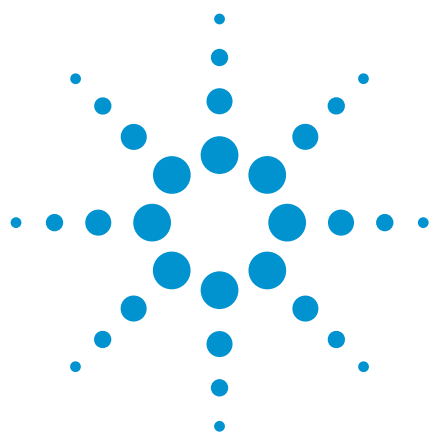




Hydrocarbons, $C_5 - C_{12}$

Naphtha and reformat analysis to ASTM D 5134

Application Note

Energy & Fuels

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil PONA CB column separates 111 components during naphtha and reformat according to ASTM D5134 in 75 minutes.

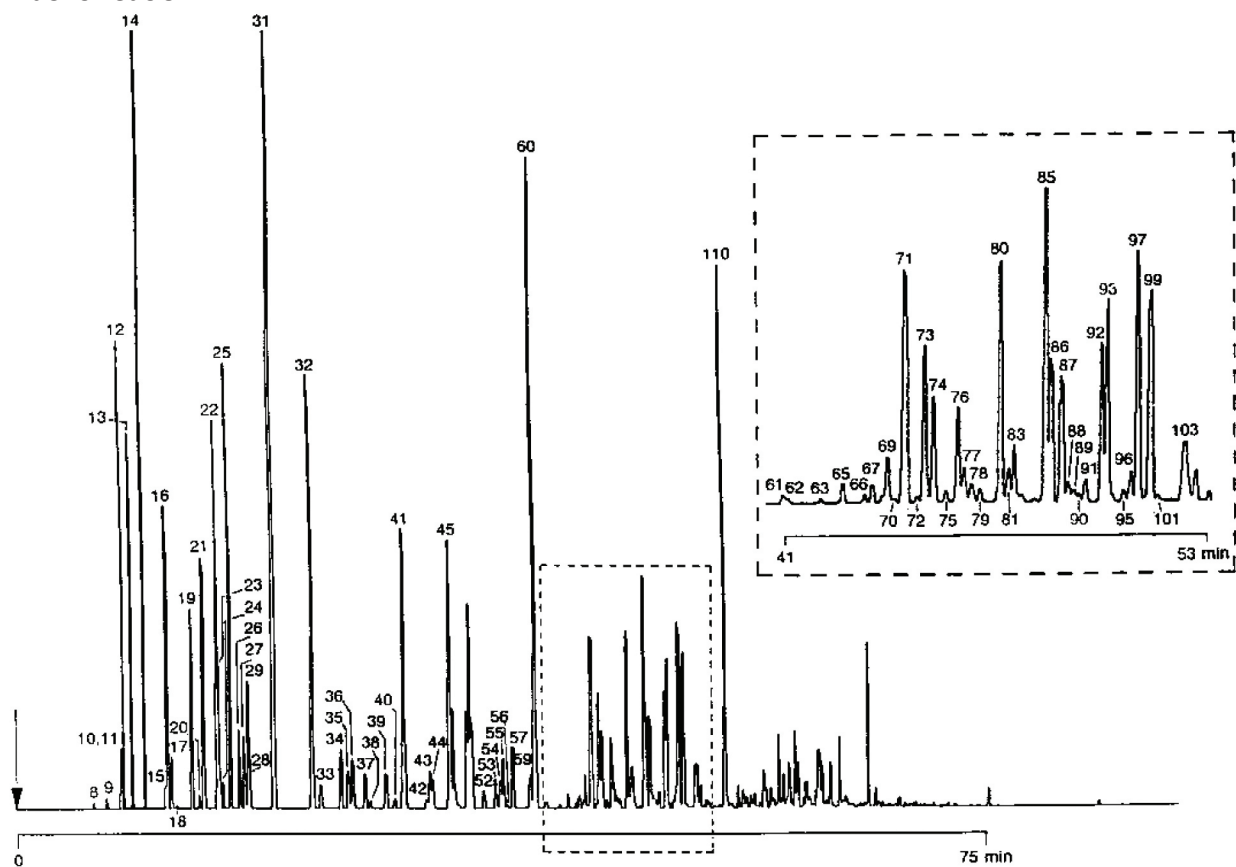


Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil PONA CB, 0.21 mm x 50 m CP-Sil
PONA CB according to ASTM D 5134
($df = 0.5 \mu m$) (Part no. CP7531)
Temperature : 35 °C (30 min) $\rightarrow$ 200 °C, 2 °C/min; 200 °C (10 min)
Carrier Gas : He, 215 kPa (2.15 bar, 30.5 psi)
Injector : Split, 200 mL/min;
T = 200 °C
Detector : FID
T = 250 °C
Sample Size : 0.1 μL
Solvent Sample : hexane

