



# Chlorinated hydrocarbons

## Analysis of chlorinated hydrocarbons to EPA 612

### Application Note

Environmental

#### Authors

Agilent Technologies, Inc.

#### Introduction

Gas chromatography with an Agilent CP-Sil 8 CB column separates nine chlorinated hydrocarbons according to EPA 612 in 48 minutes.



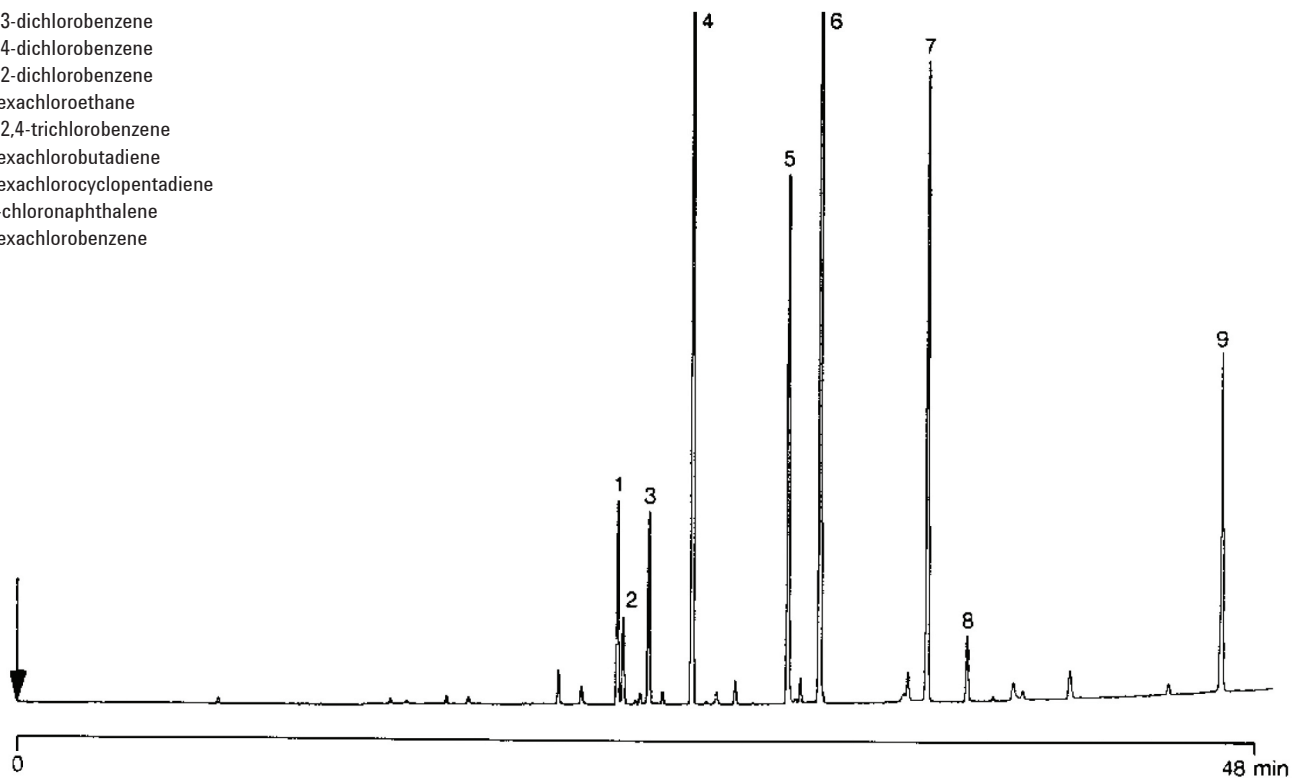
**Agilent Technologies**

## Conditions

Technique : GC-capillary  
Column : Agilent CP-Sil 8 CB, 0.32 mm x 50 m fused silica  
WCOT CP-Sil 8 CB (df = 1.2  $\mu$ m) (Part no. CP7771)  
Temperature : 50 °C (2 min)  $\rightarrow$  290 °C. 5 °C/min  
Carrier Gas : He, 90 kPa (0.9 bar, 13 psi)  
Injector : Split, 100 mL/min  
T = 265 °C  
Detector : FID  
T = 300 °C

## Peak identification

1. 1,3-dichlorobenzene
2. 1,4-dichlorobenzene
3. 1,2-dichlorobenzene
4. hexachloroethane
5. 1,2,4-trichlorobenzene
6. hexachlorobutadiene
7. hexachlorocyclopentadiene
8. 2-chloronaphthalene
9. hexachlorobenzene



[www.agilent.com/chem](http://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

A00788



**Agilent Technologies**