

Halogenated hydrocarbons, $C_1 - C_2$

Analysis of halogenated hydrocarbons in air

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 5 CB column separates eight C_1 to C_2 halogenated hydrocarbons in air in 20 minutes.



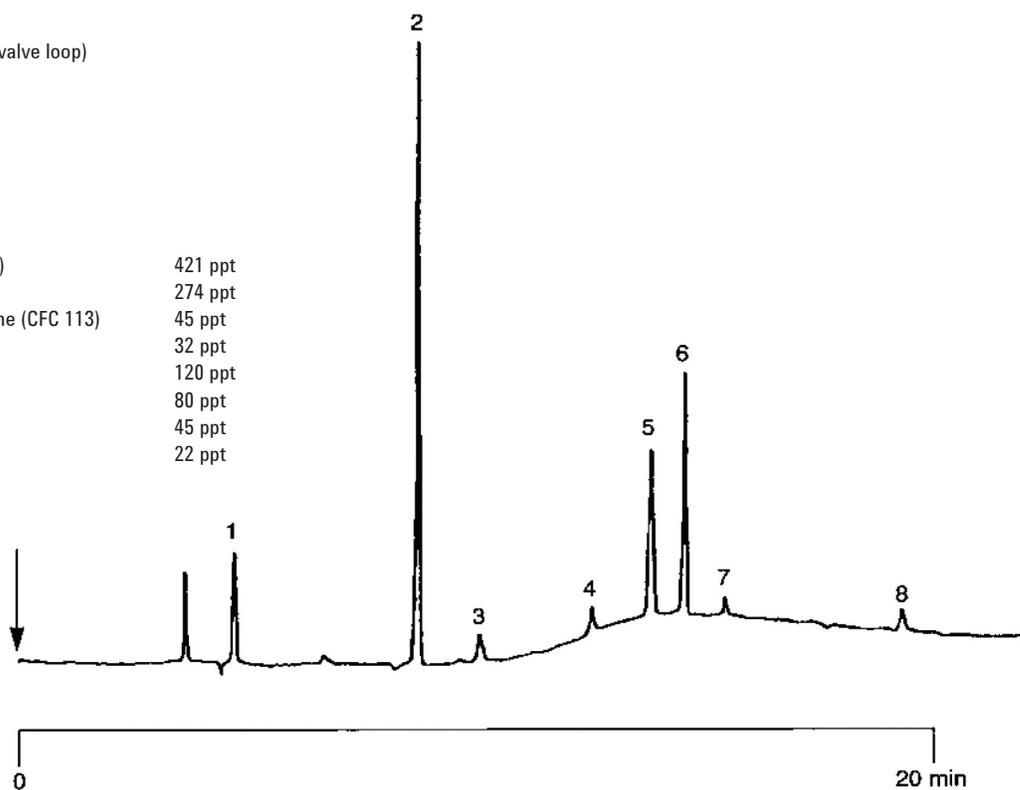
Agilent Technologies

Conditions

Technique : GC-TCT
Column : Agilent CP-Sil 5 CB, 0.32 mm x 50 m fused silica
WCOT CP-Sil 5 CB (df = 5 µm) (Part no. CP7690)
Temperature : 45 °C (3 min) → 180 °C, 10 °C/min; 180 °C (30 min)
Carrier Gas : He, 1.5 mL/min
Injector : TCT
Cold trap : Deactivated Fused-Silica
0.53 mm
Cryofocussing : -140 °C
Precool time : 15 min
Injection temp. : 100 °C
Injection time : 1 min
GC-injection block temp. : 250 °C
Detector : ECD
T = 250 °C
Sample Size : 0.5 mL (via valve loop)
Sample Temperature : ambient
Concentration Range : ppt

Peak identification

1. dichlorodifluoromethane (CFC 12)	421 ppt
2. fluorotrichloromethane (CFC 11)	274 ppt
3. 1,1,2-trifluoro-1,2,2-trichloroethane (CFC 113)	45 ppt
4. trichloromethane	32 ppt
5. 1,1,1-trichloroethane	120 ppt
6. tetrachloromethane	80 ppt
7. trichloroethylene	45 ppt
8. tetrachloroethylene	22 ppt



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

A01012



Agilent Technologies