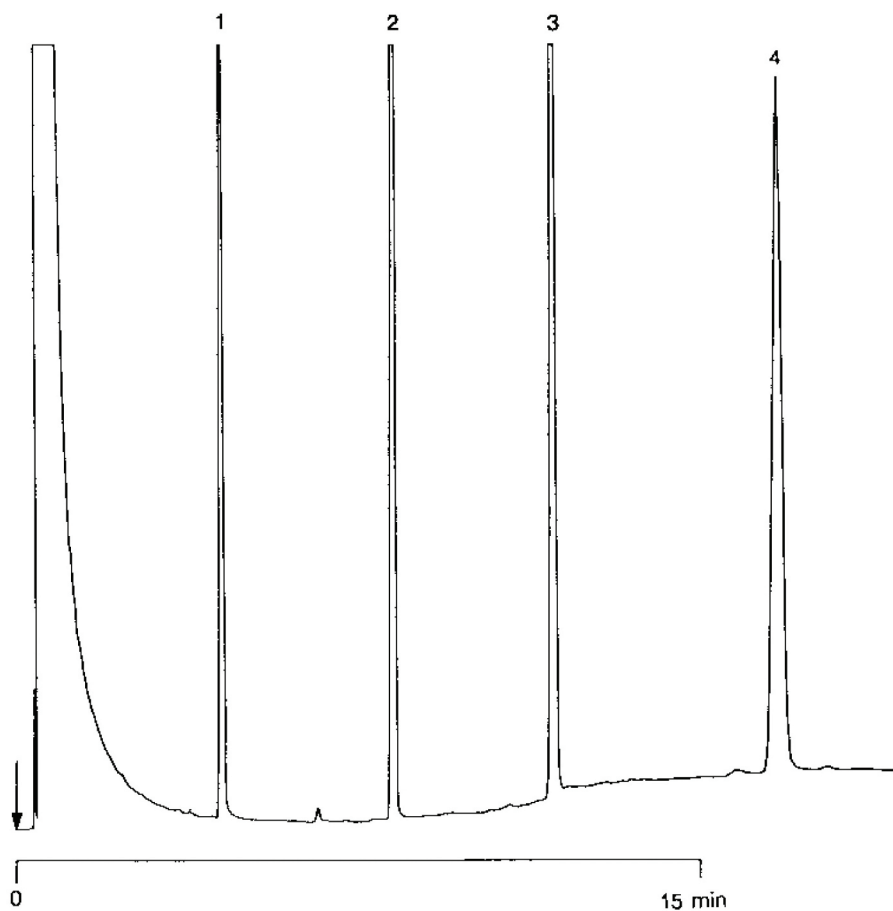
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Wax 57 CB, 0.53 mm x 25 m fused silica
WCOT CP-Wax 57 CB (df = 0.5 µm)
(Part no. CP7617)
Temperature : 100 °C → 200 °C, 10 °C/min
Carrier Gas : 50kPa (0.5 bar, 7 psi), 10 mL/min
Injector : Direct
T = 250 °C
Detector : FID, 64 x 10⁻¹² Afs
T = 250 °C
Sample Size : 0.5 µL
Concentration Range : 0.25%
Solvent Sample : ethylacetate

Peak identification

1. monoethylene glycol
2. diethylene glycol
3. triethylene glycol
4. tetraethylene 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

A00467



Agilent Technologies