

## **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 5 CB-MS column separates 19 pesticides and herbicides according to EPA 507 in 37 minutes.



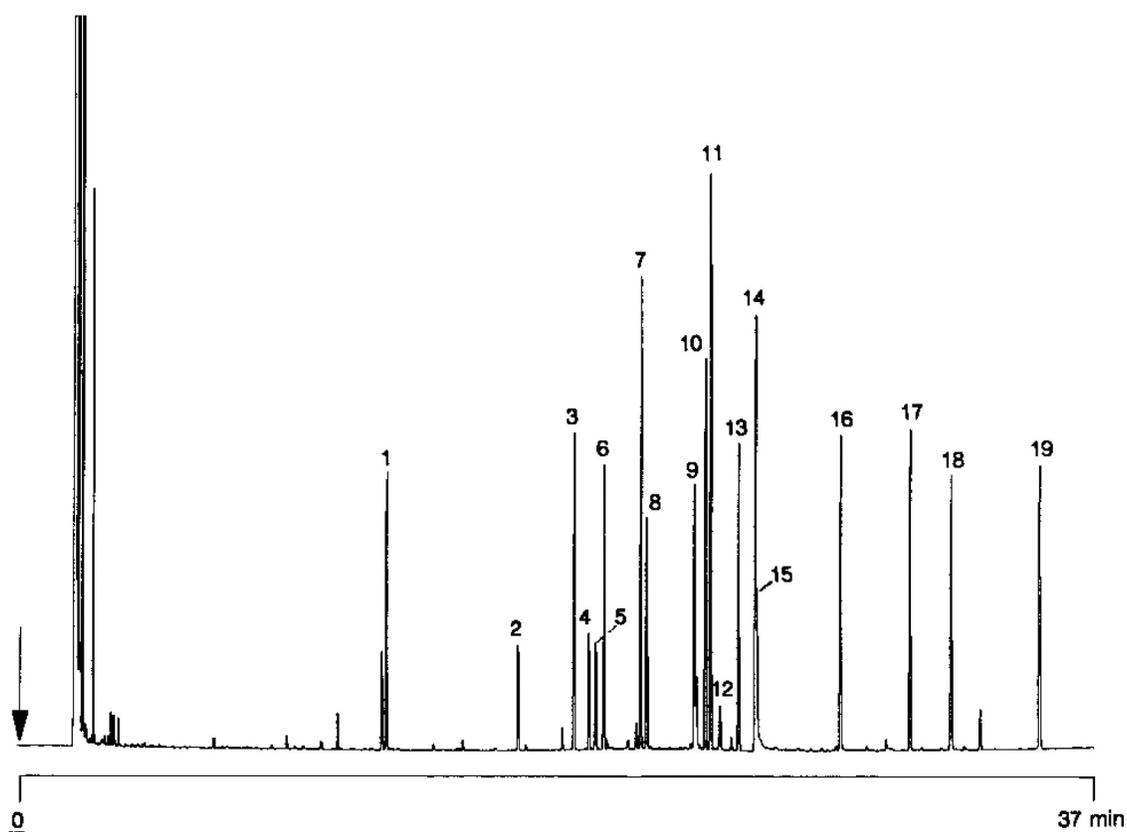
**Agilent Technologies**

## Conditions

Technique : GC-capillary  
Column : Agilent CP-Sil 5 CB-MS, 0.25 mm x 50 m fused silica  
WCOT CP-Sil 5 CB (df = 0.25 µm) (Part no. CP7843)  
Temperature : 100 °C (2 min) → 270 °C, 5 °C/min  
Carrier Gas : H<sub>2</sub>, 120 kPa (1.2 bar, 17.1 psi)  
Injector : Split, 100 mL/min  
T = 275 °C  
Detector : FID  
T = 300 °C

## Peak identification

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

A00911



**Agilent Technologies**