



## **Pesticides and herbicides**

# Analysis of pesticides and herbicides to EPA 507

## Application Note

Environmental

### **Authors**

Agilent Technologies, Inc.

### **Introduction**

Gas chromatography with an Agilent CP-Sil 8 CB Low Bleed/MS column separates 19 pesticides and herbicides according to EPA 507 in 28 minutes.



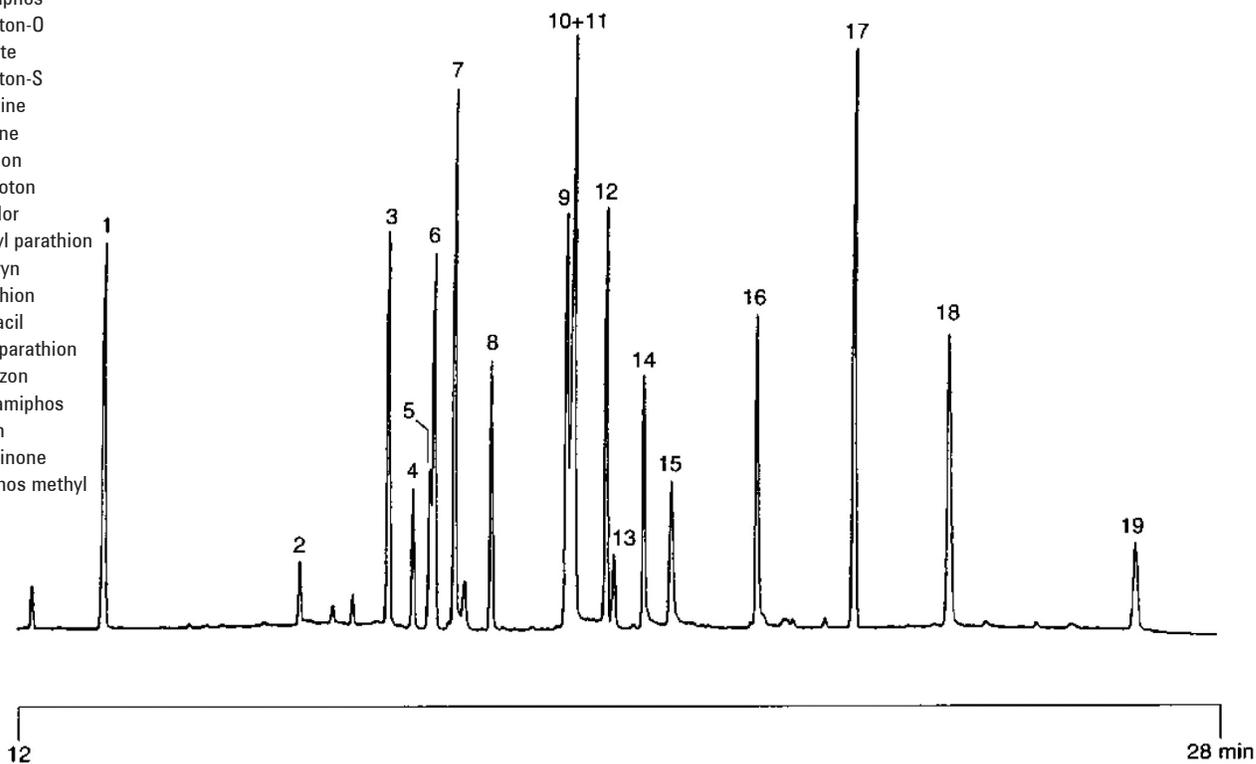
**Agilent Technologies**

## Conditions

Technique : GC-capillary  
Column : Agilent CP-Sil 8 CB low bleed/MS, 0.25 mm x 50 m  
fused silica (df = 0.25 µm) (Part no. CP5843)  
Temperature : 100 °C (2 min) → 300 °C, 10 °C/min  
Carrier Gas : He, 100 kPa (1 bar, 14 psi)  
Injector : Split, 100 mL/min  
T = 250 °C  
Detector : MSD  
T = 280 °C

## Peak identification

1. mevinphos
2. demeton-O
3. phorate
4. demeton-S
5. simazine
6. atrazine
7. diazinon
8. disulfoton
9. alachlor
10. methyl parathion
11. ametryn
12. malathion
13. bromacil
14. ethyl parathion
15. bentazon
16. phenamiphos
17. ethion
18. hexazinone
19. azinphos methyl



[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

A01002



**Agilent Technologies**