



Halogenated hydrocarbons, C₁ – C₆

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

GC/MS using an Agilent CP-Select 624 CB column separates 37 C₁ to C₆ halogenated hydrocarbons in 45 minutes.



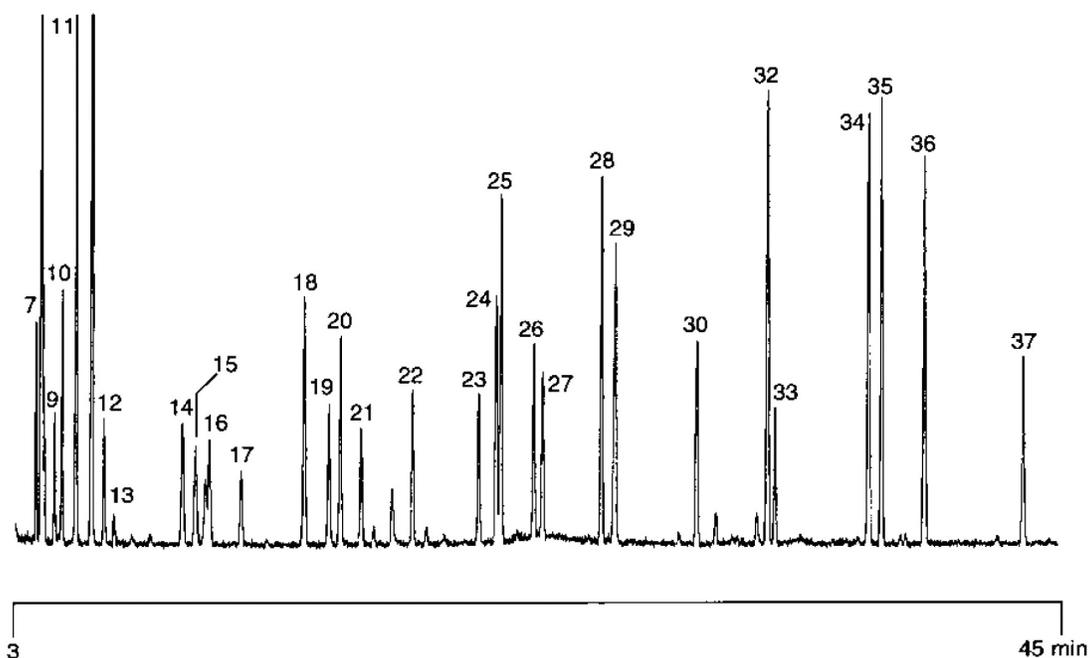
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Select 624 CB, 0.25 mm x 30 m fused silica WCOT CP-Select 624 CB (df = 1.4 µm) (Part no. CP7412)
Temperature : 35 °C (10 min) → 175 °C, 3 °C/min
Carrier Gas : He, 60 kPa (0.6 bar, 8.6 psi)
Injector : Split, 100 mL/min
T = 275 °C
Detector : MSD
T = 280 °C
Sample Size : 0.5 µL
Concentration Range : 0.1 mg/mL per compound
Solvent Sample : methanol

Peak identification

- | | |
|--------------------------------|---------------------------------|
| 1. dichlorodifluoromethane | 20. dibromomethane |
| 2. chloromethane | 21. bromodichloromethane |
| 3. vinylchloride | 22. cis-1,3-dichloropropene |
| 4. bromomethane | 23. trans-1,3-dichloropropene |
| 5. chloroethane | 24. 1,1,2-trichloroethane |
| 6. trichlorofluoromethane | 25. tetrachloroethylene |
| 7. 1,1-dichloroethylene | 26. dibromochloromethane |
| 8. methyl iodide | 27. 1,2-dibromoethane |
| 9. allyl chloride | 28. chlorobenzene |
| 10. dichloromethane | 29. 1,1,1,2-tetrachloroethane |
| 11. trans-1,2-dichloroethylene | 30. bromoform |
| 12. 1,1-dichloroethane | 31. 1,2,3-trichloropropane |
| 13. bromochloromethane | 32. 1,1,2,2-tetrachloroethane |
| 14. trichloromethane | 33. trans-1,4-dichloro-2-butene |
| 15. 1,1,1-trichloroethane | 34. 1,3-dichlorobenzene |
| 16. tetrachloromethane | 35. 1,4-dichlorobenzene |
| 17. 1,2-dichloroethane | 36. 1,2-dichlorobenzene |
| 18. trichloroethylene | 37. 1,2-dibromo-3-chloropropane |
| 19. 1,2-dichloropropane | |



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

A01232



Agilent Technologies