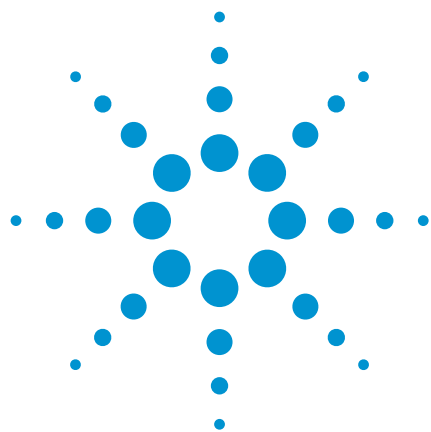




Amines

Analysis of impurities in anhydrous mono- and trimethylamine

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with a custom-made Agilent PoraPLOT Amines column separates three impurities in anhydrous mono- and trimethylamine in 4.3 minutes.



Agilent Technologies

Conditions

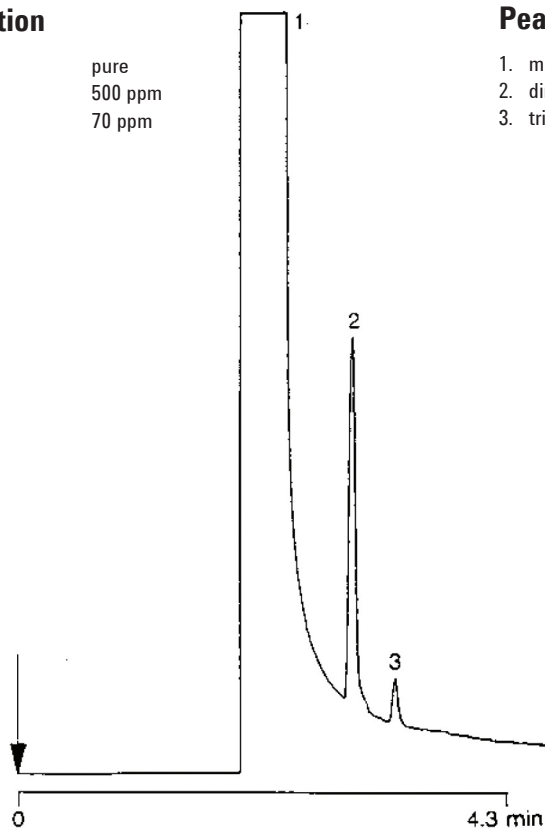
Technique : GC-capillary
Column : Agilent PoraPLOT, 0.32 mm x 25 m amines fused silica PLOT PoraPLOT (df = 10 µm)
(Part no. CP7591)
Temperature : 150 °C
Carrier Gas : He. 140 kPa (1.4 bar, 20 psi)
Injector : Splitter
Detector : FID
Sample Size : 0.1 µL

Impurities in monomethylamine

Flow: 15 mL/min

Peak identification

1. monomethylamine	pure
2. dimethylamine	500 ppm
3. trimethylamine	70 ppm

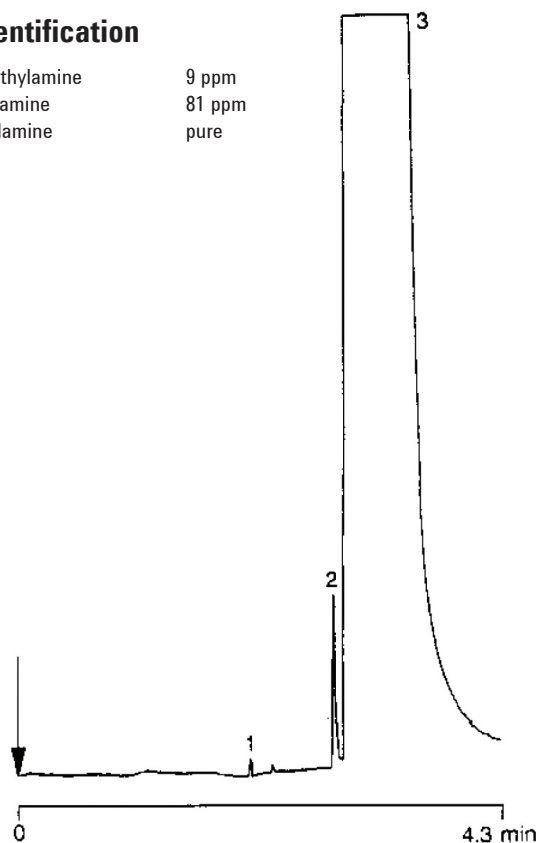


Impurities in trimethylamine

Flow: 2.5 mL/min

Peak identification

1. monomethylamine	9 ppm
2. dimethylamine	81 ppm
3. trimethylamine	pure



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

A00762



Agilent Technologies