

Application Report 391

Analysis of US EPA Method 8270D Semivolatiles Using SLB-5ms

US EPA Method 8270D specifies a procedure for the determination of semivolatiles in solid wastes and ground waters. The suitability of the SLB-5ms for the separation of 72 analytes (plus eight surrogates and six internal standards) specified in the method is demonstrated. The optimized chromatogram obtained on the SLB-5ms is presented.

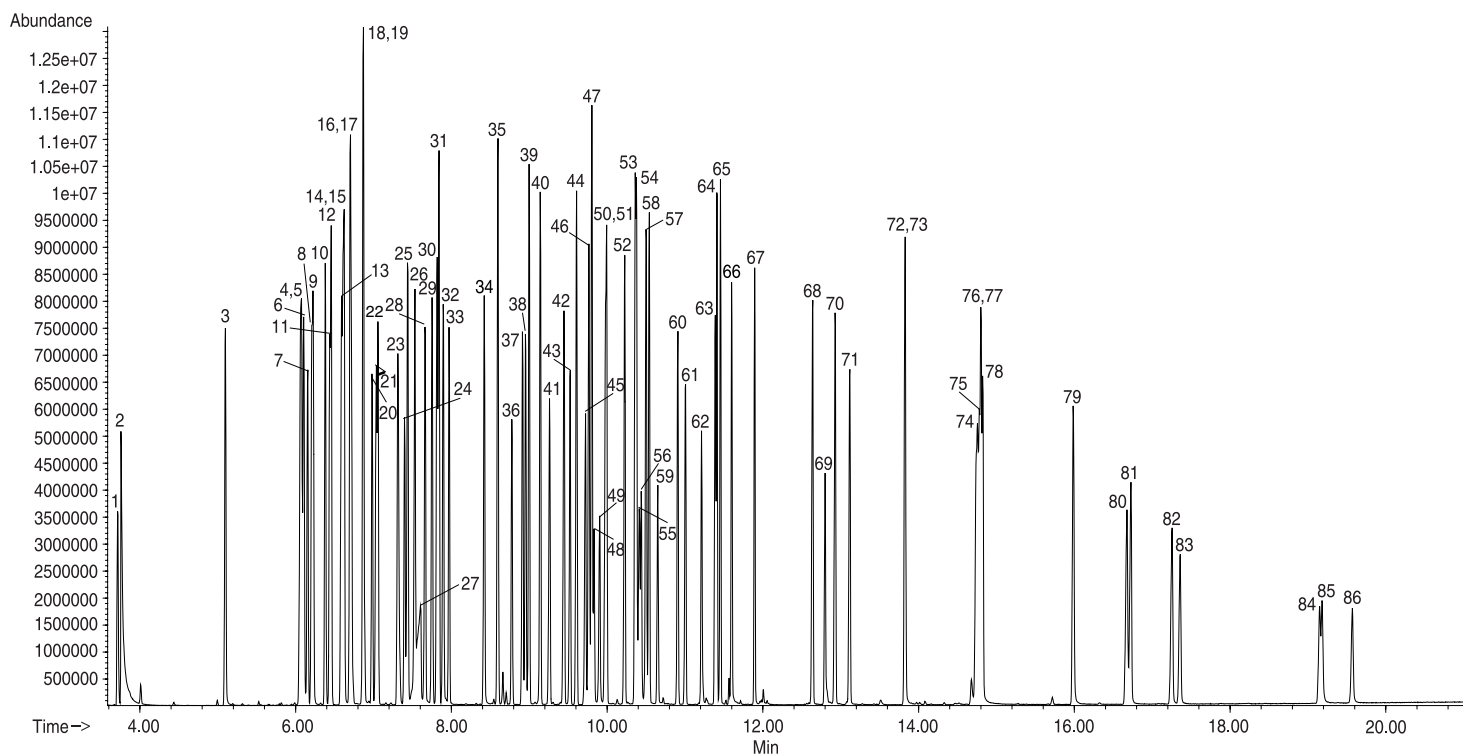
Key Words

semivolatiles, BNA, 506508, 861148, 46702-U, 48467, 861155, 46955-U, US EPA Method 8270D, RCRA, SLB-5ms, 28471-U

