

Halogenated hydrocarbons, $C_5 - C_{10}$

Analysis of halogenated hydrocarbons to EPA 8120

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 8 CB column separates 18 C_5 to C_{10} halogenated hydrocarbons according to EPA 8120 in 90 minutes.



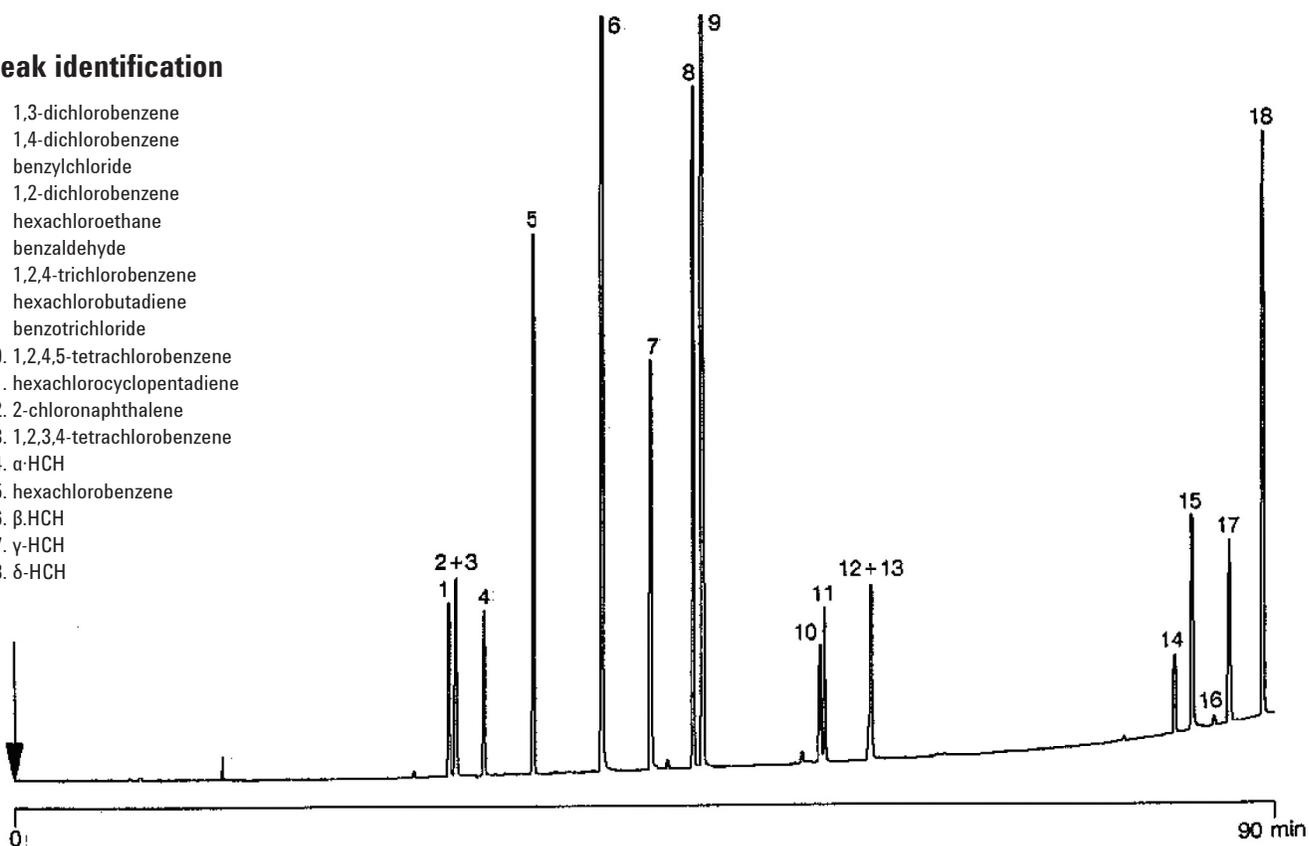
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 8 CB for pesticides, 0.32 mm x 50 m
fused silica WCOT CP-Sil 8 CB (df = 1.2 µm)
with retention gap (Part no. CP7771)
Temperature : 50 °C (2 min) → 325 °C, 3 °C/min
Carrier Gas : He, 90 kPa (0.9 bar, 13 psi)
Injector : Split, 100 mL/min
T = 265 °C
Detector : ECD
T = 300 °C

Peak identification

- 1,3-dichlorobenzene
- 1,4-dichlorobenzene
- benzylchloride
- 1,2-dichlorobenzene
- hexachloroethane
- benzaldehyde
- 1,2,4-trichlorobenzene
- hexachlorobutadiene
- benzotrichloride
- 1,2,4,5-tetrachlorobenzene
- hexachlorocyclopentadiene
- 2-chloronaphthalene
- 1,2,3,4-tetrachlorobenzene
- α-HCH
- hexachlorobenzene
- β-HCH
- γ-HCH
- δ-HCH



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

A00802



Agilent Technologies