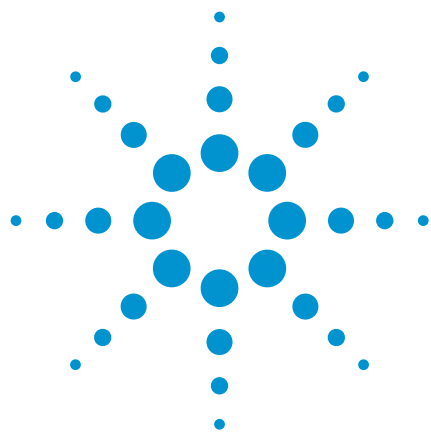




Volatile halogenated hydrocarbons

Separation of volatile halogenated hydrocarbons and other volatiles in waters

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 5 CB column separates 17 volatile halogenated hydrocarbons and other volatiles in water in 16 minutes.



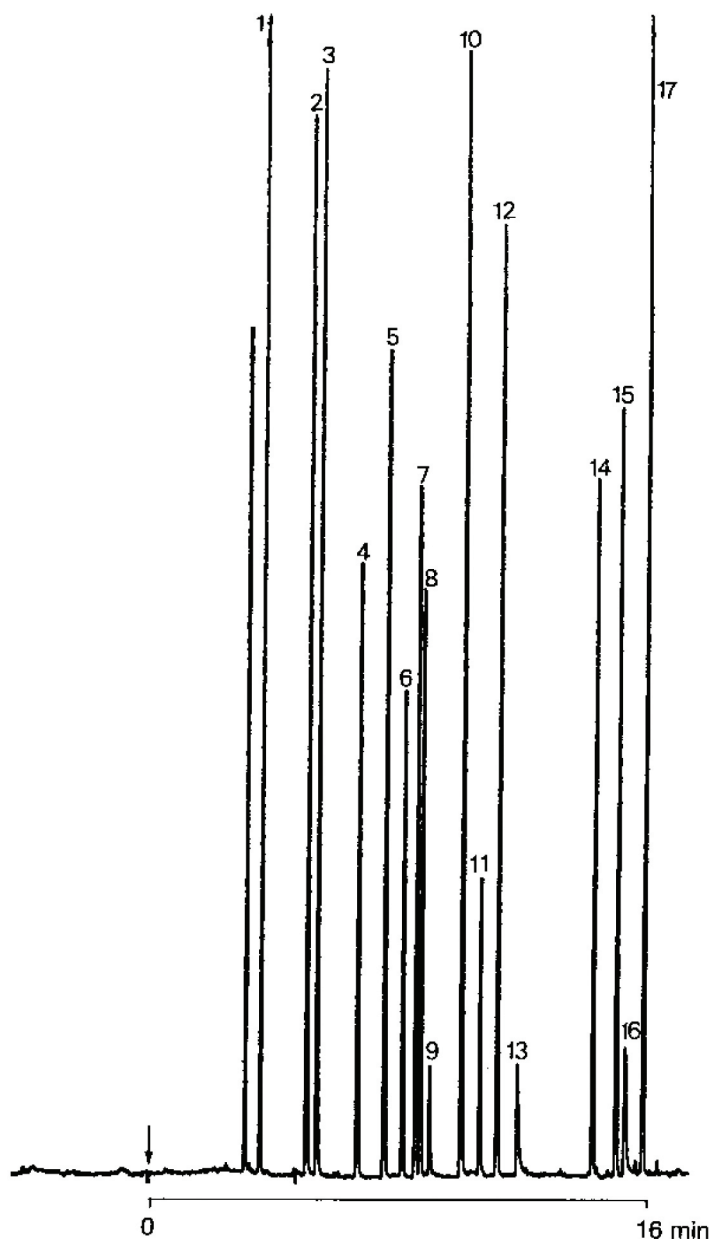
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 5 CB, 0.32 mm x 25 m fused silica
WCOT CP-Sil 5 CB (1.2 µm) (Part no. CP7760)
Temperature : 35 °C (3 min) → 220 °C, 10 °C/min
Carrier Gas : He, 55 kPa (0.55 bar, 8 psi), 18 cm/s
Injector : Splitter, 60 mL/min
T = 250 °C
Detector : FID, 4 x 10⁻¹² Afs
T = 250 °C
Sample Size : 0.1 µL
Solvent Sample : water

Peak identification

1. acetonitrile
2. butanone
3. 2-butanol
4. 1,2-dichloroethane
5. benzene
6. 1,1-dichloropropane
7. 1,2-dichloropropane
8. 2,3-dichloropropene
9. epichlorohydrin
10. methylisobutylketone
11. trans-1,3-dichloropropene
12. toluene
13. unknown
14. p-xylene
15. 1,2,3-trichloropropane
16. 2,3-dichlorohydrin
17. ethylpentylketone



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

A00119



Agilent Technologies