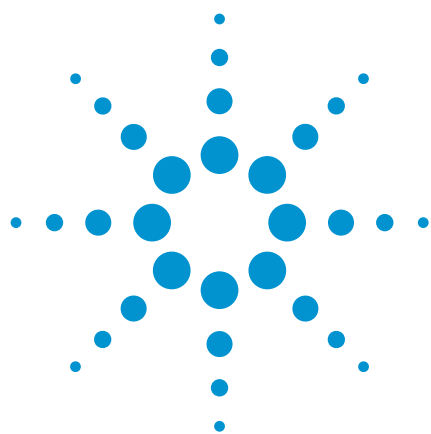




Amines, C₆ – C₁₂

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 8 CB for Amines column determines trace levels of amines in 25 minutes.



Agilent Technologies

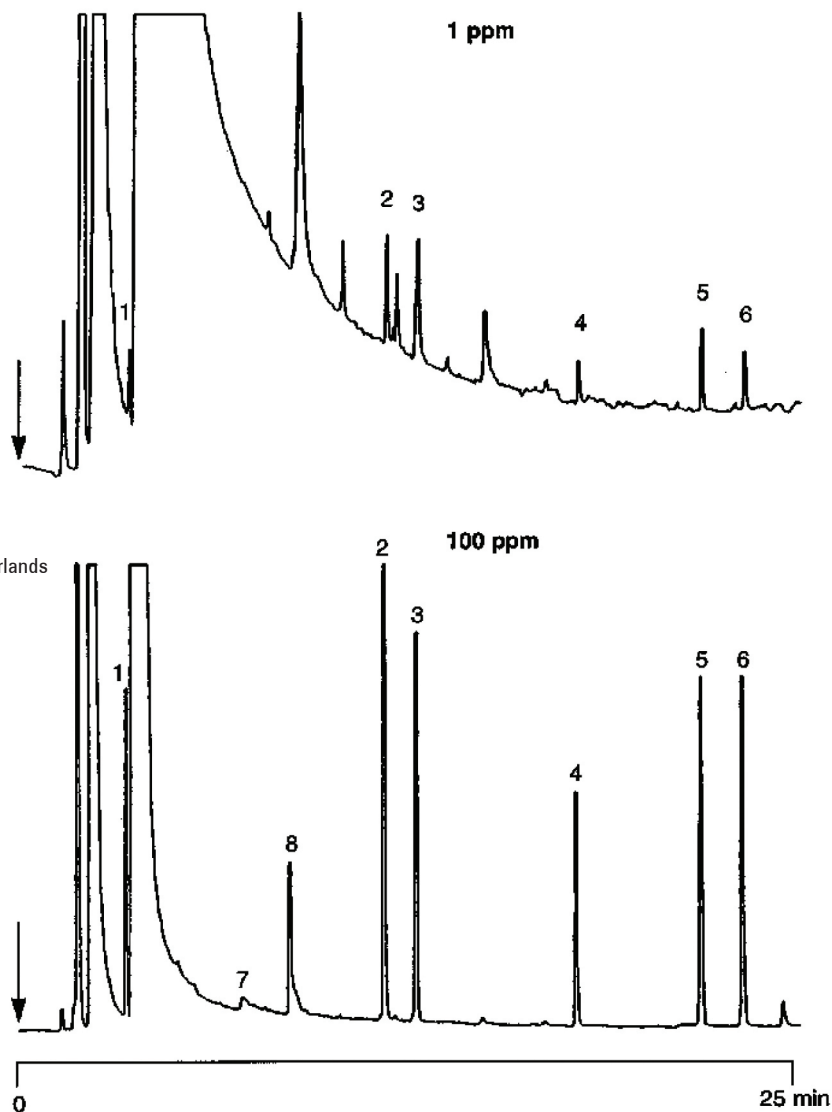
Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 8 CB for Amines,
0.32 mm x 30 m fused silica WCOT
(df = 0.5 μ m) (Custom-made)
Temperature : 40 °C (0 min) $\rightarrow$ 100 °C, 1 °C/min
Carrier Gas : He
Injector : Split, 1:23
T = 225 °C
Detector : FID
T = 300 °C
Concentration Range : 1 ppm and 100 ppm
Solvent Sample : DAB:pyridine, 1:1

Courtesy : Sjoerd van der Wal,
Brigitte Lefrere and Leo Voragen
DSM Research, Geleen, The Netherlands

Peak identification

1. cyclohexylamine
2. 2,6-dimethylaniline
3. dodecane
4. methyl dodecanoate
5. dicyclohexylamine
6. dodecylamine
7. diethylenetriamine
8. 1,6-diaminohexane



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

A01278



Agilent Technologies