Peak identification



Conditions

Technique : GC-capillary

Column : Agilent CP-Sil PONA CB, 0.21 mm x 50 m CP-Sil
PONA CB according to ASTM D 5134
(df = 0.5 µm) (Part no. CP7531)

Temperature : 35 °C (30 min) → 200 °C, 2 °C/min; 200 °C (10 min)

Carrier Gas : He, 215 kPa (2.15 bar, 30.5 psi)

Injector : Split, 200 mL/min;
T = 200 °C

Detector : FID
T = 250 °C

Sample Size : 0.1 µL

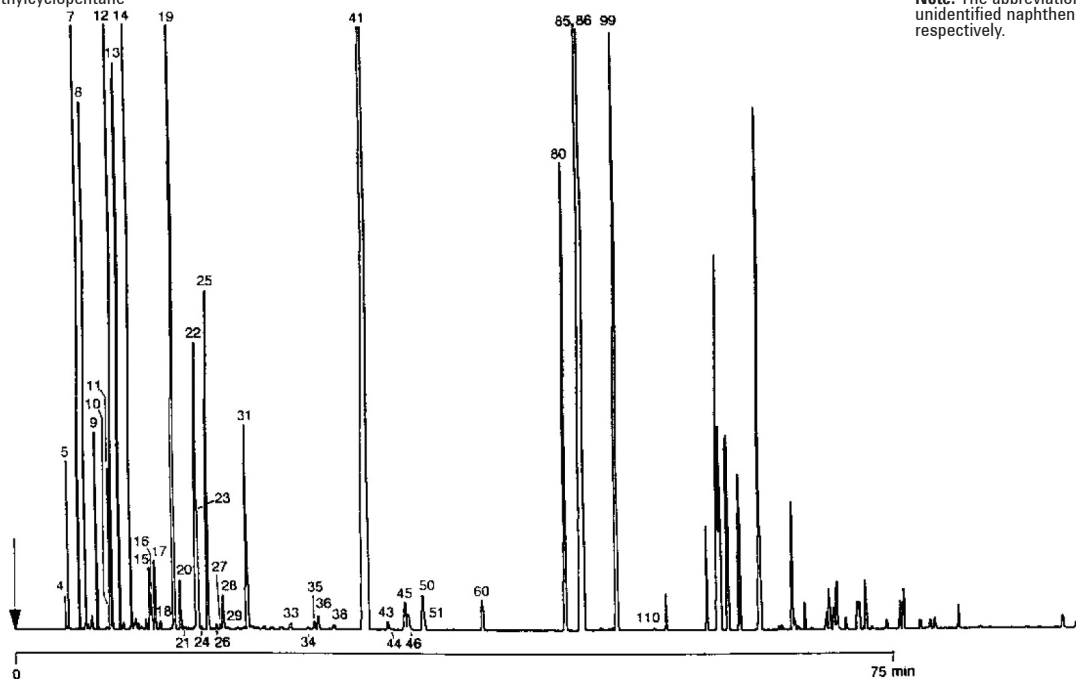
Solvent Sample : hexane

Peak identification

- | | | | |
|------------------------------------|---|--|---|
| 1. methane | 33. 1,1,3-trimethylcyclopentane + | 58. 1,cis-2,cis-3-trimethylcyclopentane | 84. unidentified naphthene + paraffin |
| 2. ethane | 2,2-dimethylhexane | 59. trans-1,3-dimethylcyclohexane + | 85. m-xylene |
| 3. propane | 34. ethylcyclopentane | cis-1,4-dimethylcyclohexane | 86. p-xylene |
| 4. isobutane | 35. 2,5-dimethylhexane + | 60. n-octane | 87. 2,3-dimethylheptane |
| 5. n-butane | 2,2,3-trimethylpentane | 61. isopropylcyclopentane + | 88. 3,4-dimethylheptane ^c + N |
| 6. 2,2-dimethylpropane | 36. 2,4-dimethylhexane | 2,4,4-trimethylhexane | 89. 3,4-dimethylheptane ^c |
| 7. isopentane | 37. 1,trans-2,cis-4-trimethylcyclopentane | 62. unidentified C ₈ -naphthene | 90. unidentified naphthene |
| 8. n-pentane | 38. 3,3-dimethylhexane | 63. unidentified C ₈ -naphthene | 91. 4-ethylheptane + N |
| 9. 2,2-dimethylbutane | 39. 1,trans-2,cis-3-trimethylcyclopentane | 64. unidentified C ₈ -naphthene | 92. 4-methyloctane |
| 10. cyclopentane | 40. 2,3,4-trimethylpentane | 65. cis-1,2-ethylmethylcyclopentane + | 93. 2-methyloctane |
| 11. 2,3-dimethylbutane | 41. toluene + 2,3,3-trimethylpentane | 2,3,5-trimethylhexane | 94. unidentified naphthene |
| 12. cyclopentane | 42. 1,1,2-trimethylcyclopentane | 66. 2,2-dimethylheptane | 95. unidentified naphthene |
| 13. 3-methylpentane | 43. 2,3-dimethylhexane | 67. cis-1,2-dimethylcyclohexane | 96. 3-ethylheptane + N |
| 14. n-hexane | 44. 2-methyl-3-ethylpentane | 68. 2,2,3-trimethylhexane + 9N | 97. 3-methyloctane |
| 15. 2,2-dimethylpentane | 45. 2-methylpentane | 69. 2,4-dimethylheptane | 98. unidentified naphthene |
