



Pesticides and PCBs

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 8 CB for Pesticides column separates 60 pesticides and polycyclic biphenyls in 24 minutes.



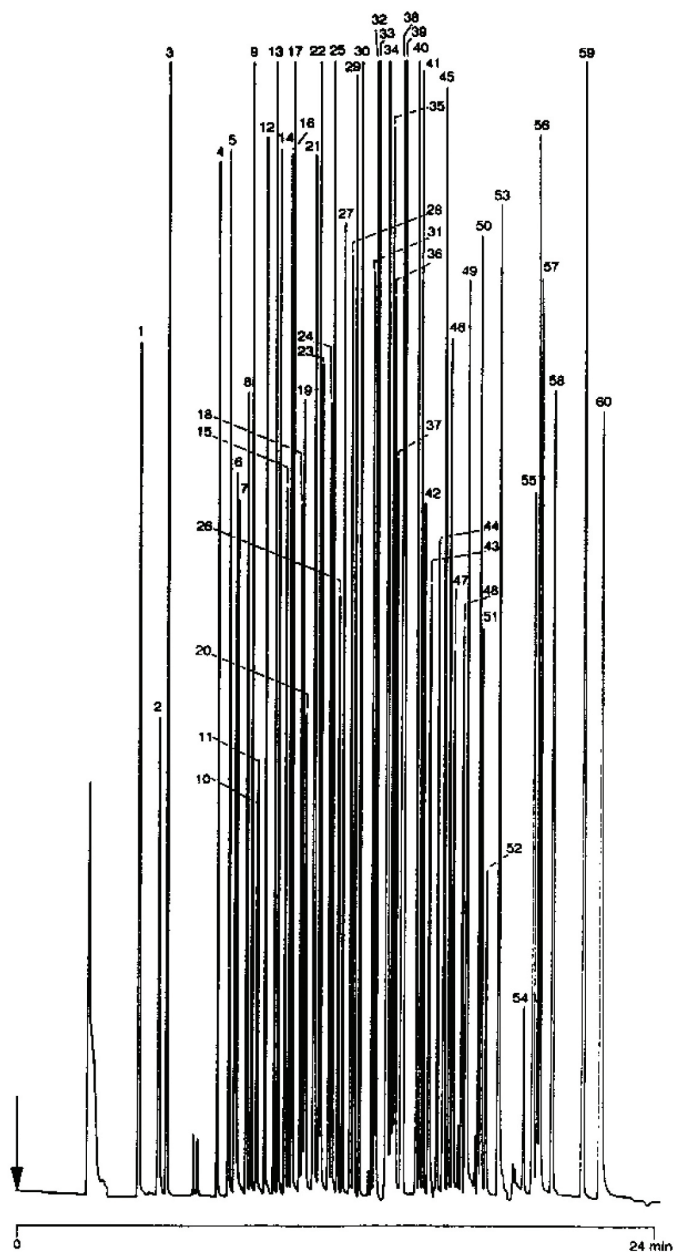
Agilent Technologies

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 : 80 °C → 220 °C, 20 °C/min;
 220 °C → 270 °C, 4 °C/min;
 Carrier Gas : H₂, 125 kPa (1.25 bar, 17.9 psi)
 Injector : On-column
 Detector : ECD
 T = 300 °C
 Sample Size : 1 µL
 Concentration range : 0.1 - 2 ng per compound
 Courtesy : Mr. Lembacher, HIPPO KG, Pfaffenhofen, Germany

Peak identification

- | | |
|-------------------------------------|-----------------------------------|
| 1. hexachloroethane | 31. octachlorostyrene |
| 2. 1,2,4-trichlorobenzene | 32. HEO |
| 3. hexachlorobutadiene | 33. modane (danthron) |
| 4. etridiazole | 34. chloroflurene |
| 5. 1,2,4-trichloro-5-nitrobenzene | 35. trans-chlordane |
| 6. α-1,2,4-trichloroacetophenone | 36. o,p'-DDE |
| 7. trichloroaniline | 37. PCB 101 |
| 8. chlorophenprop-methyl | 38. ε-endosulfan |
| 9. techazene | 39. cis-chlordane |
| 10. 2,3,5,6-tetrachloroaniline | 40. p,p'-DDE |
| 11. unknown | 41. dieldrin |
| 12. 2,3,4,5-tetrachloronitrobenzene | 42. o,p'-DDD |
| 13. α-HCH | 43. dibromoacetophenone |
| 14. HCB | 44. endrin |
| 15. β-HCH | 45. β-endosulfan |
| 16. lindane (γ-HCH) | 46. p,p'-DDD |
| 17. quintozone | 47. o,p'-DDT |
| 18. δ-HCH | 48. PCB 153 |
| 19. chlorothalonil | 49. kepone |
| 20. ε-HCH | 50. endosulfansulfate |
| 21. pentachloroaniline | 51. p,p'-DDT |
| 22. PCB 28 | 52. PCB 138 |
| 23. vinclozolin | 53. fluroxypry-1-methylhexylester |
| 24. heptachlor epoxide | 54. unknown |
| 25. S-421 | 55. hexabromobenzene |
| 26. PCB 52 | 56. methoxychlor |
| 27. dichlofluanid | 57. fenprothrin |
| 28. aldrin | 58. PCB 180 |
| 29. chlorothalonil | 59. mirex |
| 30. telofrin (isobenzan) | 60. fenarimol |



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

A00709



Agilent Technologies