



Amines

Analysis of diamines in water

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Wax for Volatile Amines and Diamines column separates five diamines in water in nine minutes.



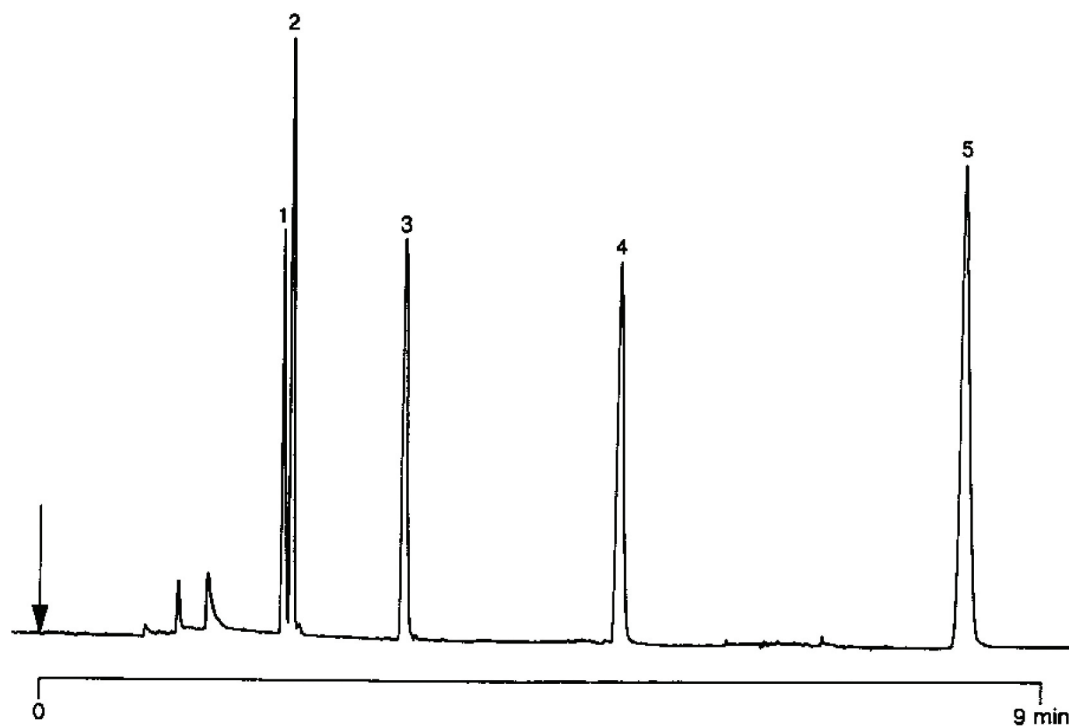
Agilent Technologies

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 : 120 °C
Carrier Gas : N₂, 40 kPa (0.4 bar, 5.7 psi)
Injector : Split, 100 mL/min
Detector : FID, 16 x 10⁻¹² Afs
Sample Size : 0.05 μ L
Concentration range : 0.03%
Solvent sample : H₂O

Peak identification

1. 1,2-diaminoethane
2. 1,2-diaminopropane
3. 1,3-diaminopropane
4. 1,4-diaminobutane
5. 1,5-diaminopentane



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

A00697



Agilent Technologies