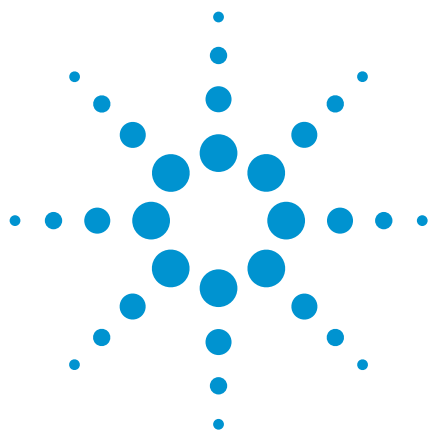




PCBs

Analysis of Aroclor 1254

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 8 CB for Pesticides column separates polychlorinated biphenyls in Aroclor 1254 in 70 minutes.



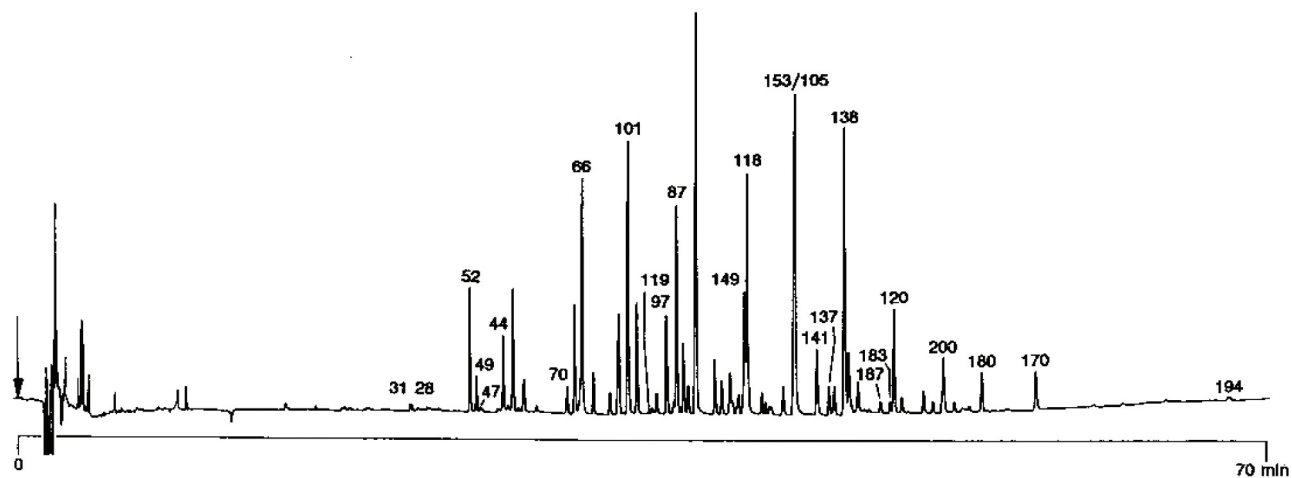
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 8 CB for pesticides, 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 (ballistic); 165 °C → 250 °C,
5° C/min
Injector : On-column
Detector : ECD
Sample Size : 1 ng/µL

Peak identification

Ballschmider 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

A00666



Agilent Technologies