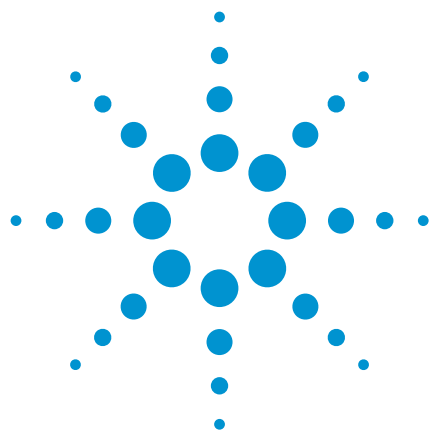




Organochlorine pesticides

Analysis of toxaphene according to EPA 8080

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 8 CB for Pesticides column separates 12 organochlorine components in toxaphene according to EPA 8080 in 58 minutes.



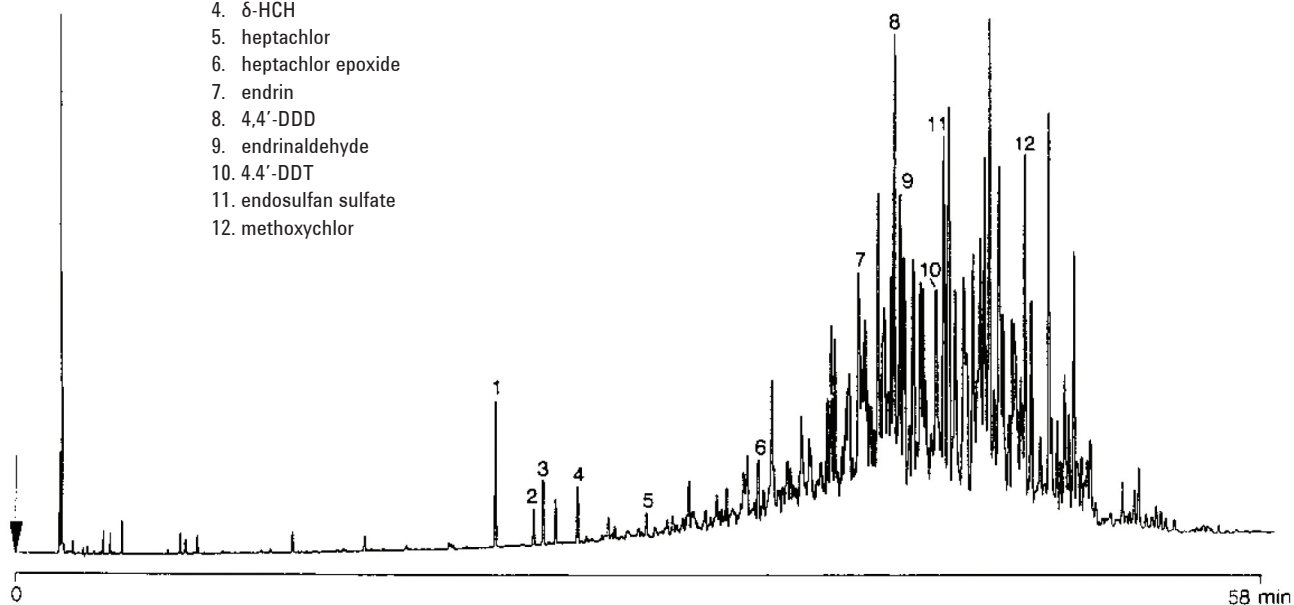
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 8 CB for Pesticides, 0.32 mm x 50 m
fused silica WCOT CP-Sil 8 CB (df = 0.12 μ m)
(Part no. CP7481)
Temperature : 100 °C (2 min) $\rightarrow$ 275 °C, 3 °C/min
Carrier Gas : H₂, 140 kPa (1.4 bar, 20 psi)
Injector : Split, 100 mL/min
T = 265 °C
Detector : ECD
T = 300 °C

Peak identification

1. α -HCH
2. β -HCH
3. γ -HCH
4. δ -HCH
5. heptachlor
6. heptachlor epoxide
7. endrin
8. 4,4'-DDD
9. endrinaldehyde
10. 4,4'-DDT
11. endosulfan sulfate
12. methoxychlor



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

A00799



Agilent Technologies