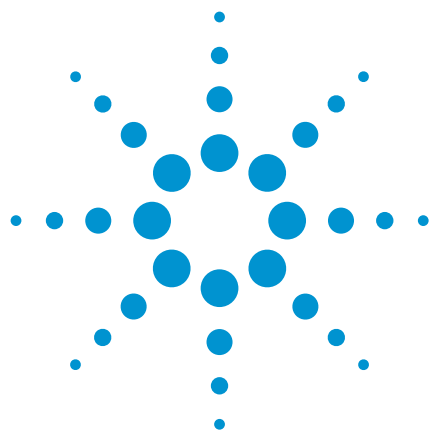




Nitroaromatics and cyclic ketones

Analysis of nitroaromatics and cyclic ketones to EPA 8080

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 13 CB column separates nine nitroaromatics and cyclic ketones according to EPA 8090 in 33 minutes.



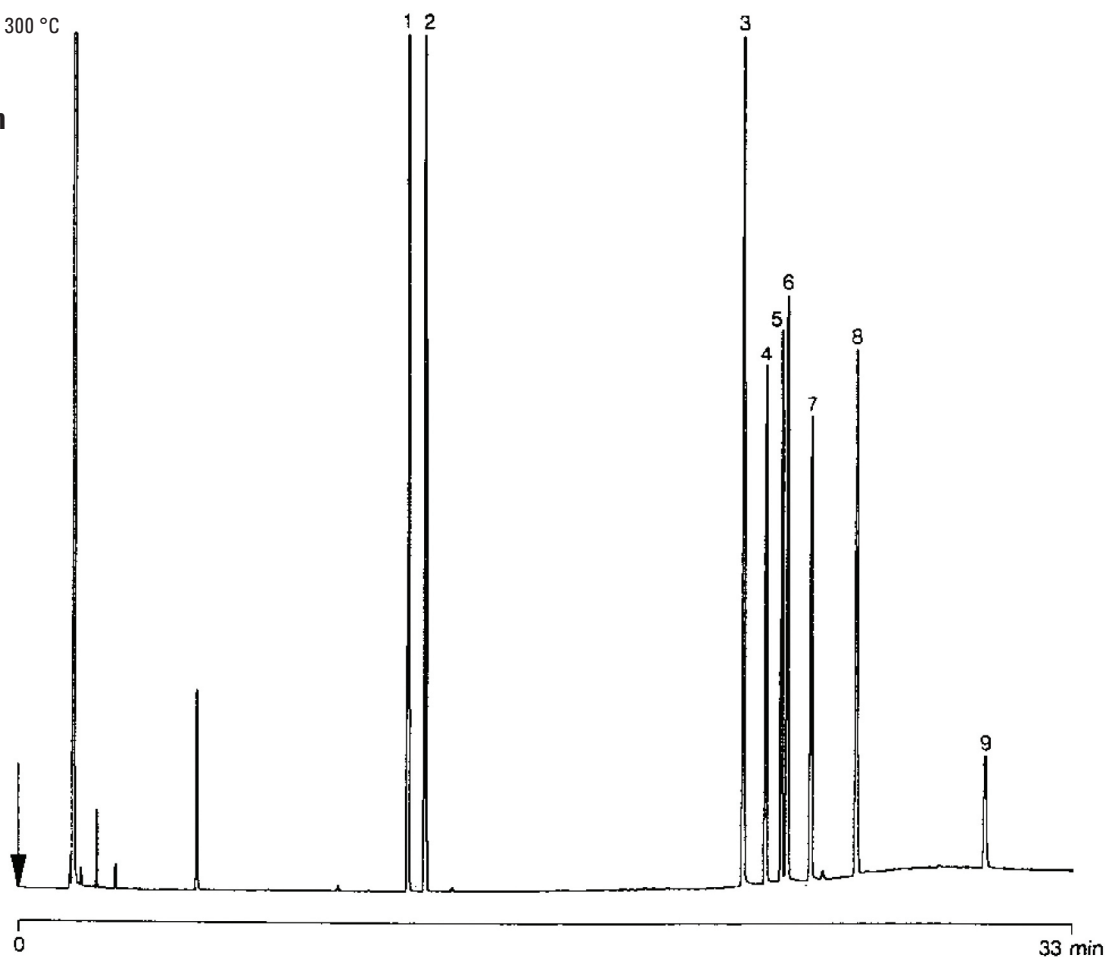
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 13 CB, 0.32 mm x 50 m fused silica
WCOT CP-Sil 13 CB (df = 1.2 µm) (Part no. CP7947)
Temperature : 80 °C → 200 °C, 5 °C/min
Carrier Gas : H₂, 100 kPa (1 bar, 14 psi)
Injector : Split, 100 mL/min
T = 265 °C
Detector : FID
T = 300 °C

Peak identification

1. nitrobenzene
2. isophorone
3. 1,4-naphthoquinone
4. 1,4-dinitrobenzene
5. 1,3-dinitrobenzene
6. 2,6-dinitrotoluene
7. 1,2-dinitrobenzene
8. 2,4-dinitrotoluene
9. 1,2-naphthoquinone



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

A00800



Agilent Technologies