| 16. methylcyclopentane | 46. 4-methylheptane + | 70. 4,4-dimethylheptane + 9N | 99. o-xylene + 1,1,2-trimethylcyclohexane |
| 17. 2,4-dimethylpentane | 3-methyl-3-ethylpentane | 71. ethylcyclohexane + | 100. unidentified naphthene + |
| 18. 2,2,3-trimethylbutane | 47. 3,4-dimethylhexane | n-propylcyclopentane | 2,4,6-trimethylheptane |
| 19. benzene | 48. 1,cis-2,trans-3-trimethylcyclopentane | 72. 2-methyl-4-ethylhexane | 101. unidentified naphthene |
| 20. 3,3-dimethylpentane | 49. cis-1,3-dimethylcyclohexane | 73. 2,6-dimethylheptane + 9N | 102. unidentified paraffin |
| 21. cyclohexane | 50. 3-methylheptane + | 74. 1,1,3-trimethylcyclohexane | 103. unidentified naphthene |
| 22. 2-methylhexane | 1,cis-2,trans-3-trimethylcyclopentane | 75. unidentified C ₈ -naphthene | 104. unidentified naphthene |
| 23. 2,3-dimethylpentane | 51. 3-ethylhexane + | 76. 2,5-dimethylheptane + 9N | 105. unidentified naphthene + paraffin |
| 24. 1,1-dimethylcyclopentane | trans-1,4-dimethylcyclohexane | 77. 3,5-dimethylheptane + N | 106. unidentified naphthene |
| 25. 3-methylhexane | 52. 1,1-dimethylcyclohexane | 78. unidentified C ₈ -naphthene | 107. unidentified naphthene |
| 26. cis-1,3-dimethylcyclopentane | 53. 2,2,5-trimethylhexane + | 79. unidentified C ₈ -naphthene | 108. unidentified naphthene |
| 27. trans-1,3-dimethylcyclopentane | trans-1,3-ethylmethylcyclopentane | 80. ethyl benzene | 109. unidentified naphthene |
| 28. 3-ethylpentane | 54. cis-1,3-ethylmethylcyclopentane | 81. unidentified C ₈ -naphthene | 110. n-nonane |
| 29. trans-1,2-dimethylcyclopentane | 55. trans-1,2-ethylmethylcyclopentane | 82. unidentified naphthene + | 111. unidentified naphthene |
| 30. 2,2,4-trimethylpentane | 56. 2,2,4-trimethylhexane + | 2,3,4-trimethylhexane | |
| 31. n-heptane | 1,1-ethylmethylcyclopentane | 83. unidentified naphthene | |
| 32. methylcyclohexane + | 57. trans-1,2-dimethylcyclohexane | | |
| cis-1,2-dimethylcyclopentane | | | |

^c - stereoisomers

Note: The abbreviations N and P refer to unidentified naphthenes and paraffins respectively.



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

A00752



Agilent Technologies