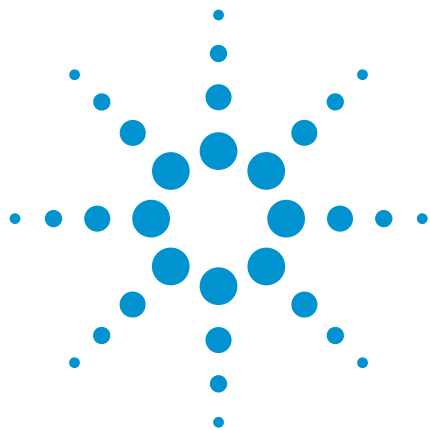




Pesticides

Analysis of α -, γ -HCH and β -, γ -PCCH

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Chirasil-DEX CB column separates four hexachlorocyclohexane pesticides in 36 minutes.



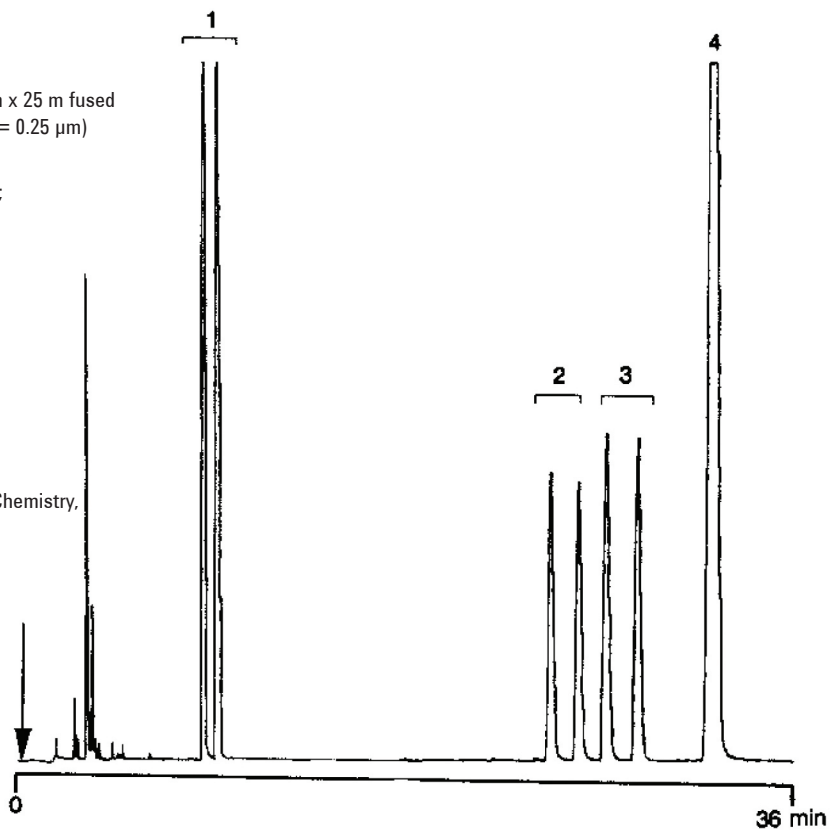
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Chirasil-DEX CB, 0.25 mm x 25 m fused silica WCOT CP-Chirasil-DEX CB (df = 0.25 µm) (Part no. CP7502)
Temperature : 150 °C (4 min) → 240 °C, 10 °C/min; 240 °C (30 min)
Carrier Gas : N₂, 80 kPa (0.8 bar, 11 psi)
Injector : Split, 100 mL/min
T = 200 °C
Detector : FID
T = 270 °C
Concentration Range : ± 25 ng/component
Courtesy : Prof. Dr. B. Luckas, Institut of Food Chemistry, University of Hohenheim, Stuttgart, Germany

Peak identification

1. γ-PCCH
2. α-HCH
3. β-PCCH
4. γ-HCH



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

A00997



Agilent Technologies