



# **Pesticides**

## **Analysis of pesticides to EPA 508**

### **Application Note**

Environmental

#### **Authors**

Agilent Technologies, Inc.

#### **Introduction**

Gas chromatography with an Agilent CP-Sil 8 CB for Pesticides column separates 29 pesticides according to EPA 508 in 64 minutes.



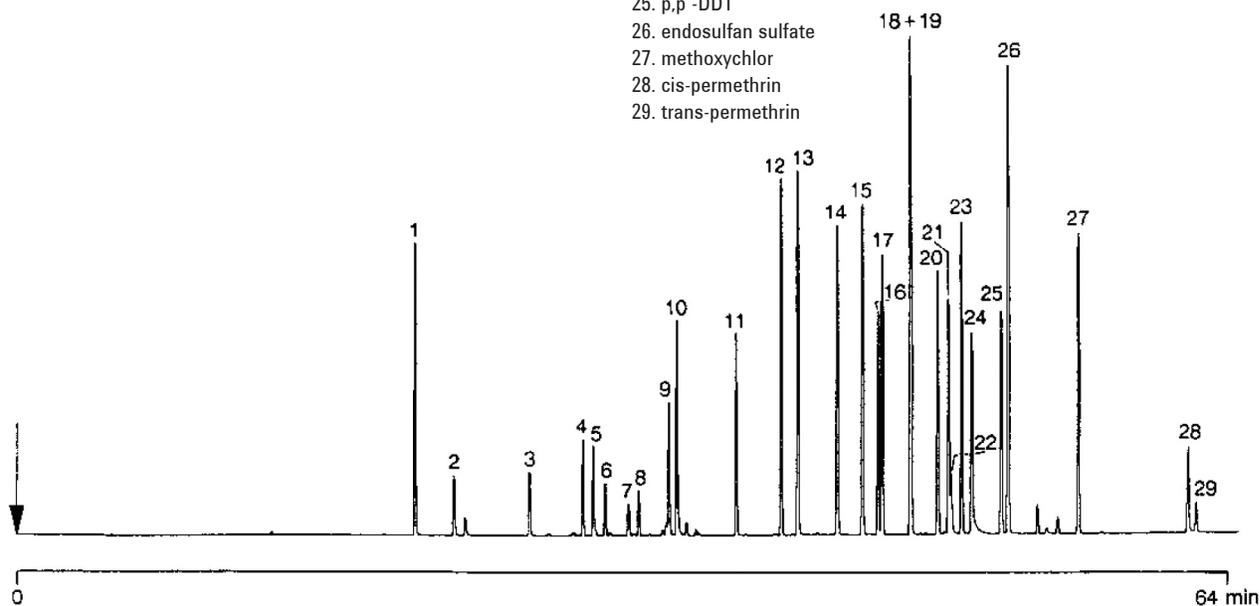
**Agilent Technologies**

## 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) with retention gap (Part no. CP7481)  
Temperature : 100 °C (2 min) → 300 °C, 3 °C/min  
Carrier Gas : He, 150 kPa (1.5 bar, 21 psi)  
Injector : Split, 100 mL/min  
T = 270 °C  
Detector : ECD  
T = 325 °C

## Peak identification

1. etridiazole
2. chlorneb
3. propachlor
4. trifluralin
5. α-HCH
6. hexachlorobenzene
7. β-HCH
8. γ-HCH (lindane)
9. δ-HCH
10. chlorothalonil
11. heptachlor
12. aldrin
13. DCPA
14. heptachlorepoxyde
15. trans-chlordane
16. endosulfan I
17. cis-chlordane
18. p,p'-DDE
19. dieldrin
20. endrin
21. endosulfan II
22. chlorobenzitate
23. p,p'-DDD
24. endrinaldehyde
25. p,p'-DDT
26. endosulfan sulfate
27. methoxychlor
28. cis-permethrin
29. trans-permethrin



[www.agilent.com/chem](http://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

A00780



**Agilent Technologies**