



Phosphorus pesticides

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 8 CB for Pesticides column separates nine organophosphate pesticides in 15 minutes.



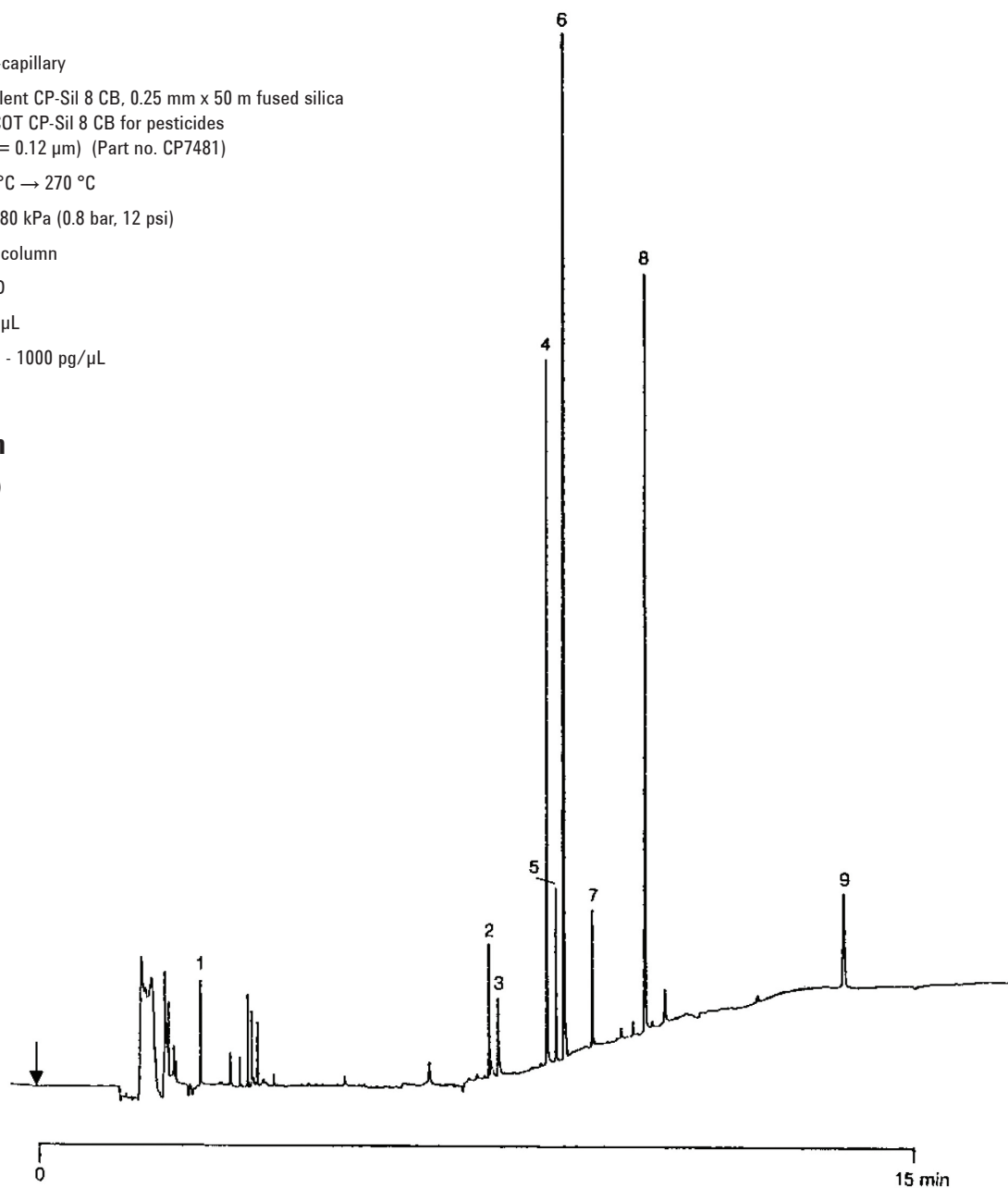
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 8 CB, 0.25 mm x 50 m fused silica
WCOT CP-Sil 8 CB for pesticides
(df = 0.12 µm) (Part no. CP7481)
Temperature : 65 °C → 270 °C
Carrier Gas : H₂, 80 kPa (0.8 bar, 12 psi)
Injector : On-column
Detector : ECD
Sample Size : 0.5 µL
Concentration range : 100 - 1000 pg/µL

Peak identification

1. dichlorvos (vapon,ddvp)
2. phorate (thimet)
3. dimethoate (cygon)
4. fenclorophos (ronnel)
5. malathion
6. ethylparathion
7. diazinon
8. ethion
9. coumaphos (coral)



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

A00358



Agilent Technologies