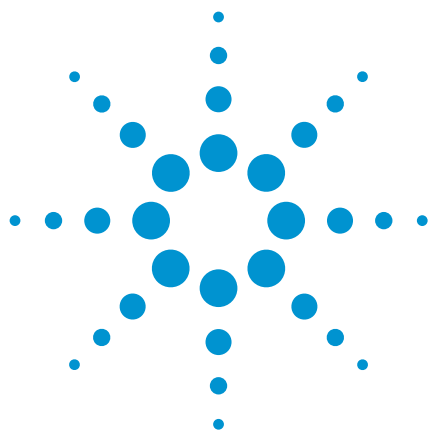




Pesticides

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

GC/MS with an Agilent CP-Sil 8 CB column separates 15 organochlorine pesticides in 29 minutes.



Agilent Technologies

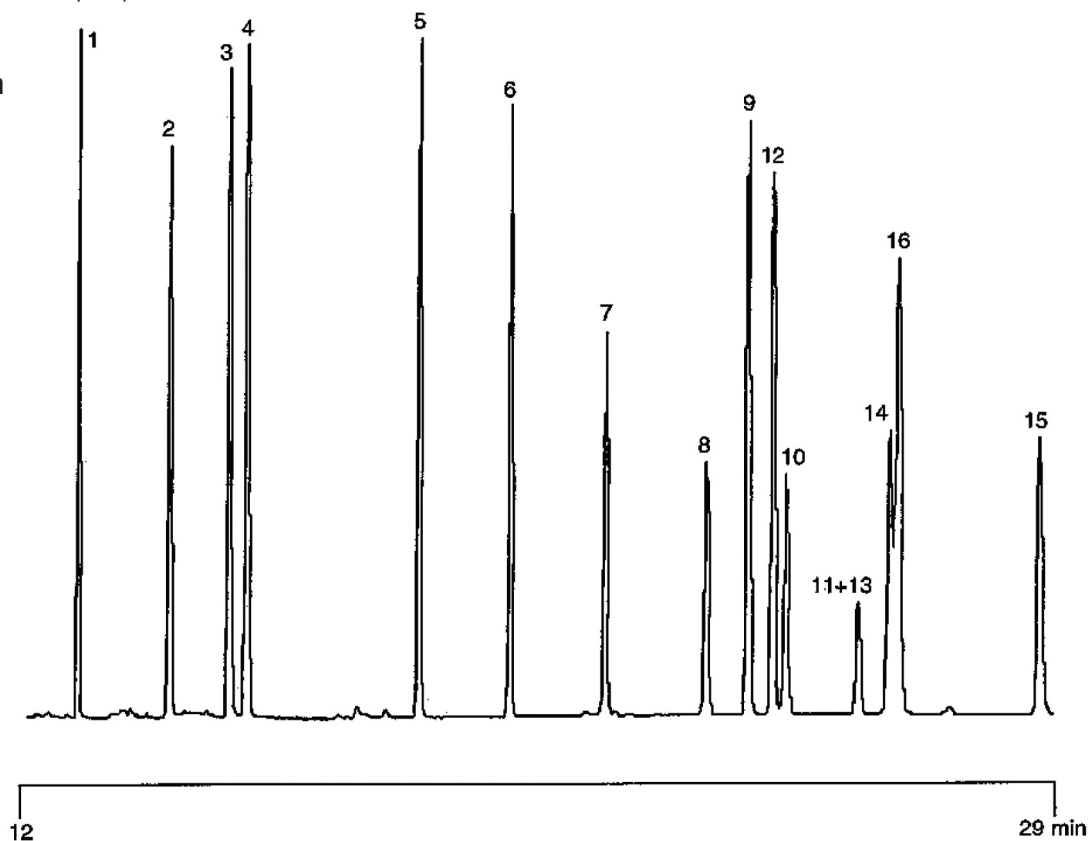
Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 8 CB Low Bleed/MS, 0.25 mm x 30 m
fused silica WCOT CP-Sil 8 CB (df = 0.5 μ m)
(Part no. CP5871)
Temperature : 50 °C $\rightarrow$ 170 °C, 25 °C/min;
170 °C $\rightarrow$ 250 °C, 4 °C/min
Carrier Gas : He
Injector : Splitless,
T = 250 °C
Detector : MS in EI mode
Injection Volume : 5 μ L

Courtesy : Mr I. Epson, North East Scientific Services,
Newcastle Upon Tyne, UK

Peak identification

1. trifluralin
2. α -HCH
3. β -HCH
4. γ -HCH
5. heptachlor
6. aldrin
7. heptachlor epoxide
8. α -endosulfan
9. p,p'-DDE
10. dieldrin
11. endrin
12. o,p'-DDD
13. β -endosulfan
14. p,p'-DDD
15. p,p'-DDT
16. endrin aldehyde



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

A01109



Agilent Technologies