



# Pollutants

## Analysis of pollutants to EPA 625

### Application Note

Environmental

#### Authors

Agilent Technologies, Inc.

#### Introduction

Gas chromatography with an Agilent CP-Sil 8 CB column separates 82 pollutants according to EPA 625 in 65 minutes.



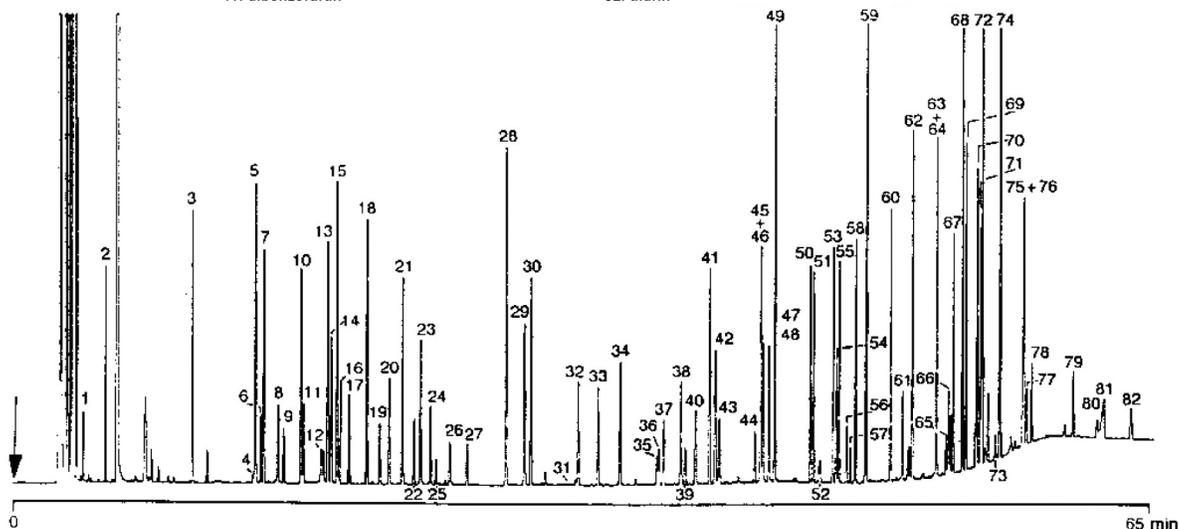
**Agilent Technologies**

## Conditions

Technique : GC-capillary  
Column : Agilent CP-Sil 8 CB, 0.25 mm x 50 m fused silica  
WCOT CP-Sil 8 CB (df = 0.25 µm) (Part no. CP7453)  
Temperature : 50 °C (2 min) → 165 °C, 3 °C/min; 165 °C → 300 °C,  
8 °C/min; 300 °C (10 min)  
Carrier Gas : H<sub>2</sub>, 80 kPa (0.8 bar, 11 psi)  
Injector : Split, 100 mL/min;  
T = 275 °C  
Detector : FID; T = 300 °C

## Peak identification

1. N-nitrosodimethylamine	21. 2,4-dimethylphenol	42. 2,4-dinitrophenol	63. heptachlorepoxyde
2. N-nitrosomethylethylamine	22. bis(2-chloroethoxy)methane	43. 2,4-dinitrotoluene	64. fluoranthene
3. N-nitrosodimethylamine	23. 2,4-dichlorophenol	44. diethylphthalate	65. benzidene
4. aniline	24. 1,2,4-trichlorobenzene	45. fluorene	66. pyrene
5. phenol	25. naphthalene	46. 4-chlorophenyl phenylether	67. endosulfan I
6. bis(2-chloroethyl)ether	26. 4-chloroaniline	47. 4-nitroaniline	68. p,p'-DDE
7. 2-chlorophenol	27. hexachlorobutadiene	48. 2-methyl-4,6-dinitrophenol	69. dieldrin
8. 1,3-dichlorobenzene	28. N-nitroso-di-n-butylamine	49. N-nitrosodiphenylamine	70. endrin
9. 1,4-dichlorobenzene	29. 4-chloro-3-methylphenol	50. 4-bromophenyl phenylether	71. endosulfan II
10. benzylalcohol	30. 2-methylnaphthalene	51. α-HCH	72. p,p'-DDD
11. 1,2-dichlorobenzene	31. hexachlorocyclopentadiene	52. hexachlorobenzene	73. butylbenzyl phthalate
12. bis(2-chloroisopropyl)ether	32. 2,4,6-trichlorophenol	53. β-HCH	74. p,p'-DDT
13. N-nitrosopyrrolidine	33. 2-chloronaphthalene	54. pentachlorophenol	75. benzo(a)anthracene
14. N-nitrosomorpholine	34. 2-nitroaniline	55. γ-HCH	76. 3,3'-dichlorobenzidine
15. N-nitroso-di-n-propylamine	35. dimethylphthalate	56. phenanthrene	77. chrysene
16. hexachloroethane	36. acenaphthylene	57. anthracene	78. bis(2-ethylhexyl)phthalate
17. nitrobenzene	37. 2,6-dinitrotoluene	58. δ-HCH	79. di-n-octyl phthalate
18. N-nitrosopiperidine	38. 3-nitroaniline	59. carbazol	80. indeno(1,2,3-cd)pyrene
19. isophorone	39. acenaphthene	60. heptachlor	81. dibenzo(a,h)anthracene
20. 2-nitrophenol	40. 4-nitrophenol	61. di-n-butyl phthalate	82. benzo(g,h,i)perylene
	41. dibenzofuran	62. aldrin	



[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

A00789



**Agilent Technologies**