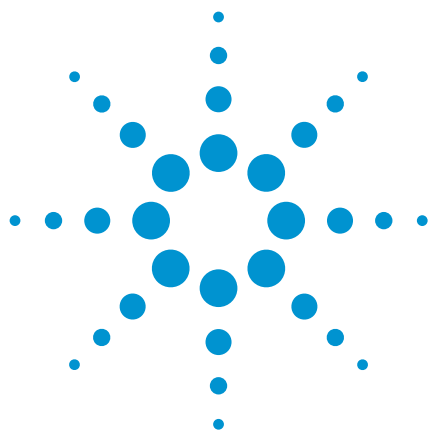




Polycyclic aromatic hydrocarbons

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

GC/MS using an Agilent CP-Sil 5 CB Low Bleed/MS column separates ten polycyclic aromatic hydrocarbons in water in 19 minutes.



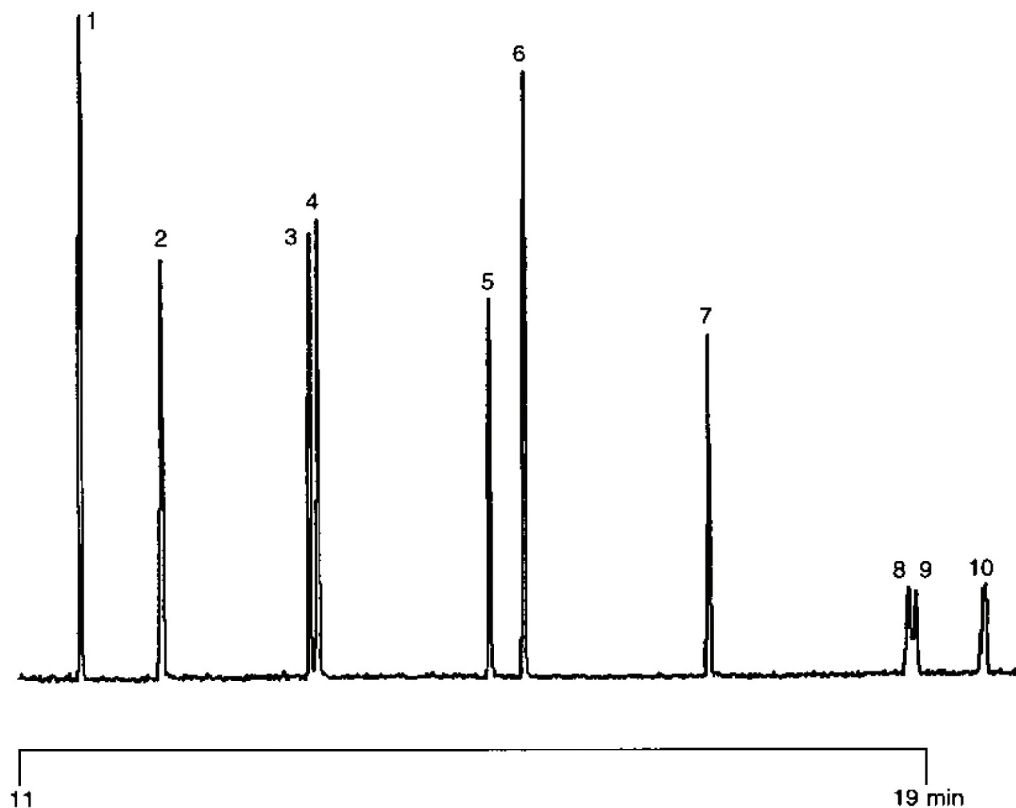
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 5 CB Low Bleed/MS, 0.25 mm x 25 m
fused silica WCOT CP-Sil 5 CB (df = 0.25 μ m)
(Part no. CP7842)
Temperature : 50 °C (5 min) $\rightarrow$ 250 °C, 20 °C/min; 250 °C (15 min)
Carrier Gas : He
Injector : Split,
T = 250 °C
Detector : MS,
T = 250 °C
Sample Size : 0.3 μ L
Concentration Range : 50 mg/L

Peak identification

1. acenaphthene
2. fluorene
3. phenanthrene
4. anthracene
5. fluoranthene
6. pyrene
7. benz[a]anthracene
8. benzo[a]pyrene
9. benz[e]acephenanthrylene
10. benzo[ghi]perylene



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

A01226



Agilent Technologies