



Aromatic hydrocarbons

Analysis of low concentrations of volatile organic components in water

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Wax 57 CB column separates six volatile aromatic hydrocarbons at low concentrations in water in 13 minutes.



Agilent Technologies

Conditions

Technique : GC-capillary

Column : Agilent CP-Wax 57 CB, 0.32 mm x 50 m fused silica
WCOT CP-Wax 57 CB (0.2µm) (Part no. CP97753)

Temperature : 55 °C

Carrier Gas : He, 70 kPa (0.7 bar, 10 psi)

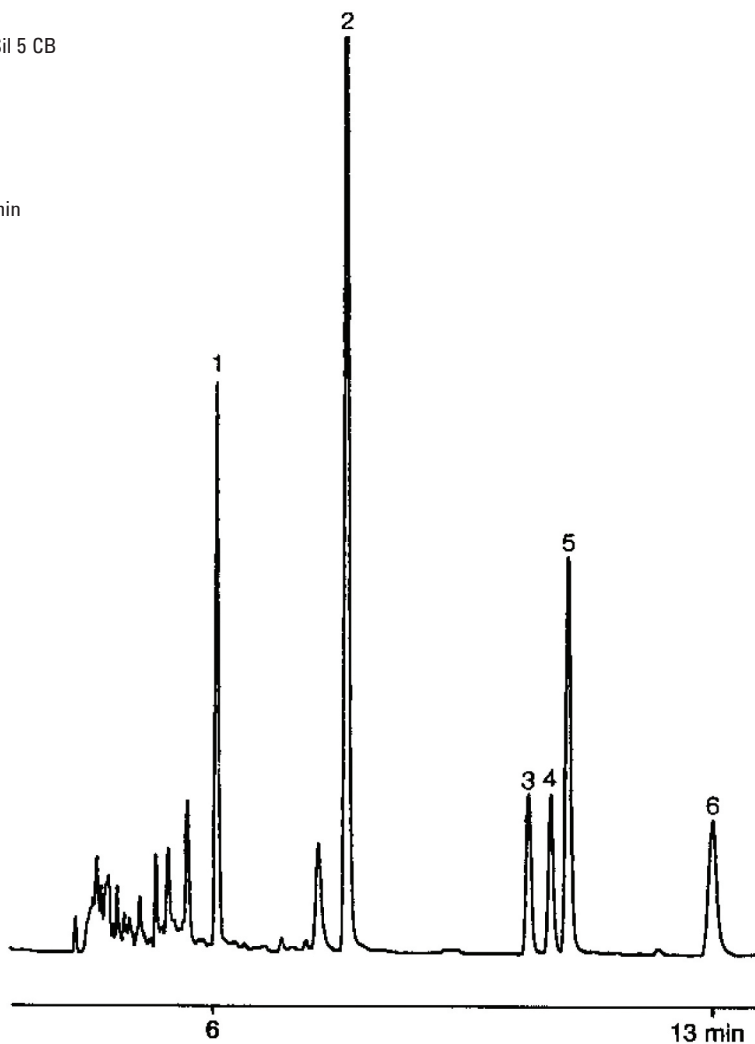
Injector : PTI
Cold Trap material 0.32 mm x 30 cm ID CP-Sil 5 CB
fused silica WCOT; (1.2 µm)
Spec.:
Cryofocussing : -120 °C
Precool time : 2 min
Purge time : 10 min
Purge flow : 10 mL/min
Desorption oven temperature : 100 °C
Condensor temperature : - 15 °C
Purge temperature : ambient
Injection temperature : 200 °C
Injection time : 5 min
GC-injection block temperature : 250 °C

Detector : FID/ECD
T= 300 °C

Sample Handling : Matrix : liquid
Sample size : 5 mL
Concentration range : 1 µg/L
Dynamic headspace : on-line
Purge (gas-stripping) : on-line

Peak identification

1. benzene
2. toluene
3. ethylbenzene
4. p-xylene
5. m-xylene
6. o-xylene



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

A00219



Agilent Technologies