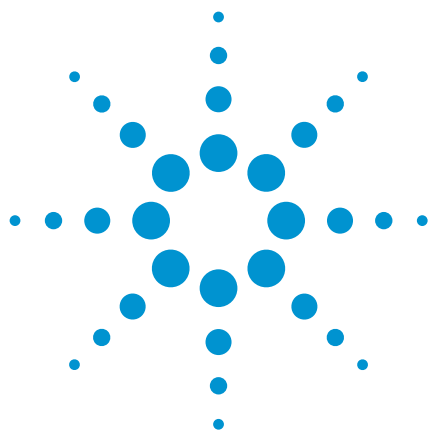




# Dioxins

## Application Note

Environmental

### Authors

Agilent Technologies, Inc.

### Introduction

GC/MS with an Agilent CP-Sil 88 for Dioxins column separates seven components in a sample of dioxin in 26 minutes.



**Agilent Technologies**

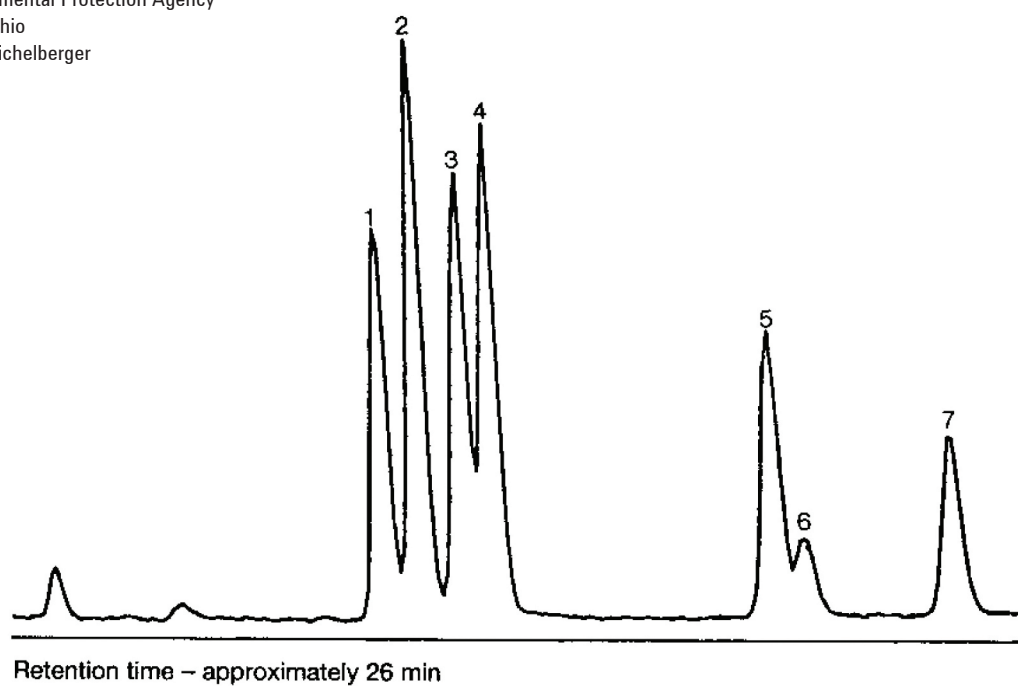
## Conditions

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 : 45 °C (3 min) → 190 °C (ballistic), → 240 °C,  
5 °C/min  
Injector : Splitless time: 1 min  
Detector : Mass Spectrometer with selected ion monitoring  
m/z = 320 and 322

Courtesy : Environment Monitoring and Support Laboratory  
U.S. Environmental Protection Agency  
Cincinnati/Ohio  
Mr James Eichelberger

## Peak identification

1. 1, 4, 7, 8 TCDD
2. 2, 3, 7, 8, TCDD
3. 1, 2, 3, 7 TCDD  
and 1, 2, 3, 8 TCDD
4. 1, 2, 3, 4, TCDD
5. 1, 2, 7, 8, TCDD
6. 1, 4, 6, 9, TCDD
7. 1, 2, 6, 7, TCDD



[www.agilent.com/chem](http://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

A00203



**Agilent Technologies**