



Pesticides

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

GC/MS with an Agilent CP-Sil 8 CB column separates 14 organophosphorus pesticides in 27 minutes.



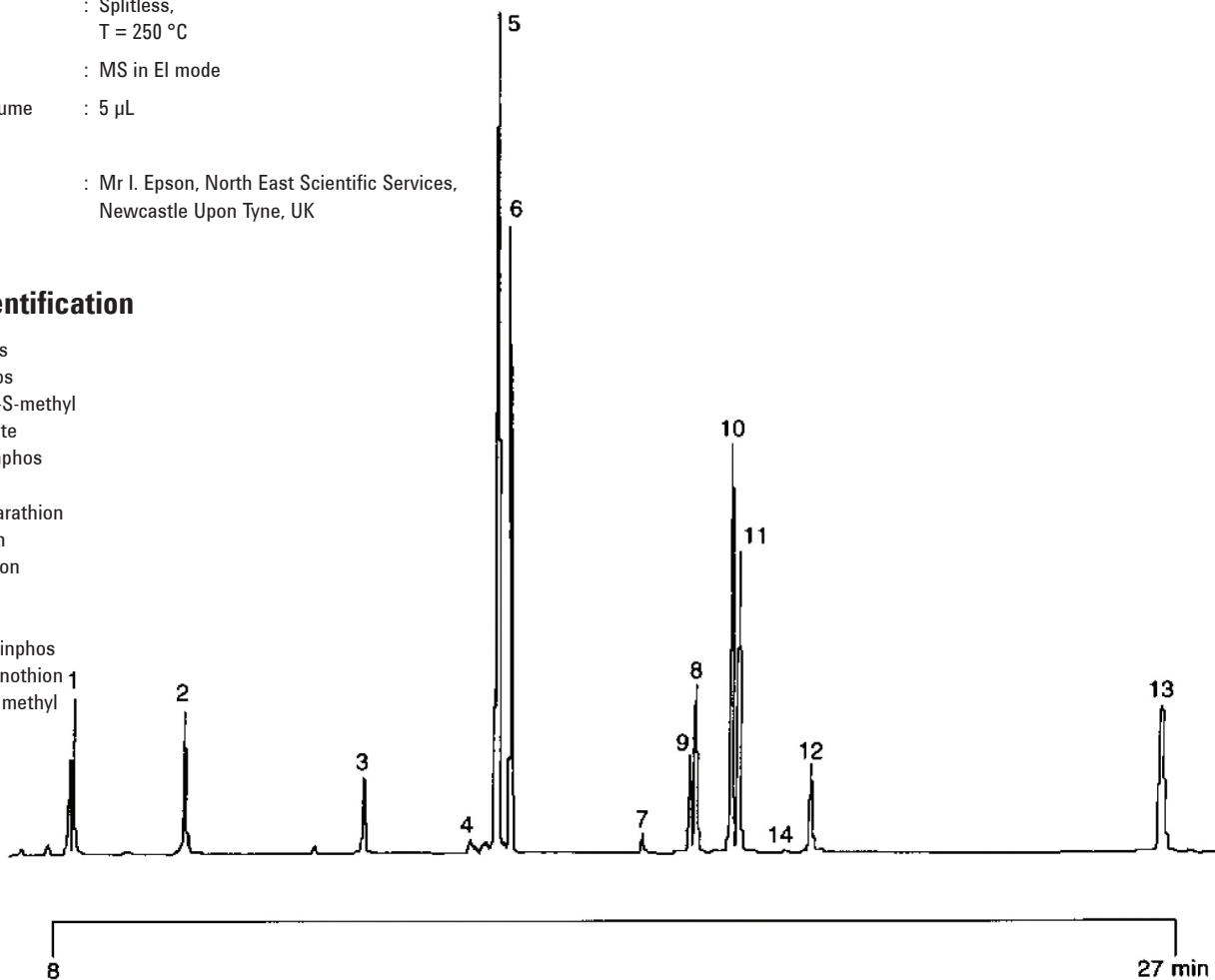
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 8 CB low bleed/MS, 0.25 mm x 30 m
fused silica WCOT CP-Sil 8 CB (df = 0.5 µm)
(Part no. CP5871)
Temperature : 50 °C → 170 °C, 25 °C/min;
170 °C → 250 °C, 4 °C/min
Carrier Gas : He
Injector : Splitless,
T = 250 °C
Detector : MS in EI mode
Injection Volume : 5 µL
Courtesy : Mr I. Epson, North East Scientific Services,
Newcastle Upon Tyne, UK

Peak identification

1. dichlorvos
2. mevinphos
3. demeton-S-methyl
4. dimethoate
5. propetamphos
6. diazinon
7. methyl parathion
8. malathion
9. fenitrothion
10. fenthion
11. parathion
12. chlorfenvinphos
13. carbophenothion
14. azinphos methyl



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

A01108



Agilent Technologies