



Polychlorodibenzofurans

Analysis of halogenated hydrocarbons

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

GC/MS with an Agilent CP-Sil 88 column analyzes pyrolysis products of a technical PCB mixture, and fly ash.



Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 88, 0.22 mm x 50 m fused silica WCOT
CP-Sil 88 (0.2 µm) (Part no. CP7588)
Temperature : 100 °C (5 min) → 190 °C, 15 °C/min → 240 °C,
2 °C/min 240 °C /40 min
Carrier Gas : He
Injector : Splitter
T =260 °C
Detector : mass selective detection (MSD) in SIM-mode
Solvent Sample : tetradecane
Courtesy : K. Ballschmiter, Department of Analytical Chemistry,
University of Ulm, D-7900 Ulm/Donau, Germany

