



## **Amines**

# **Analysis of diamines on wide-bore column**

## **Application Note**

Environmental

### **Authors**

Agilent Technologies, Inc.

### **Introduction**

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



**Agilent Technologies**

## Conditions

Technique : GC-capillary  
Column : Agilent CP-Wax for volatile amines and diamines,  
0.53 mm x 25 m fused silica WCOT CP-Wax  
(df = 1.2  $\mu$ m) (Part no. CP7422)  
Temperature : 120 °C  
Carrier Gas : N<sub>2</sub>, 10 mL/min  
Injector : Direct  
Detector : FID, 32 x 10<sup>-12</sup> Afs  
Sample Size : 0.01  $\mu$ L  
Concentration range : 0.03%  
Solvent sample : H<sub>2</sub>O

## Peak identification

1. 1,2-diaminoethane
2. 1,2-diaminopropane
3. 1,3-diaminopropane
4. 1-amino-2-propanol
5. ethanolamine
6. 1,4-diaminobutane
7. 1,5-diaminopentane



[www.agilent.com/chem](http://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

A00699



**Agilent Technologies**