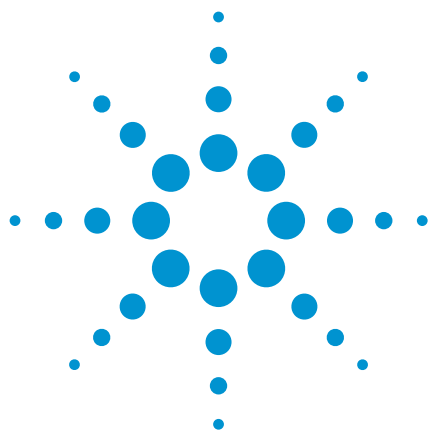




Amino alcohols, C₂ - C₅

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 8 CB for Amines column separates six C₂ to C₅ amino alcohols in 30 minutes.



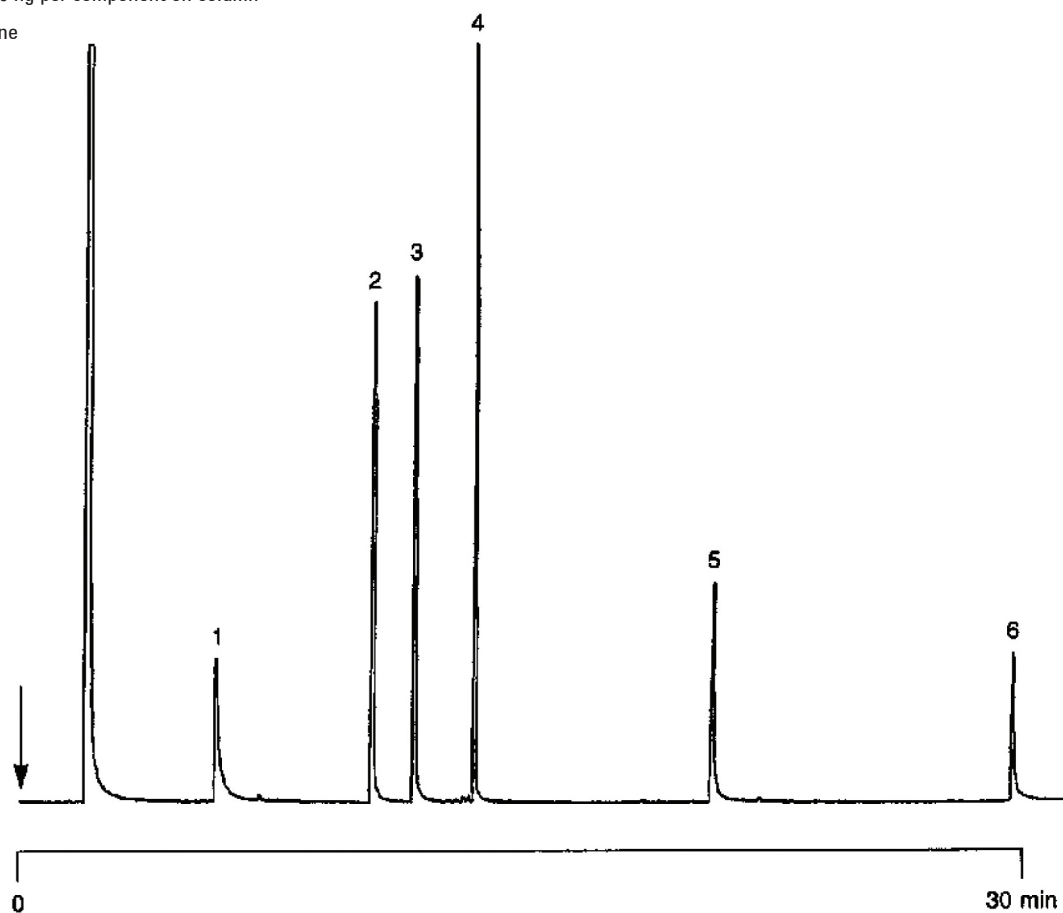
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 8 CB for Amines, 0.25 mm x 30 m
fused silica WCOT CP-Sil 8 CB (df = 0.5 µm)
(Part no. CP7595)
Temperature : 40 °C (5 min) → 200 °C, 6 °C/min
Carrier Gas : H₂, 60 kPa (0.6 bar, 8.6 psi)
Injector : Split, T = 200 °C
Detector : FID, T = 285 °C
Sample Size : 1 µL
Concentration Range : 30 - 50 ng per component on column
Solvent Sample : pyridine

Peak identification

1. ethanolamine
2. 2-ethylaminoethanol
3. 2-aminobutanol
4. diethylaminoethanol
5. diethanolamine
6. triethanolamine



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

A001254



Agilent Technologies