



Amino-alcohols

Analysis of amino-alcohols, C₃-C₅

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Wax for Volatile Amines and Diamines column separates four amino-alcohols in seven minutes.



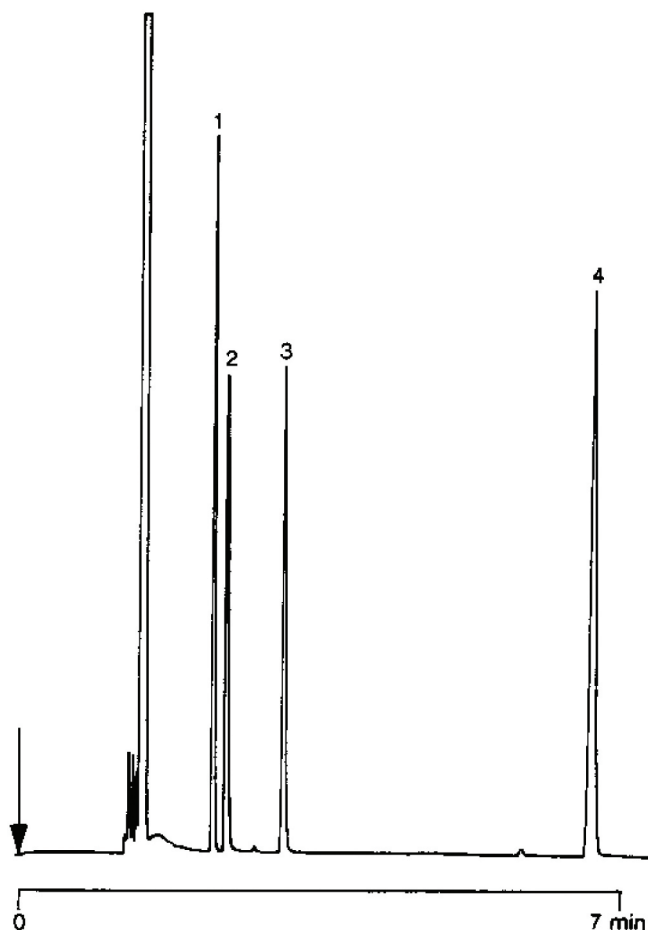
Agilent Technologies

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 µm) (Part no. CP7422)
Temperature : 170 °C
Carrier Gas : H₂, 45 kPa (0.45 bar, 6.4 psi)
Injector : Split, 100 mL
Detector : FIO, 8 x 10⁻¹² Afs
Sample Size : 0.02 µL
Concentration range : 0.03%
Solvent sample : CH₂Cl₂

Peak identification

1. 1-amino-2-propanol
2. ethanolamine
3. 2-amino-1-butanol
4. 5-amino-1-pentanol



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

A00700



Agilent Technologies