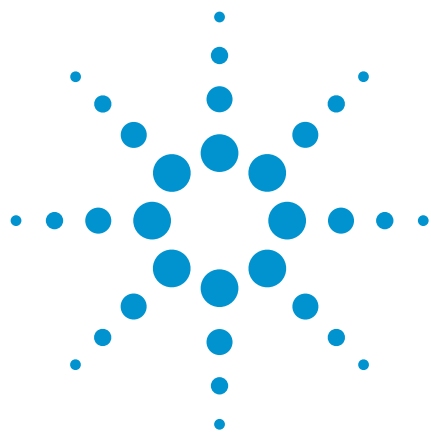




Organochlorine pesticides

Analysis of organochlorine pesticides according to EPA 8080

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 8 CB for Pesticides column separates 19 organochlorine pesticides according to EPA 8080 in 50 minutes.



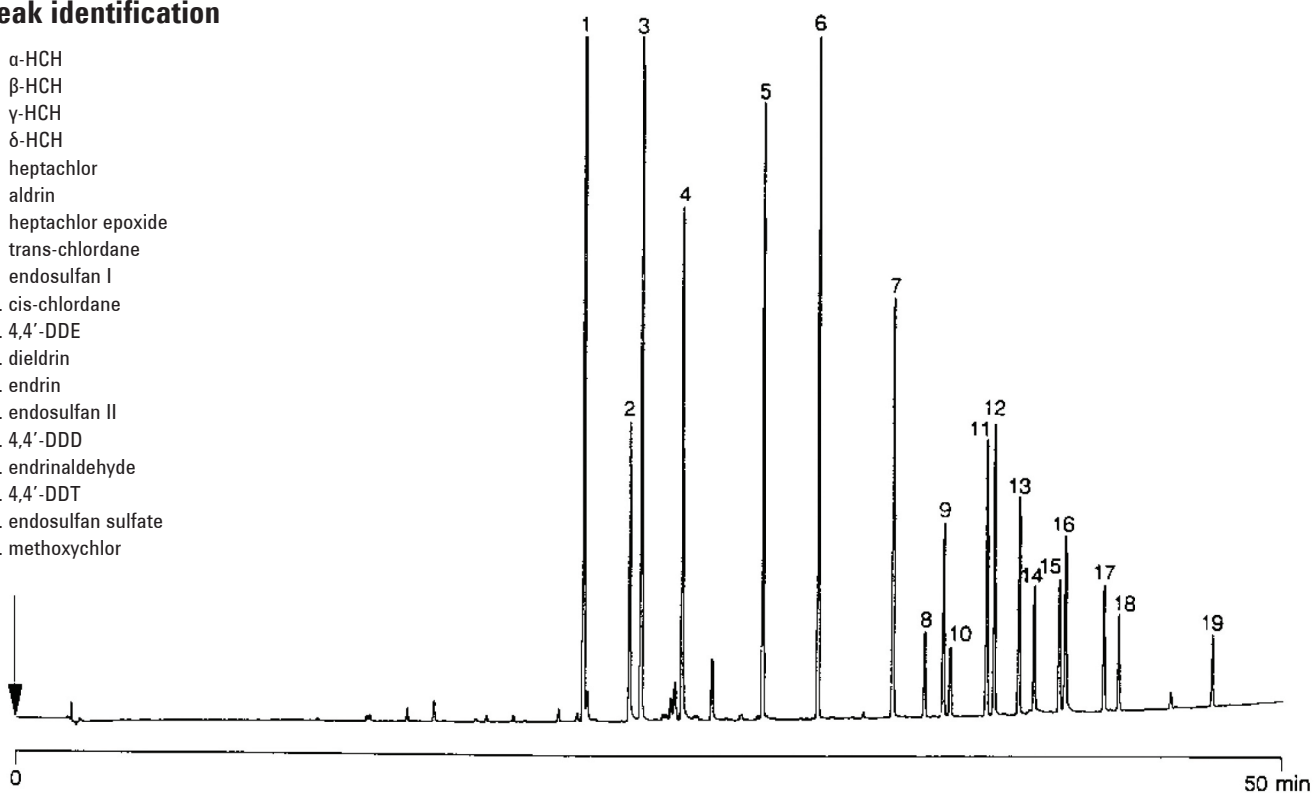
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 8 CB for Pesticides, 0.25 mm x 50 m
fused silica WCOT CP-Sil 8 CB (df = 0.12 µm)
(Part no. CP7481)
Temperature : 100 °C (2 min) → 240 °C, 3 °C/min
Carrier Gas : H₂, 140 kPa (1.4 bar, 20 psi)
Injector : Split, 100 mL/min
T = 265 °C
Detector : ECD
T = 310 °C

Peak identification

1. α-HCH
2. β-HCH
3. γ-HCH
4. δ-HCH
5. heptachlor
6. aldrin
7. heptachlor epoxide
8. trans-chlordane
9. endosulfan I
10. cis-chlordane
11. 4,4'-DDE
12. dieldrin
13. endrin
14. endosulfan II
15. 4,4'-DDD
16. endrinaldehyde
17. 4,4'-DDT
18. endosulfan sulfate
19. 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

A00796



Agilent Technologies