

Pesticides

Analysis of chlorinated pesticides to EPA 505

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 5 CB-MS column separates ten chlorinated pesticides according to EPA 505 in 37 minutes.



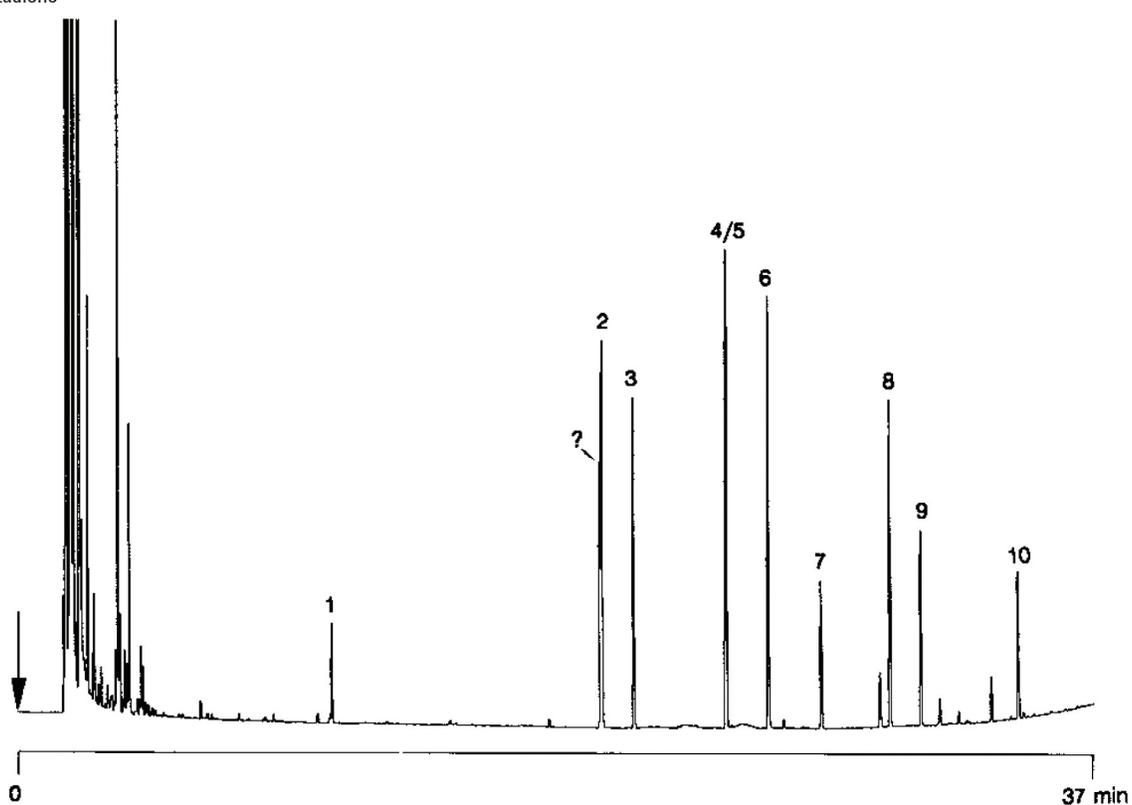
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 5 CB-MS, 0.25 mm x 50 m fused silica
WCOT CP-Sil 5 CB-MS (df = 0.25 µm)
(Part no. CP7843)
Temperature : 100 °C (2 min) → 290 °C, 5 °C/min
Carrier Gas : H₂, 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. lindane
4. heptachlor
5. alachlor
6. aldrin
7. heptachlor epoxide
8. dieldrin
9. endrin
10. methoxychlor



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

A00913



Agilent Technologies