

# Application Report 405

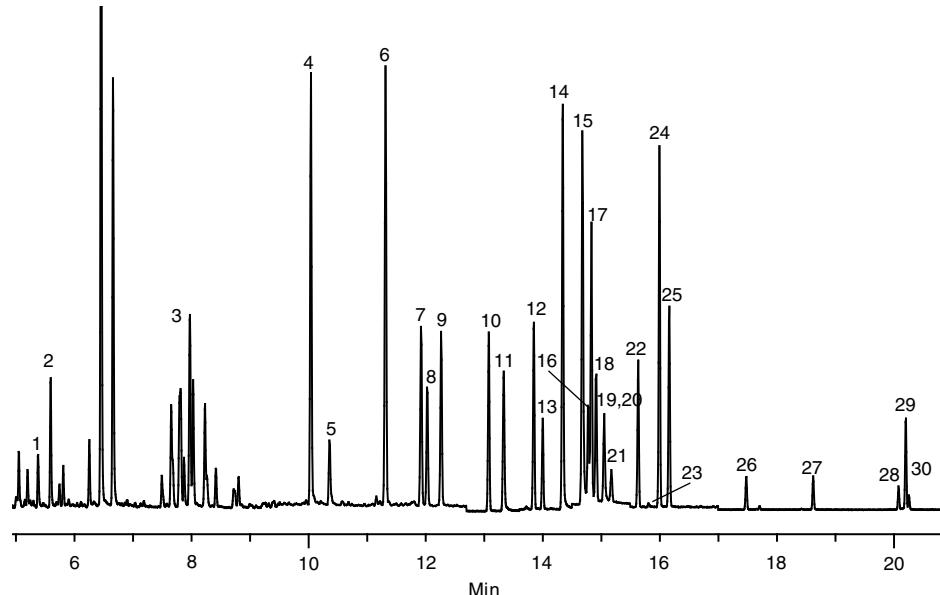
## GC-MS Analysis Agricultural Pesticides on the SLB-5ms

This application illustrates the analysis of 29 different agricultural pesticides of different classifications and functionalities. The analysis was done by GC-MS using selected ion monitoring (SIM) to allow detection of all compounds at 100 µg/L. The standard solution was made in toluene to increase its stability and enhance response of some of the pesticides. The extra peaks in the first portion of the run are due to interferences from the solvent.

**Author:** Katherine Stenerson  
**Acquisition System:** 5973 GC-MS  
**Notebook Reference:** 1569-037

### Key Words

chlorinated pesticides, organophosphorus pesticides, carbamate pesticides, agricultural pesticides, SLB-5ms, 28471-U



G003595

### Conditions

column: SLB-5ms, 30 m x 0.25 mm I.D., 0.25 µm (28471-U)  
oven: 100 °C (1 min.), 10 °C/min. to 300 °C (5 min.)  
inj.: 250 °C  
MSD interface: 300 °C  
scan range: selected ion monitoring (SIM)  
carrier gas: helium, 1 mL/min. constant flow  
injection: 1 µL, pulsed (20 psi until 0.20 min.), splitless (1.0 min.)  
liner: 4 mm I.D., single taper  
sample: 100 ppb pesticide mixture in toluene

### Peak IDs

1. Methamidiphos
2. Dichlorvos
3. Acephate
4. Propoxur
5. Ethoprophos (I.S.)
6. Hexachlorobenzene
7. γ-BHC
8. Diazinon
9. Chlorothalonil
10. Chlorpyrifos-methyl
11. Carbaryl
12. Dichlofluanid
13. Chlormpyrifos
14. p-Dichlorobenzophenone
15. Cyprodinil
16. Pencyanazole
17. Tolyfluanid
18. Heptachlor epoxide
19. Captan
20. Thia'bendazole
21. Folpet
22. cis-Chlordane
23. Imazalil
24. 4,4'-DDE
25. Dieldrin
26. Endosulfan sulfate
27. Dicofol
28. cis-Permethrin
29. trans-Permethrin
30. Coumaphos