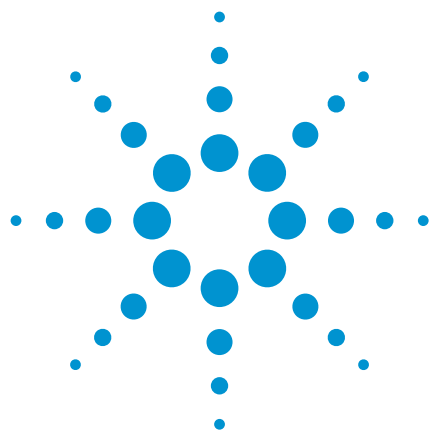




Solvents

Analysis of solvents to EPA 524.2 using GC/MS (split injector)

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Fast GC/MS with an Agilent CP-Select 624 CB column separates 60 solvents in 16 minutes according to EPA 524.2.



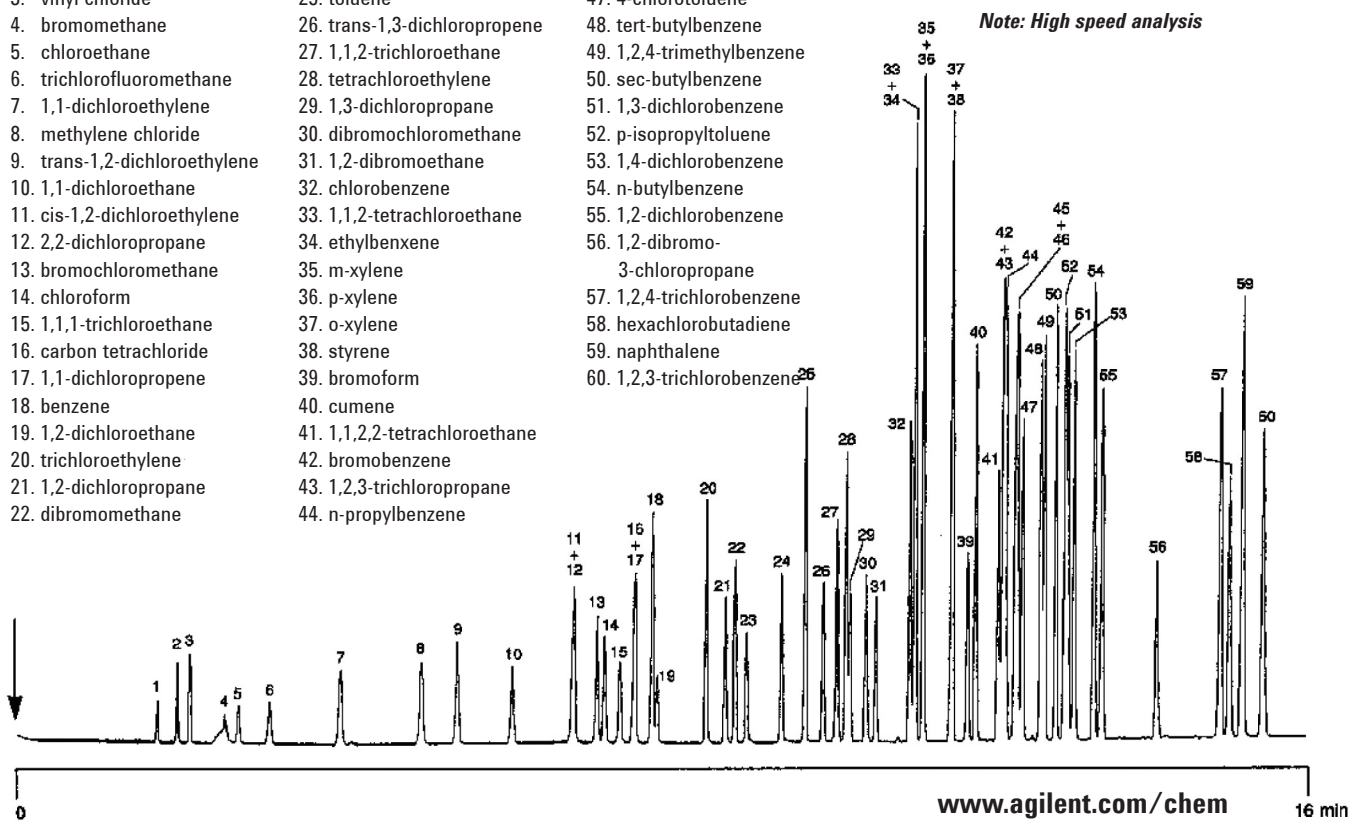
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Select 624 CB, 0.25 mm x 30 m fused silica WCOT CP-Select 624 (df = 1.4 µm) (Part no. CP7412)
Temperature : 35 °C (5 min) → 220 °C, 20 °C/min
Carrier Gas : He, 100 kPa (1.0 bar, 14.3 psi)
Injector : Splitter
Detector : MSD
Sample : 1000 ppm in methanol

Peak identification

- | | | |
|-------------------------------|-------------------------------|----------------------------|
| 1. dichlorodifluoromethane | 23. bromodichloromethane | 45. 2-chlorotoluene |
| 2. chloromethane | 24. cis-1,3-dichloropropene | 46. 1,3,5-trimethylbenzene |
| 3. vinyl chloride | 25. toluene | 47. 4-chlorotoluene |
| 4. bromomethane | 26. trans-1,3-dichloropropene | 48. tert-butylbenzene |
| 5. chloroethane | 27. 1,1,2-trichloroethane | 49. 1,2,4-trimethylbenzene |
| 6. trichlorofluoromethane | 28. tetrachloroethylene | 50. sec-butylbenzene |
| 7. 1,1-dichloroethylene | 29. 1,3-dichloropropane | 51. 1,3-dichlorobenzene |
| 8. methylene chloride | 30. dibromochloromethane | 52. p-isopropyltoluene |
| 9. trans-1,2-dichloroethylene | 31. 1,2-dibromoethane | 53. 1,4-dichlorobenzene |
| 10. 1,1-dichloroethane | 32. chlorobenzene | 54. n-butylbenzene |
| 11. cis-1,2-dichloroethylene | 33. 1,1,2-tetrachloroethane | 55. 1,2-dichlorobenzene |
| 12. 2,2-dichloropropane | 34. ethylbenzene | 56. 1,2-dibromo- |
| 13. bromochloromethane | 35. m-xylene | 3-chloropropane |
| 14. chloroform | 36. p-xylene | 57. 1,2,4-trichlorobenzene |
| 15. 1,1,1-trichloroethane | 37. o-xylene | 58. hexachlorobutadiene |
| 16. carbon tetrachloride | 38. styrene | 59. naphthalene |
| 17. 1,1-dichloropropane | 39. bromoform | 60. 1,2,3-trichlorobenzene |
| 18. benzene | 40. cumene | |
| 19. 1,2-dichloroethane | 41. 1,1,2,2-tetrachloroethane | |
| 20. trichloroethylene | 42. bromobenzene | |
| 21. 1,2-dichloropropane | 43. 1,2,3-trichloropropane | |
| 22. dibromomethane | 44. n-propylbenzene | |



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

A01255



Agilent Technologies