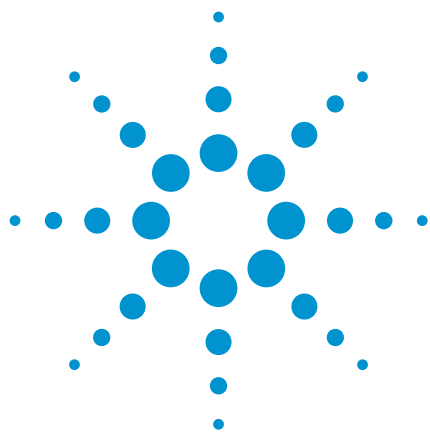




Hydrocarbons C₁ – C₉

Application Note

Environmental

Authors

Rolf Nordlinder, Olle Ramnäs and Lars-Erik Amand,
Chalmers University of Technology,
Göteborg, Sweden

Introduction

Gas chromatography using an Agilent CP-Al₂O₃/KCl column separates 35 C₁ to C₉ hydrocarbons in air in 60 minutes.



Agilent Technologies

Conditions

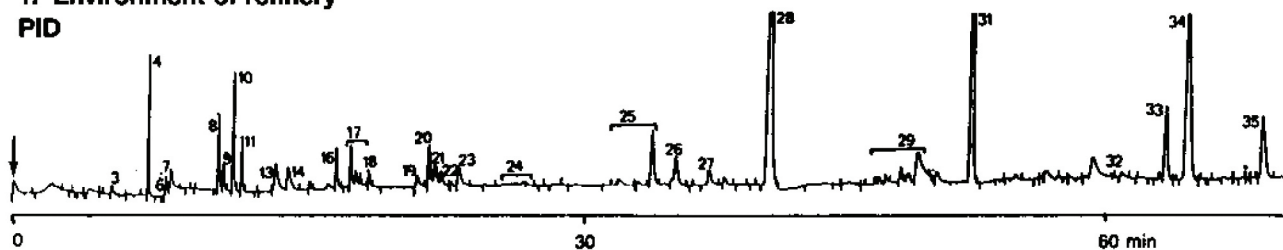
Technique : GC-capillary
 Column : Agilent CP-Al₂O₃/KCl, 0.32 mm x 50 m fused silica
 Al₂O₃/KCl PLÔT (Part no. CP7515)
 Temperature : 0 °C → 135 °C, 10 °C/min
 135 °C → 205 °C, 2 °C/min
 Carrier Gas : He, 130 kPa (1.3 bar, 19 psi)
 Injector : Air sample
 (hydrocarbons sample from 1-2 L air during 20 min)
 Detector : FID 1 x 10⁻¹² Afs
 (PID 8 x 10⁻¹² Afs (10, 2 eV lamp)
 Chromatogram 1 only)

Peak identification

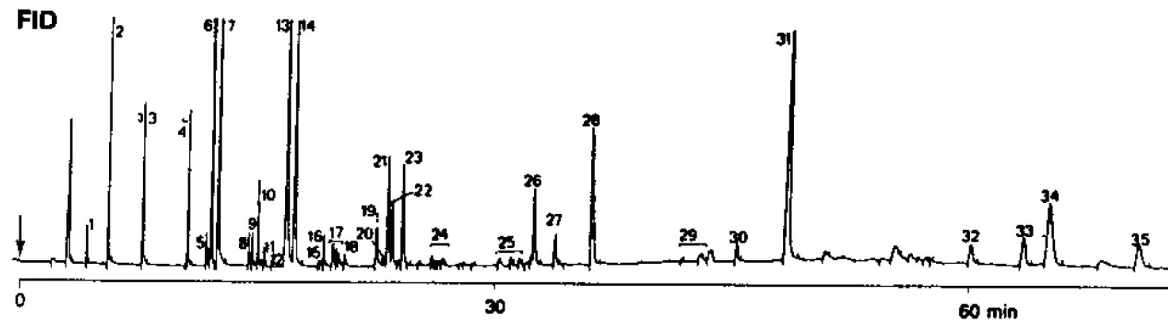
- | | |
|------------------------|---------------------------------------|
| 1. ethane | 19. methylcyclopentane |
| 2. ethene | 20. cyclohexane |
| 3. propane | 21. 2-methylpentane |
| 4. propene | 22. 3-methylpentane |
| 5. ethyne | 23. n-hexane |
| 6. methylpropane | 24. unsaturated C ₆ |
| 7. n-butane | 25. saturated branched C ₇ |
| 8. trans-2-butene | 26. 2-methylhexane |
| 9. 1-butene | 27. 3-methylhexane |
| 10. 2-methyl-1-propene | 28. n-heptane |
| 11. cis-2-butene | 29. benzene |
| 12. unknown | 30. saturated branched C ₈ |
| 13. methylbutane | 31. n-octane |
| 14. n-pentane | 32. toluene |
| 15. trans-2-pentene | 33. n-nonane |
| 16. 1-pentene | 34. ethylbenzene |
| 17. methylbutenes | 35. m-,p-xylene |
| 18. cis-2-pentene | 36. o-xylene |

