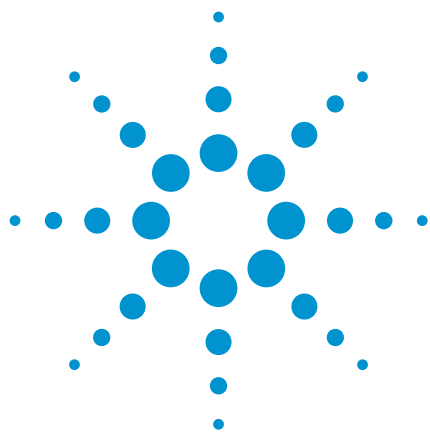




Pesticides and PCBs

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 8 CB column separates 23 pesticides and polychlorinated biphenyls in a standard mixture in 26 minutes.



Agilent Technologies

Conditions

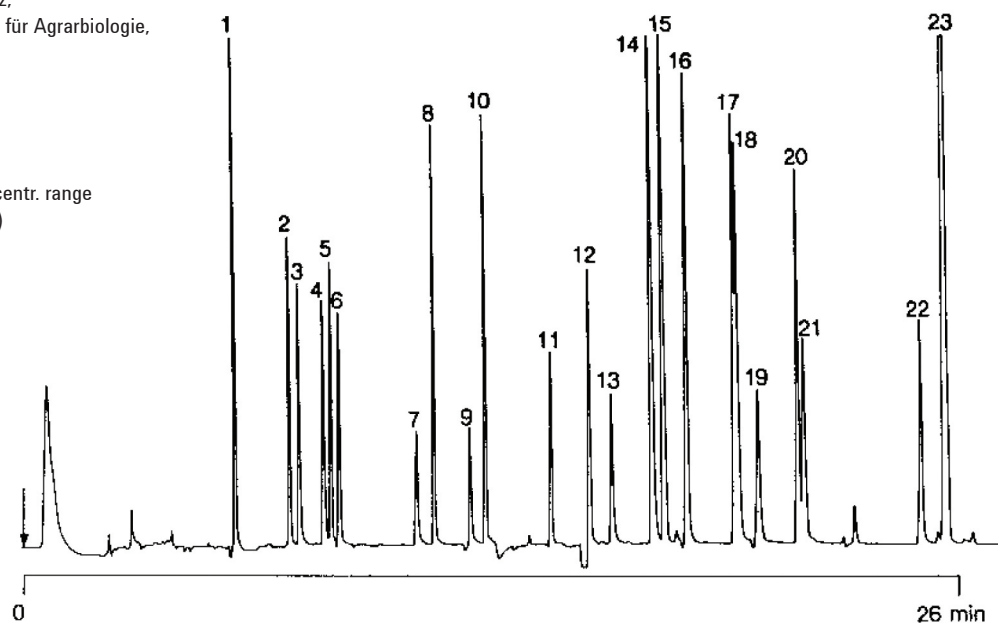
Technique : GC-capillary
Column : Agilent CP-Sil 8 CB, 0.32 mm x 50 m fused silica
WCOT CP-Sil 8 CB (df = 0.12 µm) (Part no. CP7751)
Temperature : 140 °C (2.5 min) → 244 °C,
4 °C/min (0 min) → 290 °C, 20 °C/min (3 min)
Carrier Gas : He, 140 kPa (1.4 bar, 20 psi)
Injector : Splitless (140 sec)
T = 210 °C
Detector : ECD,
T = 300 °C
Sample Size : 5.0 µL
Solvent Sample : n-decane

Courtesy : Ing. W Brodacz,
Bundesanstalt für Agrarbiologie,
Linz, Austria

Peak Identification :

- | | |
|-------------------------|----|
| 1. technazene (TCNB) | 10 |
| 2. α-HCH | 5 |
| 3. HCB | 5 |
| 4. β-HCH | 10 |
| 5. lindane | 5 |
| 8. quintozone (PCNB) | 5 |
| 7. PCB 28 | 10 |
| 8. heptachlor | 10 |
| 9. PCB 52 | 10 |
| 10. aldrin | 10 |
| 11. β-heptachlorepoxide | 5 |
| 12. γ-chlordane | 10 |
| 13. PCB 101 | 10 |
| 14. dieldrin | 20 |
| 15. p,p'-DDE | 20 |
| 16. endrin | 20 |
| 17. p,p'-DDD (TDE) | 20 |
| 18. o,p'-DDT | 20 |
| 19. PCB 153 | 10 |
| 20. p,p'-DDT | 20 |
| 21. PCB 138 | 10 |
| 22. PCB 180 | 10 |
| 23. mirex | 60 |

Concentr. range
(ppb)



www.agilent.com/chem

This information is subject to change without notice.

© Agilent Technologies, Inc. 2011

Printed in the USA

31 October, 2011

Published in UK, August 03, 2010

A00478



Agilent Technologies