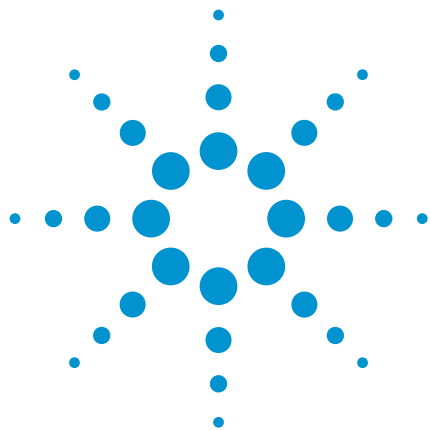




Nitrosamines

Analysis of nitrosoamines to EPA 607

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Wax 51 for Amines column separates nine nitrosamines according to EPA 607 in 18 minutes.



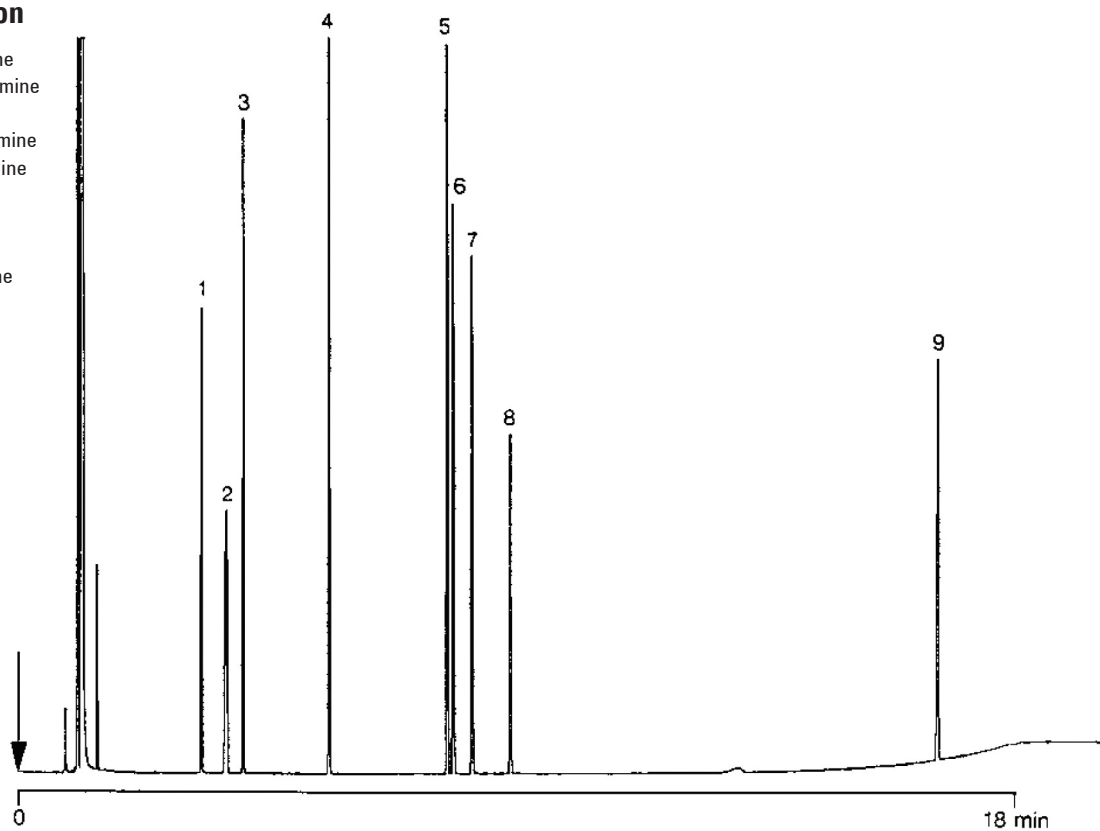
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Wax 51 for Amines, 0.25 mm x 25 m
fused silica WCOT CP-Wax 51 (df = 0.2 µm)
(Part no. CP7405)
Temperature : 60 °C → 240 °C, 10 °C/min
Carrier Gas : H₂, 75 kPa (0.75 bar, 10.5 psi)
Injector : Split, 100 mL/min
T = 265 °C
Detector : FID
T = 300 °C

Peak identification

1. N-nitrosodimethylamine
2. N-nitrosomethylethylamine
3. N-nitrosodiethylamine
4. N-nitroso di-n-propylamine
5. N-nitroso di-n-butylamine
6. N-nitrosopiperidine
7. N-nitrosopyrrolidine
8. N-nitrosomorpholine
9. N-nitrosodiphenylamine



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

A00785



Agilent Technologies