



Halogenated hydrocarbons, C₂

Analysis of halogenated hydrocarbons in water

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Wax 52 CB column separates three C₂ halogenated hydrocarbons in water in 40 minutes.



Agilent Technologies

Conditions

Technique : GC-PTI

Column : Agilent CP-Wax 52 CB, 0.25 mm x 60 m fused silica
WCOT CP-Wax 52 CB (df = 0.5 µm)
(Part no. CP8748)

Temperature : 30 °C (1 min) → 200 °C, 5 °C/min

Injector : TCT
Cold trap : CP-Sil 8 CB
0.53 mm; df = 5 µm
Precool temperature : -140 °C
Precool time : 3 min
Desorb time : 4 min
Desorb flow : 8.5 mL/min
Injection temp. : 220 °C
Injection time : 1 min
Backflush flow : 40 mL/min

Detector : ECD
T = 275 °C

Sample Size : 5 µL

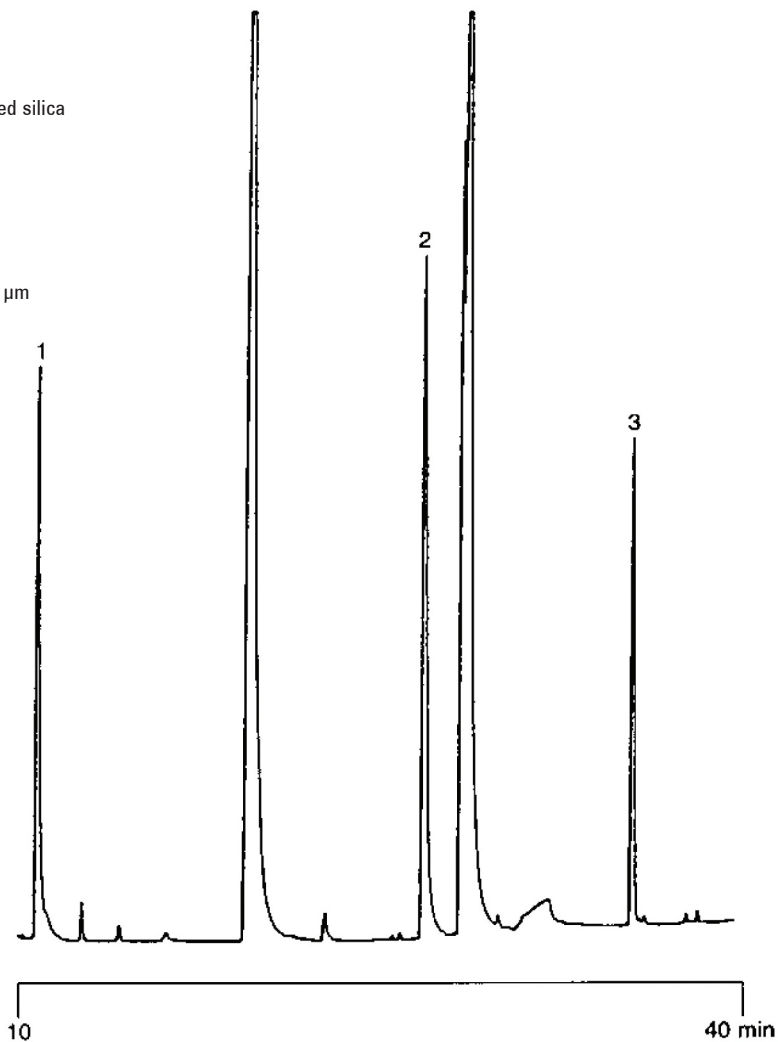
Sample Temperature : ambient

Concentration Range : 10 µg/L (10 ppb)

Courtesy : GL Sciences, Japan

Peak identification

1. 1,1,1-trichloroethane
2. trichloroethylene
3. tetrachloroethylene



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

A01018



Agilent Technologies