



Pesticides

Analysis of organochlorine in pesticides in water

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 19 CB column separates 30 organochlorine pesticides in 59 minutes.



Agilent Technologies

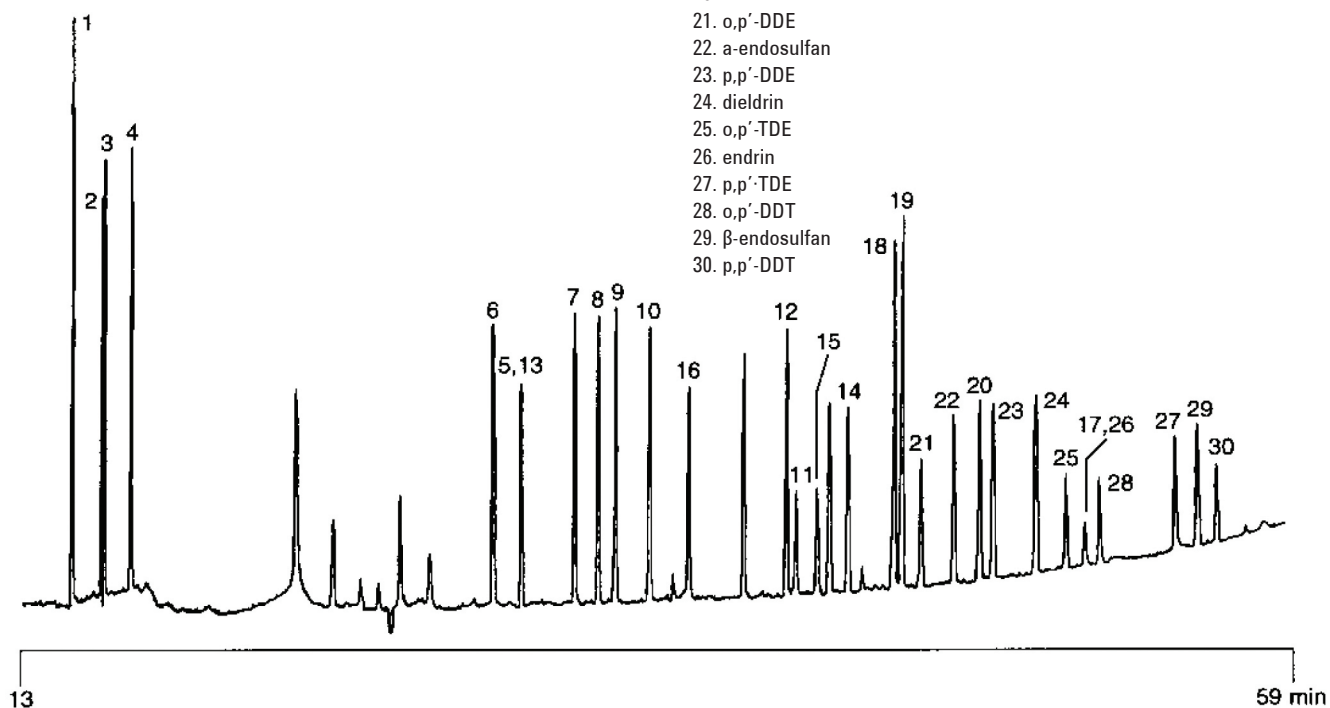
Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 19 CB for pesticides, 0.25 mm x 50 m
fused silica WCOT CP-Sil 19 CB (df = 0.2 µm)
(Part no. CP7407)
Temperature : 45 °C (4 min) → 150 °C, 15 °C/min
Carrier Gas : He
Injector : On-column
Detector : ECD
T = 300 °C

Courtesy : Mr W. Civil, NRA, Exeter, UK

Peak identification

1. 1,3,5-TCB
2. HCB
3. 1,2,4-TCB
4. 1,2,3-TCB
5. trifluralin
6. HCB
7. α-HCB
8. pentachloronitrobenzene (I.S.)
9. triallale
10. γ-HCH
11. β-HCH
12. vinclozolin
13. heptachlor
14. δ-HCH
15. ε-HCH
16. aldrin
17. isodrin
18. cis-HCE
19. trans-HCE
20. cis-chlordane
21. o,p'-DDE
22. a-endosulfan
23. p,p'-DDE
24. dieldrin
25. o,p'-TDE
26. endrin
27. p,p'-TDE
28. o,p'-DDT
29. β-endosulfan
30. p,p'-DDT



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

A01094



Agilent Technologies