



PCBs

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 5 CB column separates 30 polychlorinated biphenyls in 65 minutes.



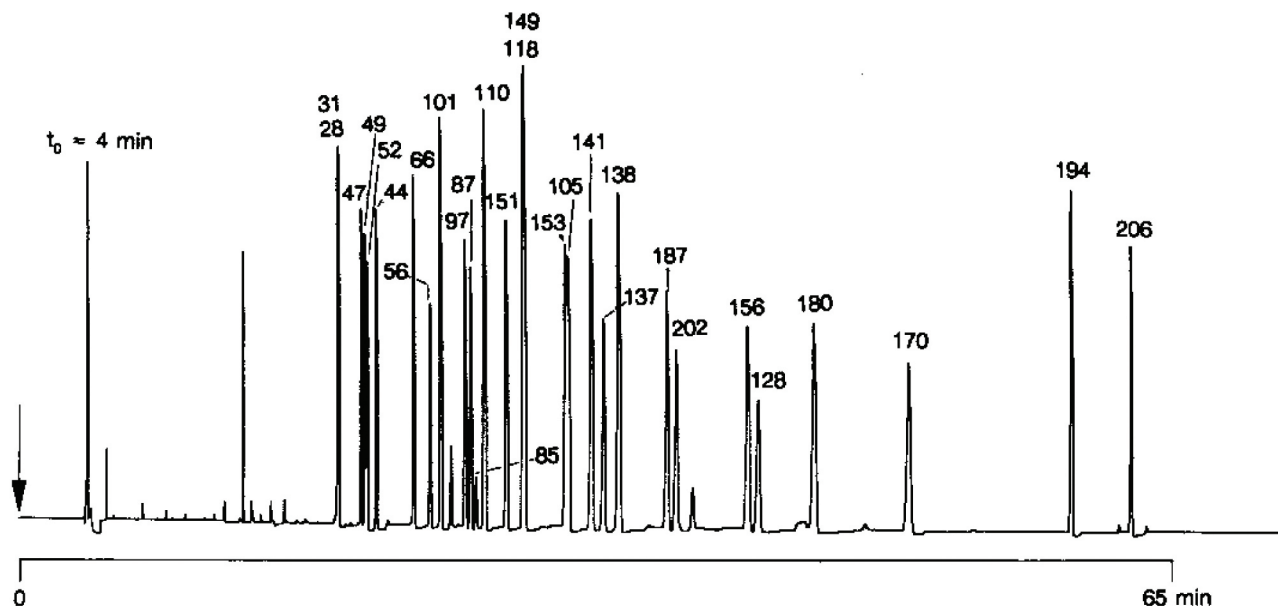
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 5 CB, 0.25 mm x 50 m fused silica
WCOT CP-Sil 5 CB (df = 0.12 µm)
(Part no. CP7740)
Temperature : 90 °C (3 min) → 215 °C, 15 °C/min;
215 °C (40 min) → 270 °C, 5 °C/min;
270 °C (60 min)
Carrier Gas : He, 200 kPa (2 bar, 29 psi)
Injector : Splitter
T = 270 °C
Detector : ECD, attn 32
T = 315 °C
Sample Size : 0.2 µL
Concentration Range : 1 ppb

Peak identification

PCB numbers according to Ballschmiter



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

A00916



Agilent Technologies