



Volatile amino alcohols, C₂ – C₅

Analysis of amino alcohols in water

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography using an Agilent CP-Sil 5 CB column separates five C₂ to C₅ volatile amino alcohols in water in ten minutes.



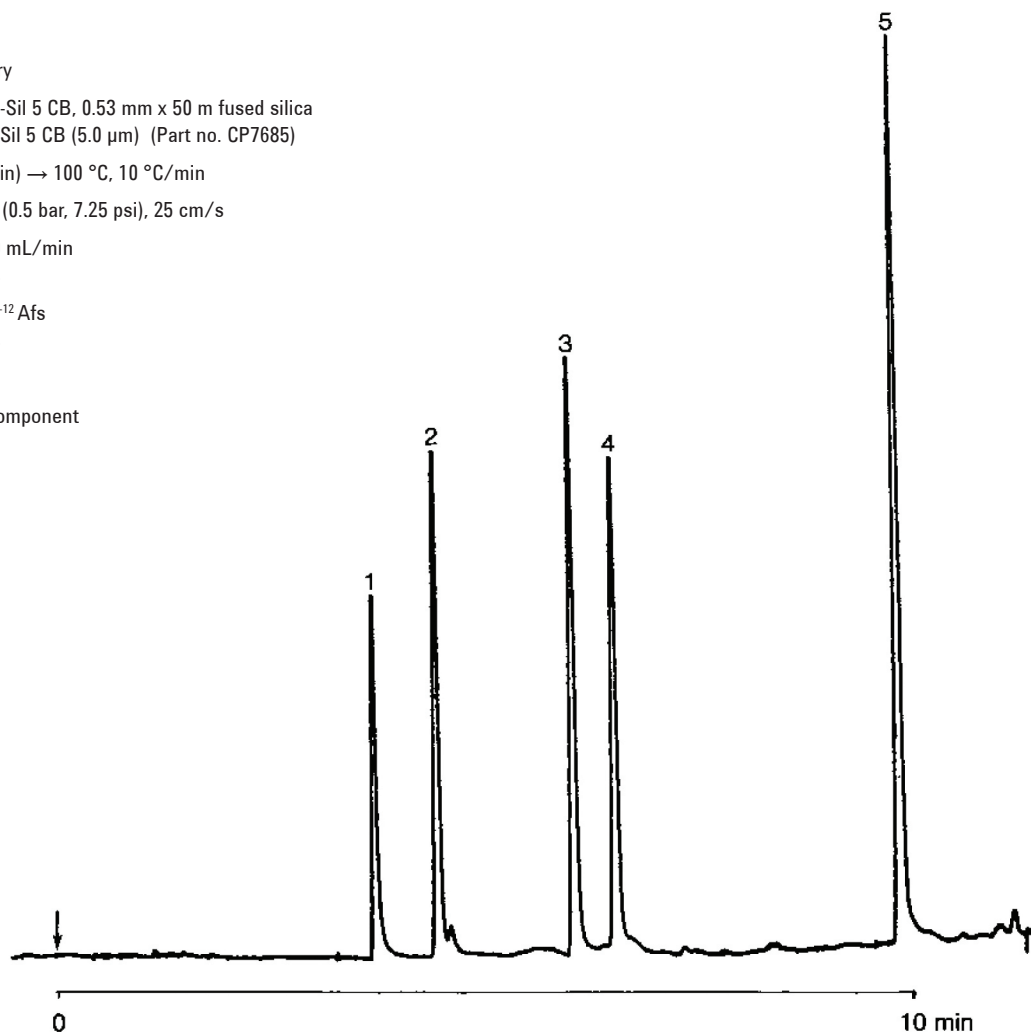
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 5 CB, 0.53 mm x 50 m fused silica
WCOT CP-Sil 5 CB (5.0 μ m) (Part no. CP7685)
Temperature : 65 °C (6 min) $\rightarrow$ 100 °C, 10 °C/min
Carrier Gas : H₂, 50 kPa (0.5 bar, 7.25 psi), 25 cm/s
Injector : Splitter, 50 mL/min
T = 240 °C
Detector : FID, 8 x 10⁻¹² Afs
T = 250 °C
Sample Size : 0.2 μ L
Concentration Range : $\approx$ 0.1 %/component
Solvent Sample : water

Peak identification

1. 2-amino-ethanol
2. 1-amino-2-propanol
3. 2-(ethylamino)-ethanol
4. 2-amino-1-butanol
5. 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

A00163



Agilent Technologies