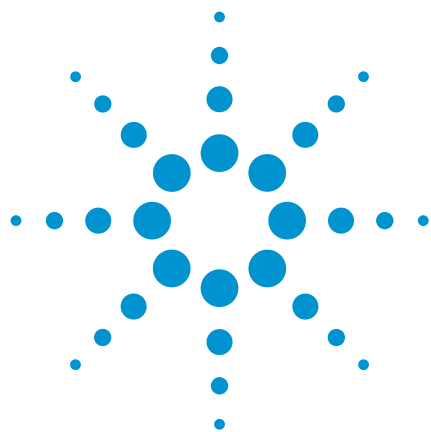




Amines, C₆ – C₁₂

Analysis of alkylamines and aromatic amines

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography using an Agilent CP-Wax 51 for Amines column separates six alkylamines and aromatic amines in six minutes.



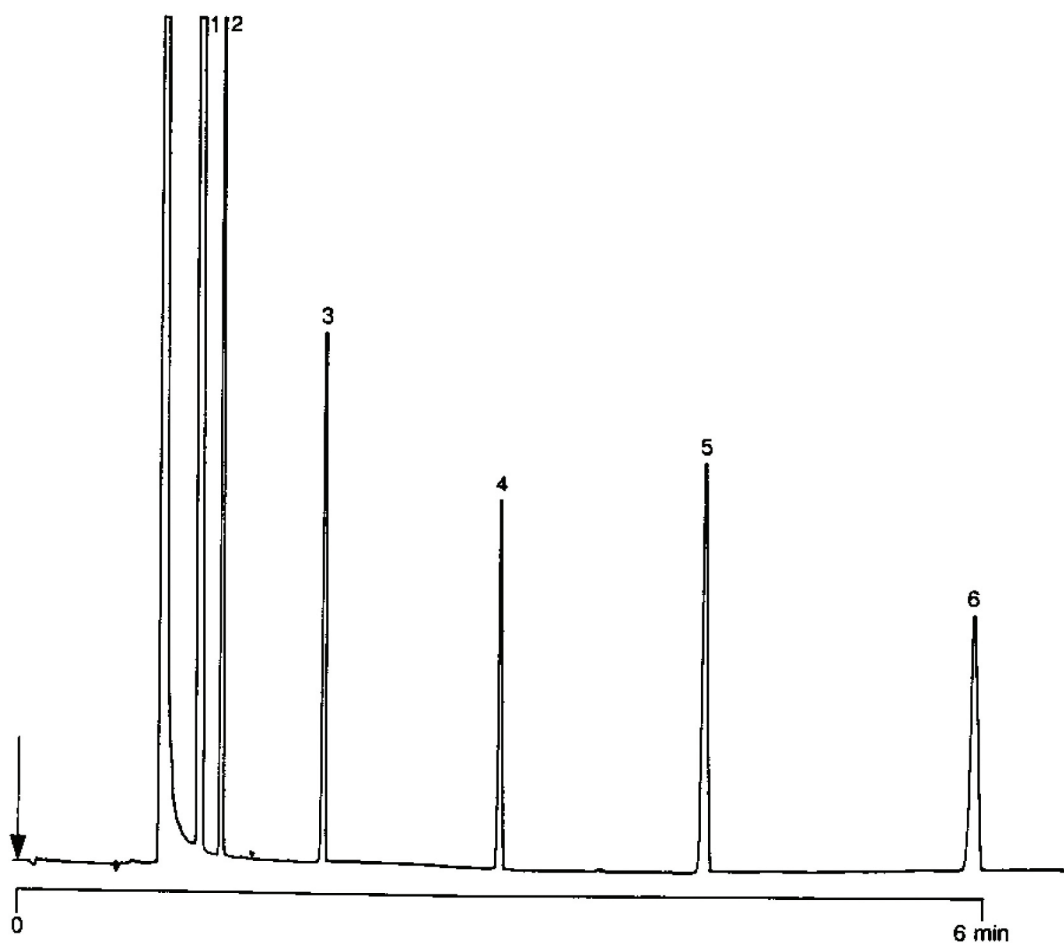
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Wax 51 for Amines, 0.25 mm x 25 m
fused silica WCOT CP-Wax
(df = 0.2 µm) (Part no. CP7405)
Temperature : 130 °C
Carrier Gas : N₂, 120 kPa (1.2 bar, 17 psi)
Injector : Splitter, 100 mL/min
Detector : FID, 8 x 10⁻¹² Afs

Peak identification

1. 2-aminooctane
2. 1-aminooctane
3. 1-aminodecane
4. dicyclohexylamine
5. aniline
6. 2,6-dimethylaniline



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

A00698



Agilent Technologies