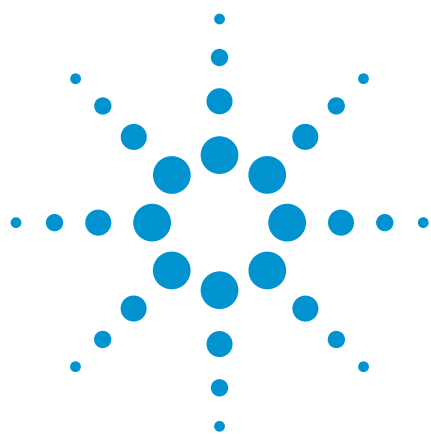




Organic water pollutants

Analysis of organic water pollutants to EPA 525

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 8 CB column separates 41 organic water pollutants according to EPA 525 in 50 minutes.



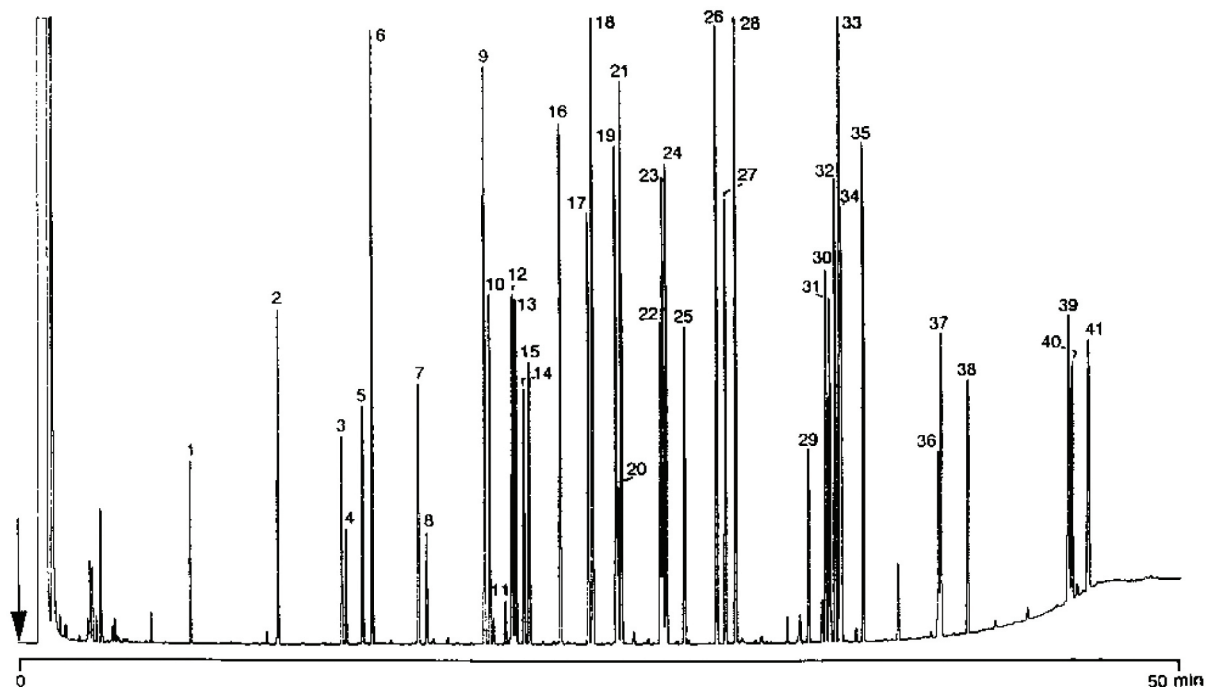
Agilent Technologies

Conditions

Technique : GC-capillary
 Column : Agilent CP-Sil 8 CB fused, 0.25 mm x 25 m silica
 WCOT CP-Sil 8 CB (df = 0.25 µm) (Part no. CP7451)
 Temperature : 80 °C (2 min) → 300 °C, 5 °C/min
 Carrier Gas : H₂, 60 kPa (0.6 bar, 8.6 psi)
 Injector : Split, 100 mL/min
 T = 275 °C
 Detector : FID
 T = 300 °C

Peak identification

1. naphthalene	11. simazine	21. aldrin	31. chrysene
2. hexachlorocyclopentadiene	12. altrazine	22. fluoranthene	32. 2,2',3,3',4,4',6-heptachlorobiphenyl
3. acenaphthylene	13. lindane	23. heptachlorepoxyde	33. methoxychlor
4. dimethylphthalate	14. phenanthrene	24. 2,2',3',4,6-pentachlorobiphenyl	34. 2,2',3,3',4,5',6,6-octachlorobiphenyl
5. acenaphthene	15. anthracene	25. pyrene	35. bis(2-ethylhexyl)phthalate
6. 2-chlorobiphenyl	16. 2,4,5-trichlorobiphenyl	26. dieldrin	36. benzo(b)fluoranthene
7. fluorene	17. heptachlor	27. 2,2',4,4',5,6'-hexachlorobiphenyl	37. benzo(k)fluoranthene
8. diethylphthalate	18. alachlor	28. endrin	38. benzo(a)pyrene
9. 2,3-dichlorobiphenyl	19. 2,2',4,4'-tetrachlorobiphenyl	29. bis(2-ethylhexyl)adipate	39. indeno(1,2,3-cd)pyrene
10. hexachlorobenzene	20. di-n-butylphthalate	30. benzo(a)anthracene	40. dibenzo(a,h)anthracene
			41. benzo(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

A00781



Agilent Technologies