

Alcohols and Glycols

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

The high polarity (CP-Index 57) of the Agilent CP-Wax 57 CB for Alcohols and Glycols polyethylene glycol phase, its inertness and resistance to water make it possible to analyze both alcohols and glycols in aqueous samples.



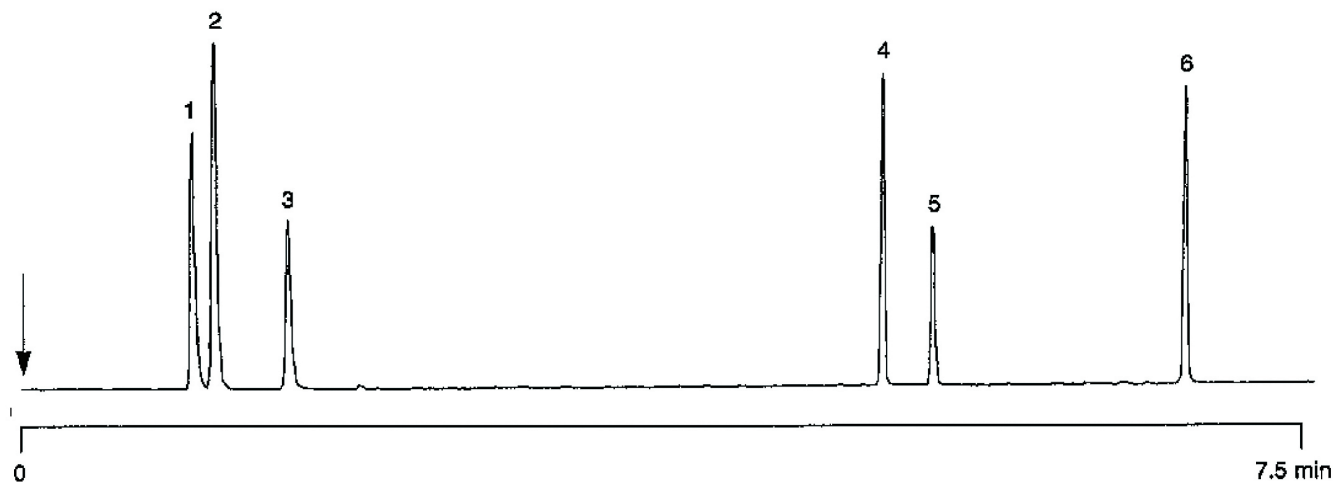
Agilent Technologies

Conditions

Technique : GC-wide-bore
Column : Agilent CP-Wax 57 CB for Alcohols and Glycols,
0.53 mm × 25 m, 1.2 μm (p/n CP7617)
Temperature : 45 °C (0.75 min) → 180 °C, 20 °C/min
Carrier Gas : He, 10 mL/min, 50 kPa (0.5 bar, 7.2 psi)
Injector : Direct,
T = 220 °C
Detector : FID
T = 230 °C
Sample Size : 0.2 μL,
fast injection, no air plug
Solvent Sample : Water

Peak identification

1. Methanol
2. Ethanol
3. 1-Propanol
4. Ethylene glycol
5. Trichloroethanol
6. Diethylene glycol



www.agilent.com/chem

This information is subject to change without notice.

© Agilent Technologies, Inc. 2015

Printed in the USA

November 30, 2015

First published prior to 11 May, 2010

A01330



Agilent Technologies