



# Chlorinated pesticides

## Chlorinated pesticides to EPA 505

### Application Note

Environmental

#### Authors

Agilent Technologies, Inc.

#### Introduction

Gas chromatography with an Agilent CP-Sil 8 CB for PCBs column separates 12 chlorinated pesticides according to EPA 505 in 23 minutes.



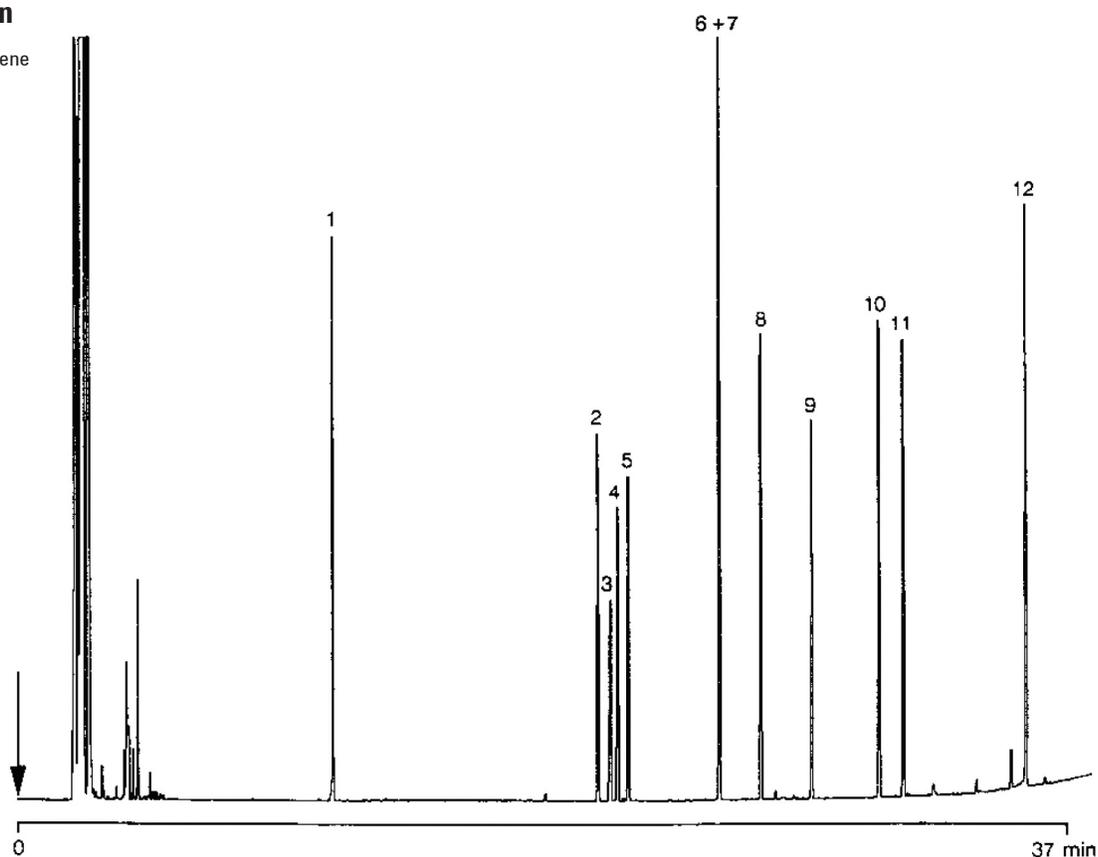
**Agilent Technologies**

## Conditions

Technique : GC-capillary  
Column : Agilent CP-Sil 8 CB for PCBs, 0.25 mm x 50 m fused silica WCOT CP-Sil 8 CB (df = 0.25 µm)  
(Part no. CP7482)  
Temperature : 100 °C (2 min) → 290 °C, 5 °C/min  
Carrier Gas : H<sub>2</sub>, 150 kPa (1.5 bar, 21 psi)  
Injector : Split, 100 mL/min  
T = 265 °C  
Detector : FID  
T = 300 °C

## Peak identification

1. hexachlorocyclopentadiene
2. hexachlorobenzene
3. simazine
4. atrazine
5. lindane
6. heptachlor
7. alachlor
8. aldrin
9. heptachlor epoxide
10. dieldrin
11. endrin
12. methoxychlor



[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

A00778



**Agilent Technologies**