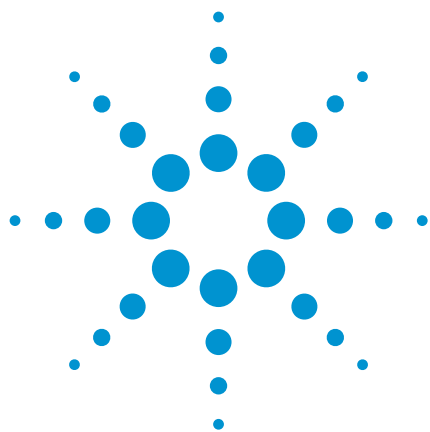




Amines, $C_3 - C_{14}$

Analysis of alkyl and aromatic amines

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Wax for Volatile Amines and Diamines column separates 15 alkyl and aromatic amines in 12 minutes.



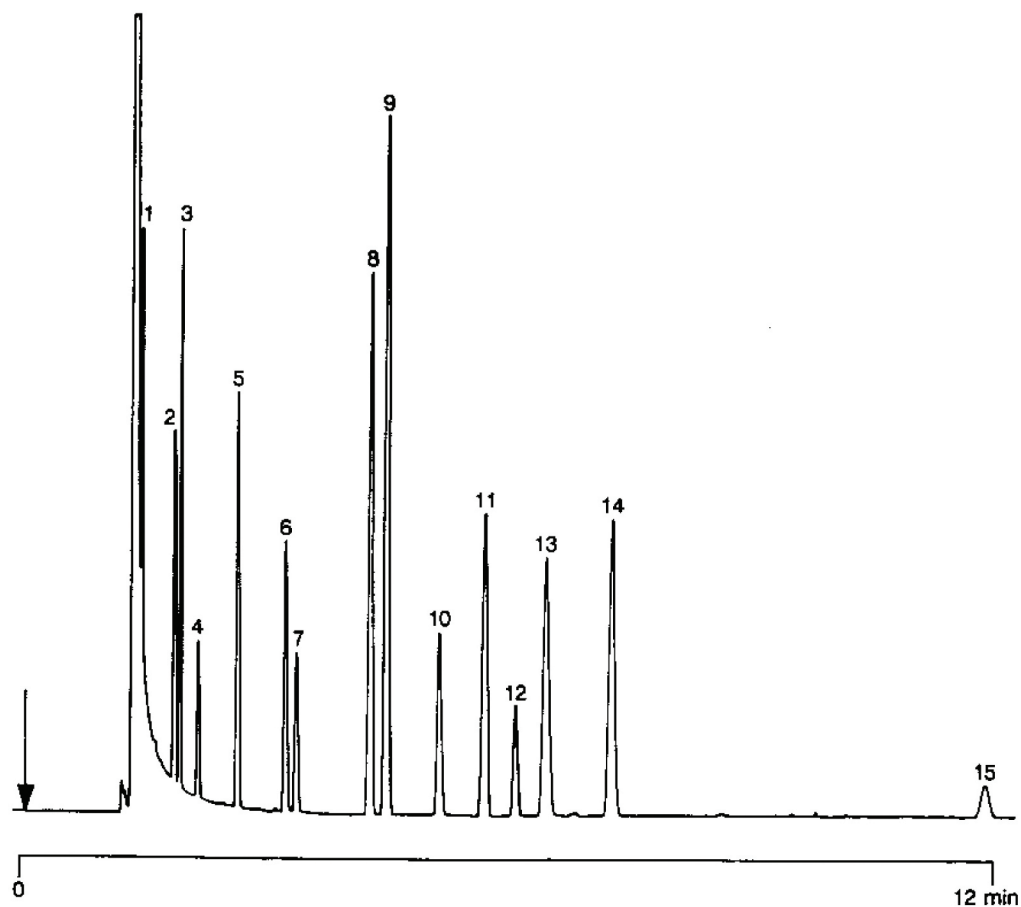
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Wax for volatile amines and diamines,
0.32 mm x 25 m fused silica WCOT CP-Wax
(df = 1.2 μ m) (Part no. CP7422)
Temperature : 170 °C
Carrier Gas : H₂, 40 kPa (0.4 bar, 5.7 psi)
Injector : Split, 50mL
Detector : FID, 16 x 10⁻¹² Afs
Sample Size : 0.02 μ L
Concentration Range : 0.1 %
Solvent Sample : methanol

Peak identification

1. tri-n-propylamine
2. tri-n-butylamine
3. pyridine
4. octylamine
5. N,N-dimethylbenzylamine
6. di-n-hexylamine
7. decylamine
8. N,N-dimethylaniline
9. benzylamine
10. N,N-diethylaniline
11. dicyclohexylamine
12. dodecylamine
13. tri-n-hexylamine
14. aniline
15. tetradecylamine



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

A00701



Agilent Technologies