1. Environment of refinery

PID



FID



Analysis of C₁-C₉ hydrocarbons in environmental air

Sampling and analysis of mixtures of light hydrocarbons (C₁-C₉) in environmental air has been a problem for several years. Gas chromatographic systems for the separation of the hydrocarbons working with two or three columns have been used. The new Al₂O₃/KCl PLOT column has now made it possible to separate complex mixtures of light hydrocarbons on one column. The column has been tested in a system for the analysis of hydrocarbons in environmental air. The system is based on a Carlo Erba 2900 gas chromatograph and has a thermal desorption oven followed by a liquid nitrogen cooling trap. The effluent from the column is split to three different detectors, one FID (Carlo Erba FID 20), one ECD (Carlo Erba AT-25) and one PID (HNU P1-52). Glass tubes filled with different layers of adsorbents and a personal monitoring pump were used for the sampling of hydrocarbons in air. A more detailed description will be published elsewhere.

We have used the column daily for more than 6 months and the number of analyses is over 1000. The separation of the hydrocarbons C₁-C₅ was excellent. Separation of hydrocarbons up to C₉ was possible with temperature programmed analysis. The column has been run both at low (< 20 °C) and high (205 °C) temperatures. Conditioning of the column at 200 °C over night has been used frequently. Water in the air samples has been the only problem.

In our case it was necessary to have a moisture filter between the desorption oven and the cooling trap. We have only used the column for air samples and no liquid samples have been injected.

The analytical system described has mainly been developed for the estimation of fugitive hydrocarbon emissions from chemical plants. The ECD is used for detection of SF₆ when tracer technique is applied. The system is also suitable for the analysis of halogenated hydrocarbons in environmental air.

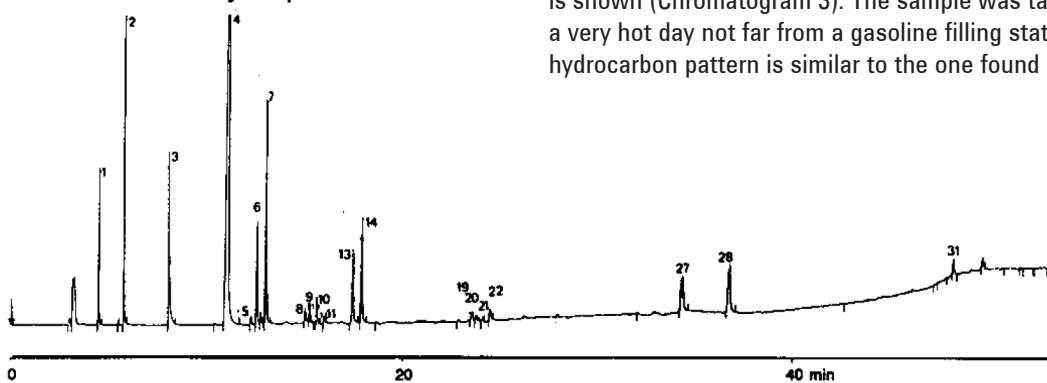
The first application is shown in Chromatogram 1. The sample was taken about 200 m downwind from a petrochemical refinery. The chromatogram shows a complex mixture of hydrocarbons where saturated aliphatics up to C₆ dominate. Aromatic hydrocarbons from the reforming unit of the refinery can also be seen.

The concentration of e.g. benzene was about 10 µg/m³. In the second application a hydrocarbon pattern different from the first sample is found (Chromatogram 2). The sample was taken downwind from an ethylene plant.

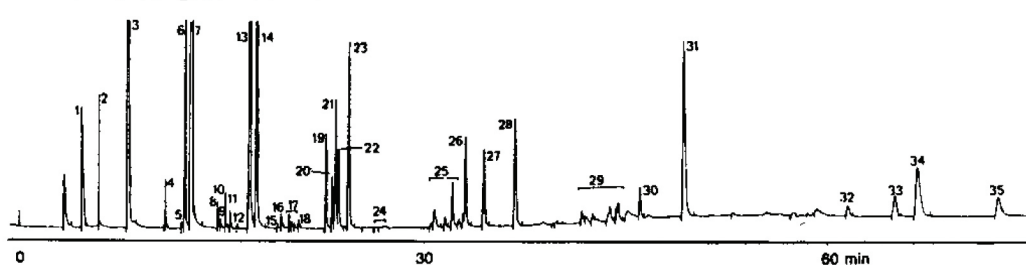
Unsaturated aliphatic hydrocarbons dominate here. The high concentration of propene probably indicates a leakage in the plant area. Aromatic hydrocarbons, mainly from the steamed cracked naphtha unit of the plant, are also determined.

Finally, an air sample from a street in the city of Göteborg is shown (Chromatogram 3). The sample was taken on a very hot day not far from a gasoline filling station. The hydrocarbon pattern is similar to the one found in gasoline.

2. Environment of ethylene plant



3. Environment of gasoline station



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

A00056



Agilent Technologies