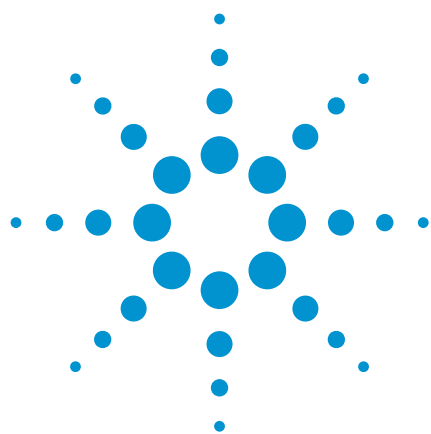




Organophosphorus pesticides

Analysis of organophosphorus pesticides with thioether groups

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 13 CB column separates 17 organophosphorus pesticides with thioether groups in 24 minutes.



Agilent Technologies

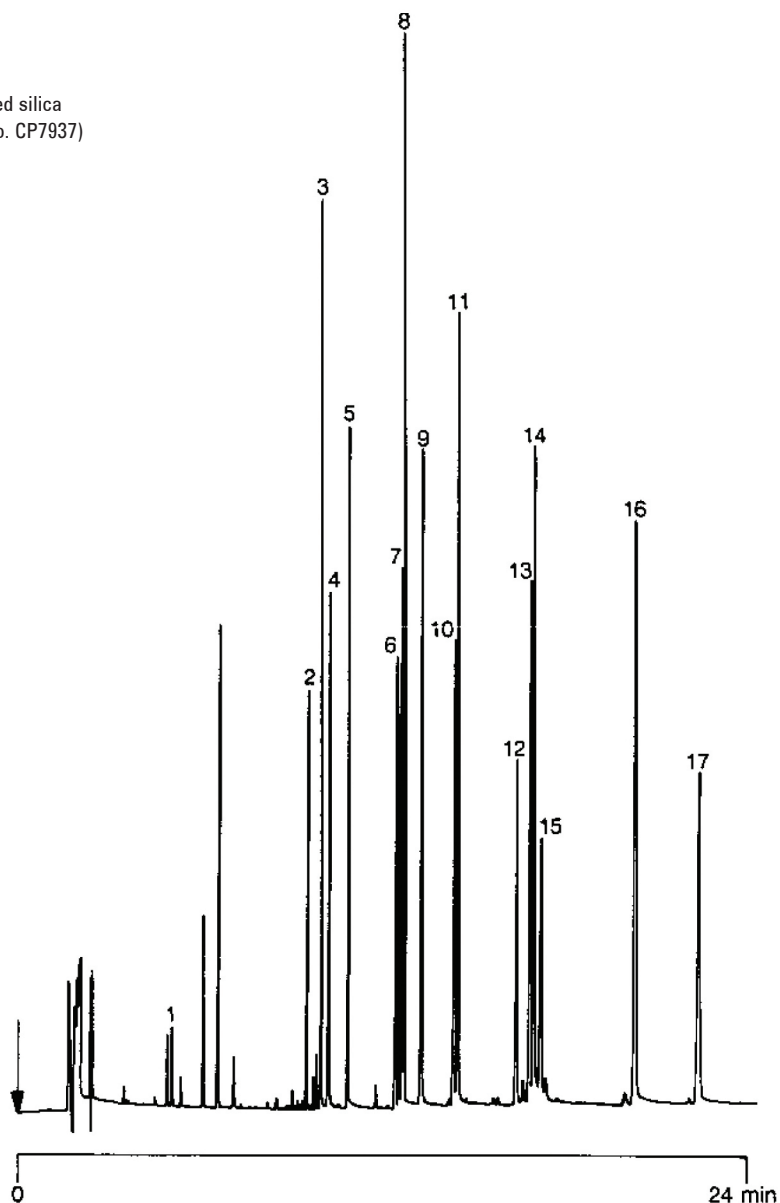
Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 13 CB, 0.32 mm x 50 m fused silica
WCOT CP-Sil 13 CB (df = 0.4 µm) (Part no. CP7937)
Temperature : 80 °C → 270 °C, 20 °C/min
Carrier Gas : H₂, 100 kPa (1.0 bar, 14 psi)
Injector : On-column
Detector : NPD
T = 300 °C
Sample Size : 1 µL
Concentration Range : 0.25 - 1 ng/µL

Courtesy : Mr Lembacher,
Hipp K. G., Pfaffenhofen,
Germany

Peak identification

	Concentration range (ng/µL)
1. triethyl phosphate	0.25
2. demeton-S-methyl	0.5
3. phorate	0.5
4. demeton	0.5
5. disulfoton	0.5
6. demeton-S-methylsulfon	1.0
7. fenthion	0.5
8. phorate-sulfon	1.0
9. demeton-sulfon	1.0
10. fenamiphos	0.5
11. disulfoton-sulfon	1.0
12. fensulfothion	0.5
13. fenthion-sulfon	1.0
14. fenamiphos-sulfon	1.0
15. carbophenothion	0.5
16. fensulfothion-sulfon	1.0
17. carbophenothion-sulfon	1.0



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

A00614



Agilent Technologies