



Polycyclic aromatic hydrocarbons

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with a custom-made Agilent CP-Sil 8 column separates 20 polycyclic aromatic hydrocarbons in 20 minutes.



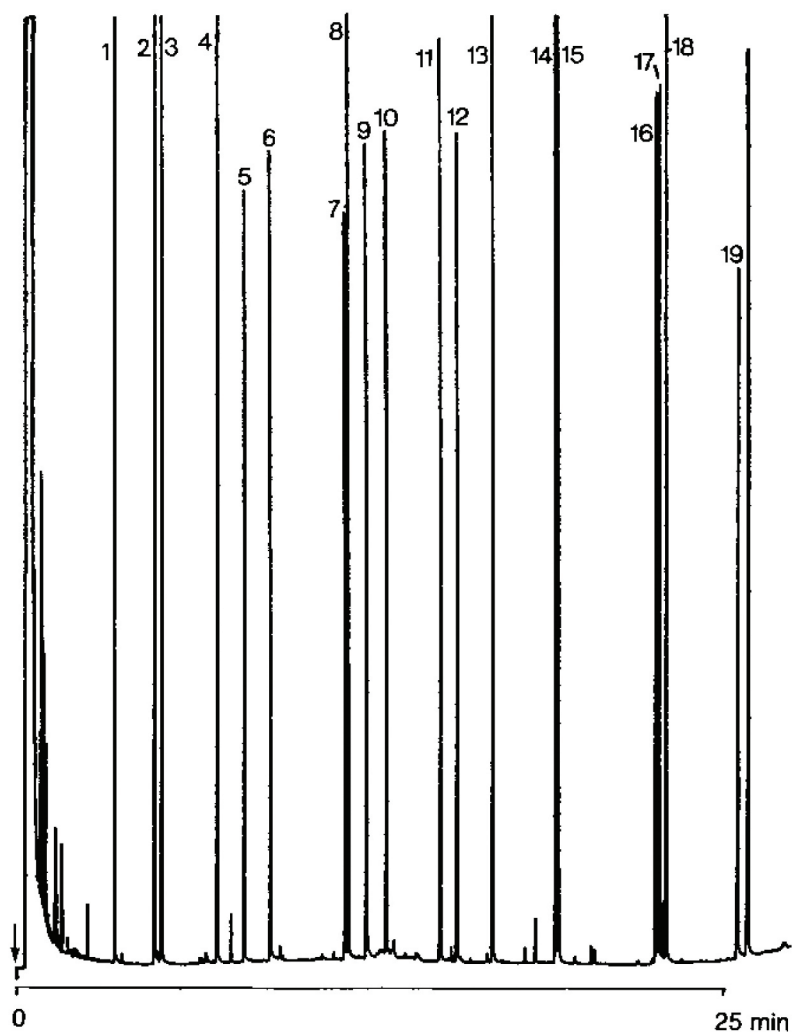
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 8, 0.32 mm x 25 m fused silica WCOT
CP-Sil 8 (0.12 μ m) (Part no. CP6030)
Temperature : 65 °C $\rightarrow$ 325 °C, 6 °C/min
Carrier Gas : H₂, 35 kPa (0.35 bar, 5.1 psi), 60 cm/s
Injector : On-column
Detector : FID, 8 x 10⁻¹² Afs
T = 325 °C
Sample Size : 0.5 μ L

Peak identification

1. naphthalene
2. 2-methylnaphthalene
3. 1-methylnaphthalene
4. acenaphthalene
5. dibenzofuran
6. fluorene
7. phenanthrene
8. anthracene
9. carbazole
10. 2-methylphenanthrene
11. fluoranthene
12. pyrene
13. 2,3-dibenzofluorene
14. benz (a) anthracene
15. chrysene
16. benz (e) pyrene
17. benz (a) pyrene
18. perylene
19. dibenzanthracene
20. benz (g,h,i) 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

A00125



Agilent Technologies