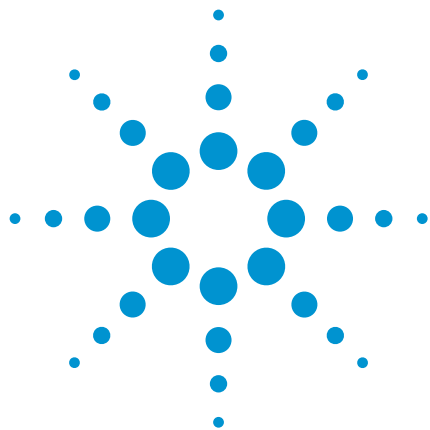




Sulfur pesticides

Analysis of miticides

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 8 CB for Pesticides column separates eight sulfur miticides in 22 minutes.



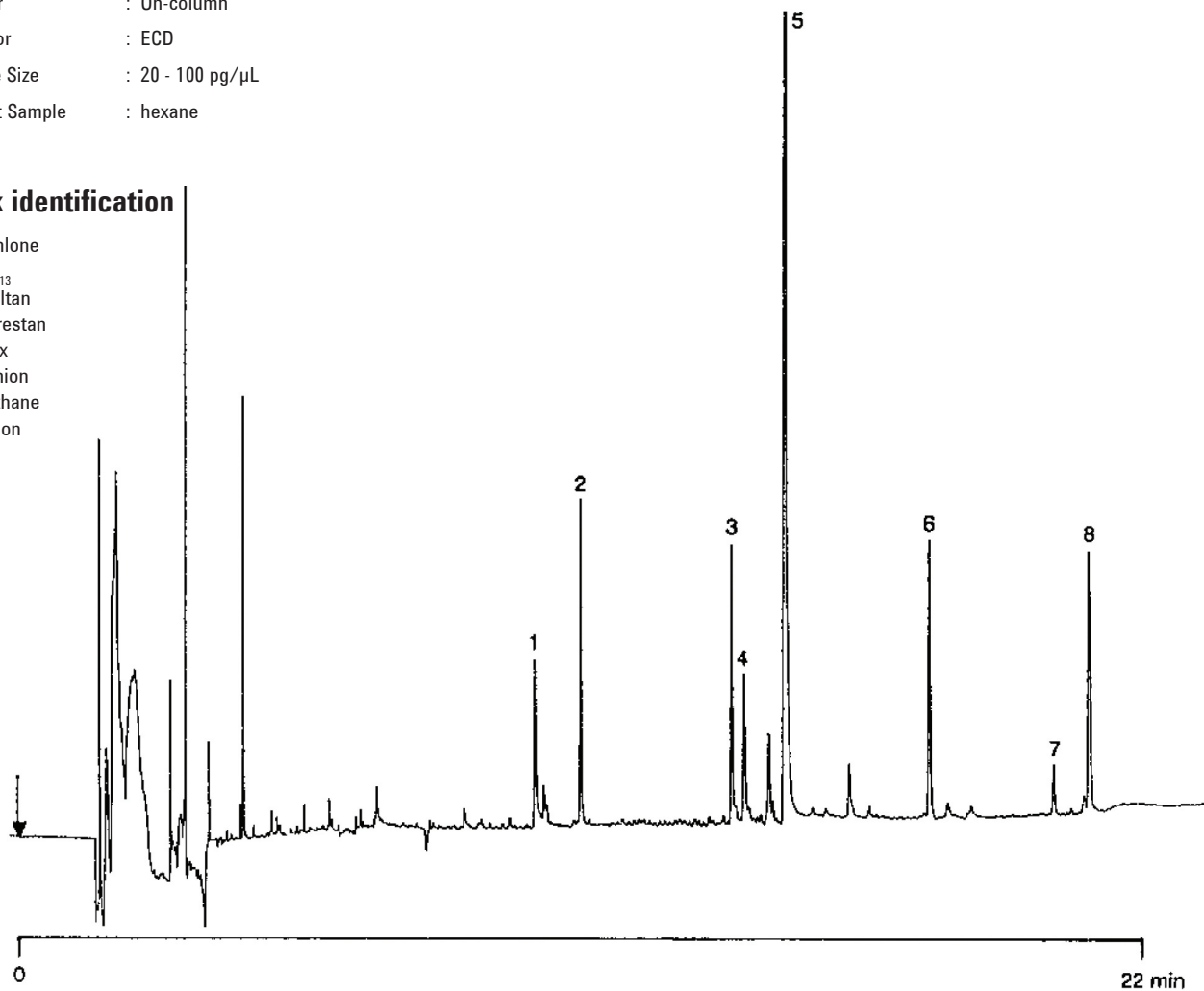
Agilent Technologies

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 → 165 °C + 5 °C/min → 270 °C
Carrier Gas : H₂, 80 kPa (0.8 bar, 12 psi)
Injector : On-column
Detector : ECD
Sample Size : 20 - 100 pg/µL
Solvent Sample : hexane

Peak identification

1. dichloro
2. VC₁₃
3. phaltan
4. morestan
5. ovex
6. trithion
7. kelthane
8. tedion



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

A00359



Agilent Technologies