

# Application Report 388

## US EPA Method 608, 8081, OLM04.2 Organochlorine Pesticides on the SPB-608

This is the analysis of a 22-component standard containing 20 pesticides and 2 surrogate compounds commonly analyzed by US EPA Method 8081. The run conditions shown in this application were also used in a separate application to separate the same compounds on a 30 m x 0.25 mm I.D. x 0.25  $\mu$ m SLB-5ms. Between both the SPB-608 and SLB-5ms, all 22 compounds were resolved, making the two columns a possible combination for the dual column analysis of pesticides by US EPA Method 8081.

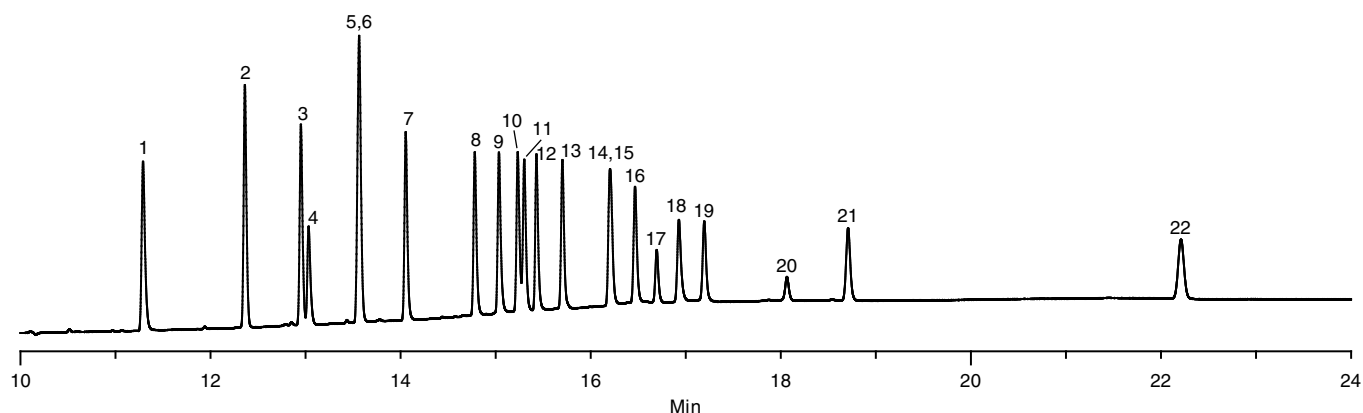
### Key Words

organochlorine pesticides, US EPA Method 8081, SPB-608, 46845-U, 24103-U

Author: Katherine Stenerson

Acquisition System: 6249

Notebook Reference: 1569-016



G003530

### Conditions

column: SPB-608, 30 m x 0.25 mm I.D., 0.25  $\mu$ m (24103-U)  
oven: 100 °C (2 min.), 15 °C/min. to 300 °C (5 min.)  
inj.: 250 °C  
det.: ECD, 300 °C  
carrier gas: helium, 0.9 mL/min, constant flow  
injection: 2.0  $\mu$ L, splitless (0.75 min.)  
liner: 4 mm I.D., single taper  
sample: chlorinated pesticide standard (46845-U), diluted to 50 ppb in n-hexane

### Peak IDs

1. 2,4,5,6-tetrachloro-m-xylene (surr.)	12. 4,4'-DDE
2. $\alpha$ -BHC	13. Dieldrin
3. $\gamma$ -BHC	14. Endrin
4. $\beta$ -BHC	15. 4,4'-DDD
5. $\delta$ -BHC	16. Endosulfan II
6. Heptachlor	17. 4,4'-DDT
7. Aldrin	18. Endrin aldehyde
8. Heptachlor epoxide	19. Endosulfan sulfate
9. $\gamma$ -chlordane	20. Methoxychlor
10. $\alpha$ -chlordane	21. Endrin ketone
11. Endosulfan I	22. Decachlorobiphenyl (surr.)