Author: Katherine Stenerson

Acquisition System: 6890 GC / 5973 MSD

Notebook Reference: 1569-10



G003533

Conditions

column: SLB-5ms, 30 m x 0.25 mm I.D., 0.25 μ m (28471-U)
oven: 40 $^{\circ}$ C (2 min.), 22 $^{\circ}$ C/min. to 240 $^{\circ}$ C, 10 $^{\circ}$ C/min. to 330 $^{\circ}$ C, (1 min.)
inj.: 250 $^{\circ}$ C
MSD interface: 330 $^{\circ}$ C
scan range: 40-450 m/z
carrier gas: helium, 1.0 mL/min. (11 min.), 10 mL/min² to 1.5 mL/min. (hold remainder of run)
injection: 0.5 μ L, pulsed (15 psi until 0.10 min.) splitless (0.50 min.)
liner: 2 mm I.D., straight
sample: 50 ng on-column of a 72 component semivolatile standard and 8 surrogate compounds, plus 6 internal standards (at 40 ng on-column)

27. Benzoic acid
28. 2,4-dichlorophenol
29. 1,2,4-trichlorobenzene
30. Naphthalene-d₈ (I.S.)
31. Naphthalene
32. 4-chloroaniline
33. Hexachlorobutadiene
34. 4-chloro-3-methylphenol
35. 2-methylnaphthalene
36. Hexachlorocyclopentadiene
37. 2,4,6-trichlorophenol
38. 2,4,5-trichlorophenol
39. 2-fluorobiphenyl (surr.)
40. 2-chloronaphthalene
41. 2-nitroaniline
42. Dimethyl phthalate
43. 2,6-dinitrotoluene
44. Acenaphthylene
45. 3-nitroaniline
46. Acenaphthene-d₁₀ (I.S.)
47. Acenaphthene
48. 2,4-dinitrophenol
49. 4-nitrophenol
50. Dibenzofuran
51. 2,4-dinitrotoluene
52. Diethyl phthalate
53. 4-chlorophenyl phenyl ether
54. Fluorene
55. 4-nitroaniline
56. 2-methyl-4,6-dinitrophenol
57. N-nitrosodiphenylamine
58. Azobenzene

59. 2,4,6-tribromophenol (surr.)
60. 4-bromophenyl phenyl ether
61. Hexachlorobenzene
62. Pentachlorophenol
63. Phenanthrene-d₁₀ (I.S.)
64. Phenanthrene
65. Anthracene
66. Carbazole
67. Di-n-butyl phthalate
68. Fluoranthene
69. Benzidine
70. Pyrene
71. Terphenyl-d₁₄ (surr.)
72. 3,3'-dimethylbenzidine
73. Butylbenzyl phthalate
74. 3,3'-dichlorobenzidine
75. Benzo(a)anthracene
76. Bis(2-ethylhexyl)phthalate
77. Chrysene-d₁₂ (I.S.)
78. Chrysene
79. Di-n-octyl phthalate
80. Benzo(b)fluoranthene
81. Benzo(k)fluoranthene
82. Benzo(a)pyrene
83. Perylene-d₁₂ (I.S.)
84. Indeno(1,2,3-cd)pyrene
85. Dibenzo(a,h)anthracene
86. Benzo(g,h,i)perylene

Peak IDs

1. N-nitrosodimethylamine	14. 1,2-dichlorobenzene-d ₄ (surr.)
2. Pyridine	15. 1,2-dichlorobenzene
3. 2-fluorophenol (surr.)	16. 2-methylphenol
4. Phenol-d ₆ (surr.)	17. Bis(2-chloroisopropyl)ether
5. Phenol	18. N-nitroso-di-n-propylamine
6. Aniline	19. 4-methylphenol
7. Bis(2-chloroethyl)ether	20. Hexachloroethane
8. 2-chlorophenol-d ₄ (surr.)	21. Nitrobenzene-d ₅ (surr.)
9. 2-chlorophenol	22. Nitrobenzene
10. 1,3-dichlorobenzene	23. Isophorone
11. 1,4-dichlorobenzene-d ₄ (I.S.)	24. 2-nitrophenol
12. 1,4-dichlorobenzene	25. 2,4-dimethylphenol
13. Benzyl alcohol	26. Bis(2-chloroethoxy)methane