



PCBs

Analysis of Aroclor

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

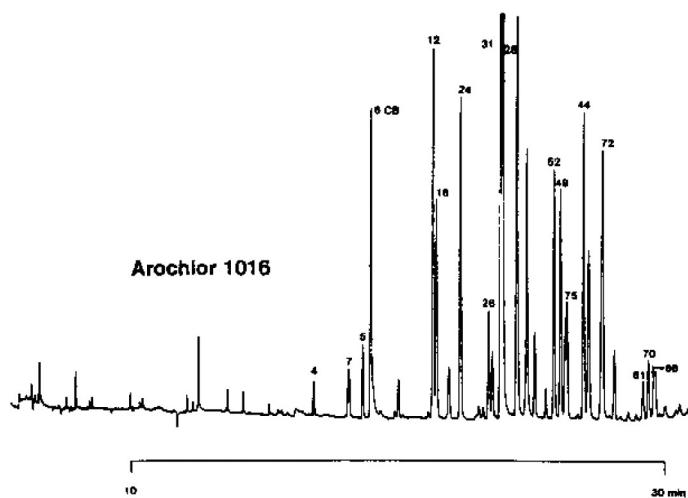
Gas chromatography with an Agilent CP-Sil 8 CB for Pesticides column separates polychlorinated biphenyls in six Aroclor PCB mixtures in 30 to 45 minutes.



Agilent Technologies

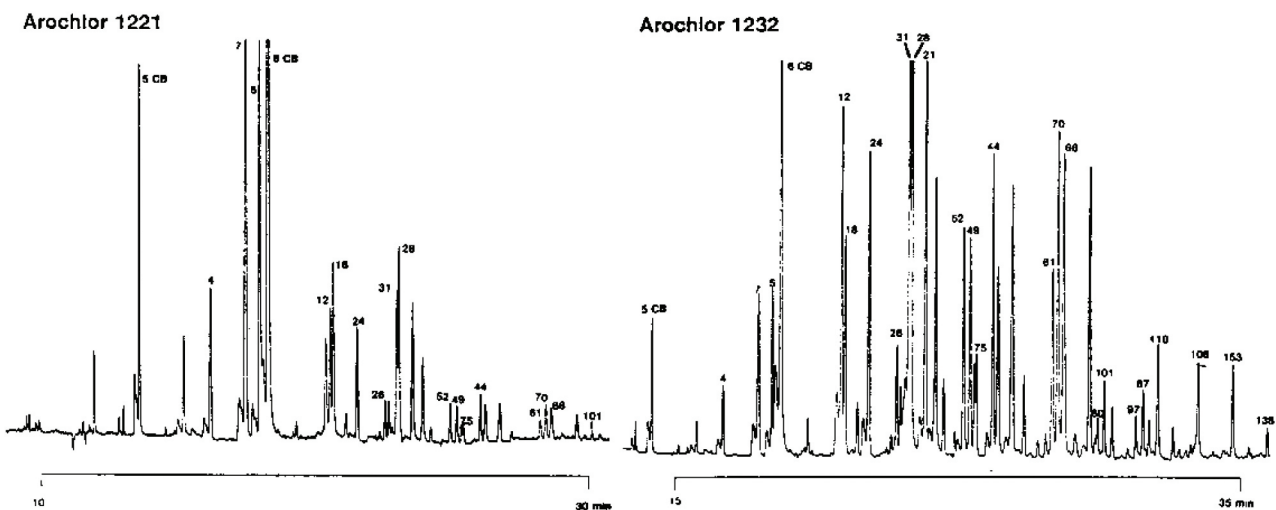
Conditions

Technique : GC-capillary
 Column : Agilent CP-Sil 8 CB for PCBs, 0.25 mm x 50 m fused silica WCOT CP-Sil 8 CB (df = 0.25 µm) (Part no. CP7482)
 Temperature : 60 °C (2 min) → 180 °C, 15 °C/min; 180 °C (6 min) → 220 °C, 4 °C/min; 220 °C (2 min) → 275 °C, 5 °C/min; 275 °C (15 min)
 Carrier Gas : H₂, 100 kPa (1 bar, 14 psi)
 Injector : Splitless, 60 s, 1.0 µL
 Detector : ECD
 Sample : 1 - 30 pg/compound/µL



