

Aromatic hydrocarbons, $C_{16} - C_{20}$

Analysis of alkylbenzenes

Application Note

Food Testing & Agriculture

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 8 CB column separates 19 alkylbenzenes in 18 minutes.



Agilent Technologies

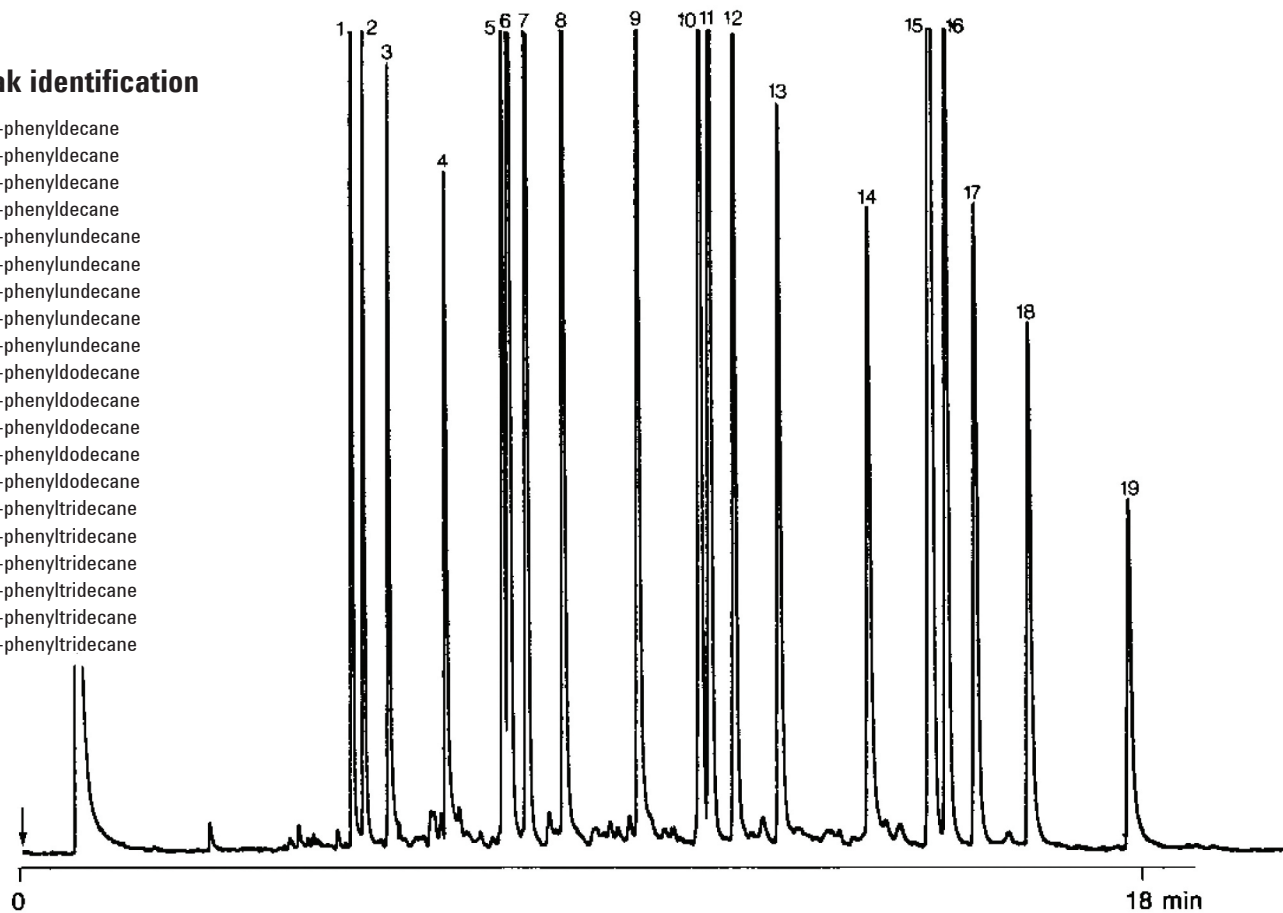
Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 8 CB, 0.22 mm x 25 m fused silica
WCOT CP-Sil 8 CB, (0.12 µm) (Part no. CP7711)
Temperature : 120 °C → 150 °C, 2 °C/min
Carrier Gas : H₂, 130 kPa (1.3 bar, 19 psi)
Injector : Splitter, 1 mL/min
T = 275 °C
Detector : FID
T = 275 °C

Courtesy : Dr. C. Mariani, Stazione sperimentale degli oli e dei grassi, Via Colombo 79, Milano, Italy

Peak identification

1. 5-phenyldecane
2. 4-phenyldecane
3. 3-phenyldecane
4. 2-phenyldecane
5. 6-phenylundecane
6. 5-phenylundecane
7. 4-phenylundecane
8. 3-phenylundecane
9. 2-phenylundecane
10. 6-phenyldodecane
11. 5-phenyldodecane
12. 4-phenyldodecane
13. 3-phenyldodecane
14. 2-phenyldodecane
15. 7-phenyltridecane
16. 6-phenyltridecane
17. 5-phenyltridecane
18. 4-phenyltridecane
19. 3-phenyltridecane
19. 2-phenyltridecane



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

A00301



Agilent Technologies