



Pesticides

Application Note

Food Testing & Agriculture

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with a custom-made Agilent CP-Sil 5 column separates 11 pesticides in 17 minutes.



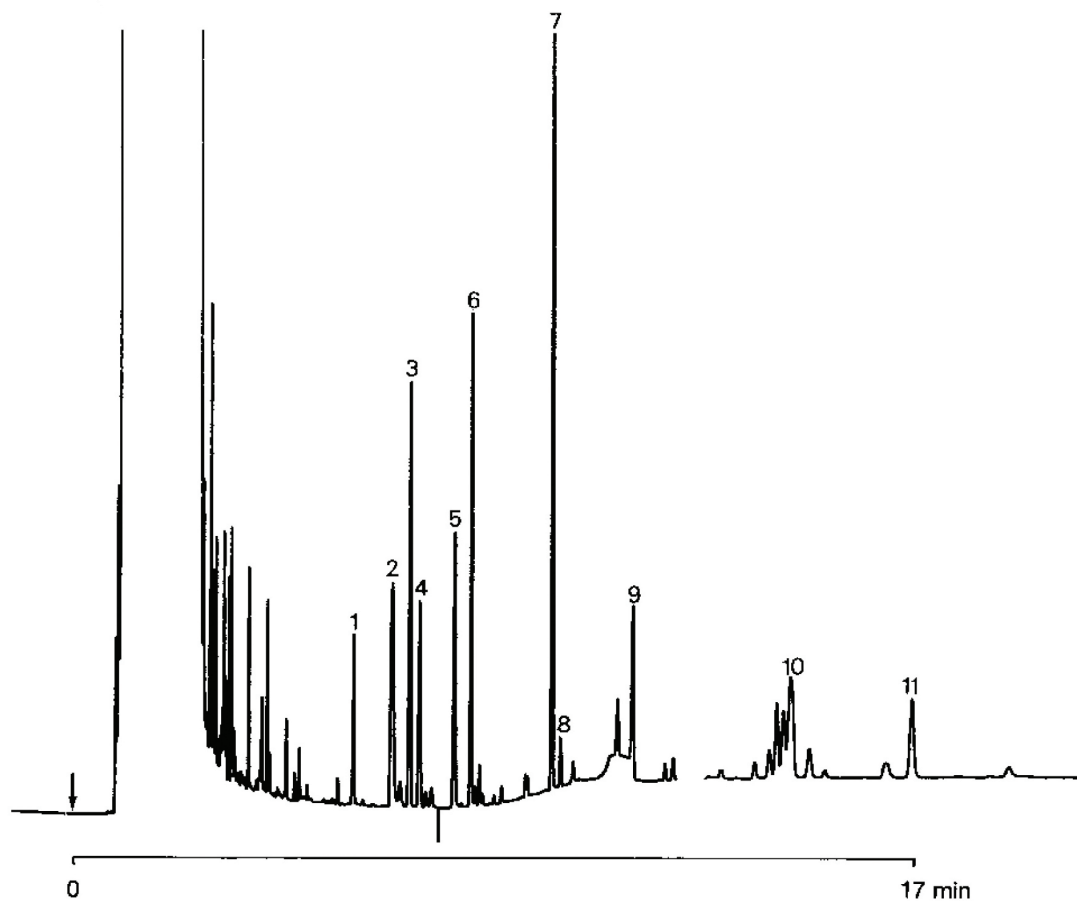
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 5, 0.32 mm x 25 m fused silica WCOT
CP-Sil 5 (0.12 μ m) (Part no. CP7470)
Temperature : 190 °C (6.5 min) $\rightarrow$ 275 °C, 10 °C/min
Carrier Gas : He, 50 kPa (0.5 bar, 7.3 psi), 35 cm/s
Injector : On-column
Detector : FID, 8×10^{-12} Afs
T = 300 °C
Sample Size : 2 μ L
Concentration range :

Peak identification

1. TCNB
2. dichloran
3. lindane
4. PCNB
5. pentachloroaniline
6. ronilane
7. antor
8. pp'-DDE
9. rovril
10. cypermethrin
11. decamethrin



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

A00123



Agilent Technologies