Peak identification

Ballschmitter no: s

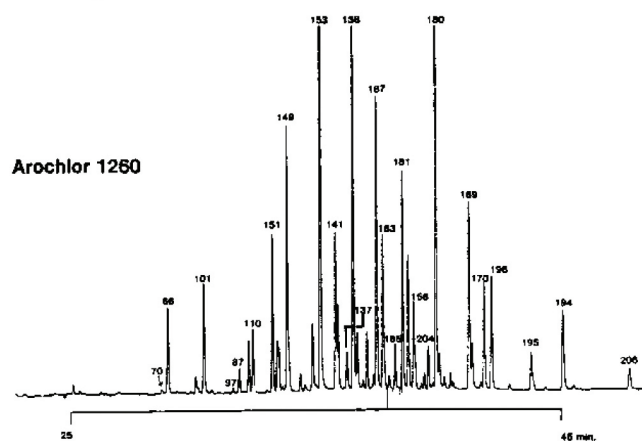
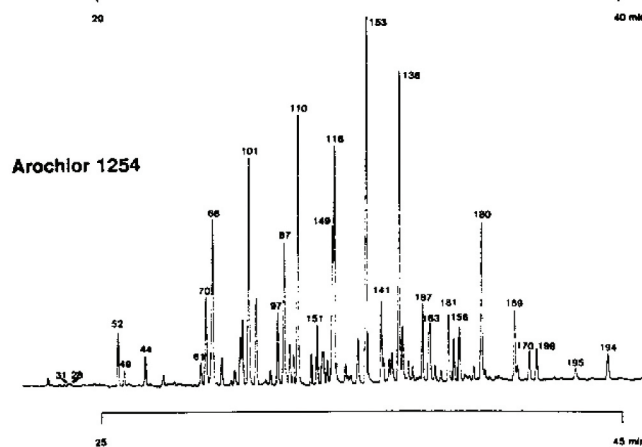
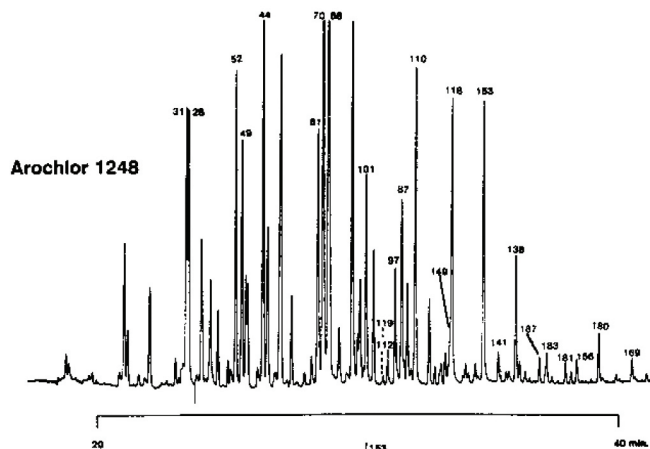


Conditions

Technique	: GC-capillary
Column	: Agilent CP-Sil 8 CB for PCBs, 0,25 mm x 50 m fused silica WCOT CP-Sil 8 CB (df = 0.25 µm) (Part no. CP7482)
Temperature	: 60 °C (2 min) → 180 °C, 15 °C/min; 180 °C (6 min) → 220 °C, 4 °C/min; 220 °C (2 min) → 275 °C, 5 °C/min; 275 °C (15 min)
Carrier Gas	: H ₂ , 100 kPa (1 bar, 14 psi)
Injector	: Splitless, 60 s, 1.0 µL
Detector	: ECD
Sample	: 1 - 30 pg/compound/µL

Peak identification

Ballschmitter no.'s



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

A00672



Agilent Technologies