



Nitroaromatics and cyclic ketones

Analysis of nitro aromatics and cyclic ketones to EPA 8090

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

GC/MS analysis of eight nitroaromatics and cyclic ketones according to EPA 8090, with an Agilent FactorFour VF-200ms column, is achieved in less than 20 minutes.



Agilent Technologies

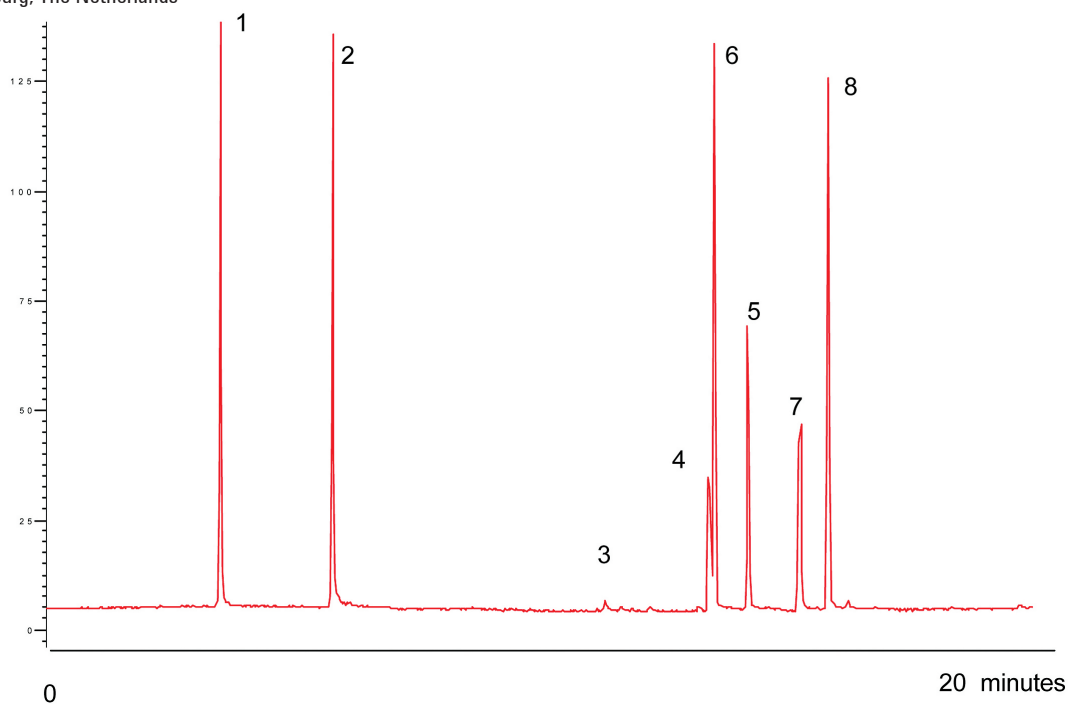
Conditions

Technique : GC
Column : Agilent FactorFour VF-200ms, 0.25 mm x 30 m
(df = 0.25 μ m) (Part no. CP8858)
Temperature : 45 °C, 10 °C/min \rightarrow 325 °C
Carrier Gas : Helium, ca. 1.0 mL/min
Pressure program : 60 kPa
Injector : Split/Splitless, in split mode, 1:100
Detector : FID
Sample Size : 1 μ L
Solvent : methylene chloride, 2000 μ g/mL

Courtesy : Jan Peene, Agilent Application Laboratory,
Middelburg, The Netherlands

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



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

A02097



Agilent Technologies