



# Anilines

## Application Note

Environmental

### Authors

Agilent Technologies, Inc.

### Introduction

HPLC analysis of anilines is achieved with an Agilent Polaris C18-A column in less than 12 minutes.



**Agilent Technologies**

## Conditions

Column : Agilent Polaris C18-A, 4.6 x 150 mm, 5  $\mu$ m  
Part no. A2000150X046

Eluent : H<sub>2</sub>O 30:70

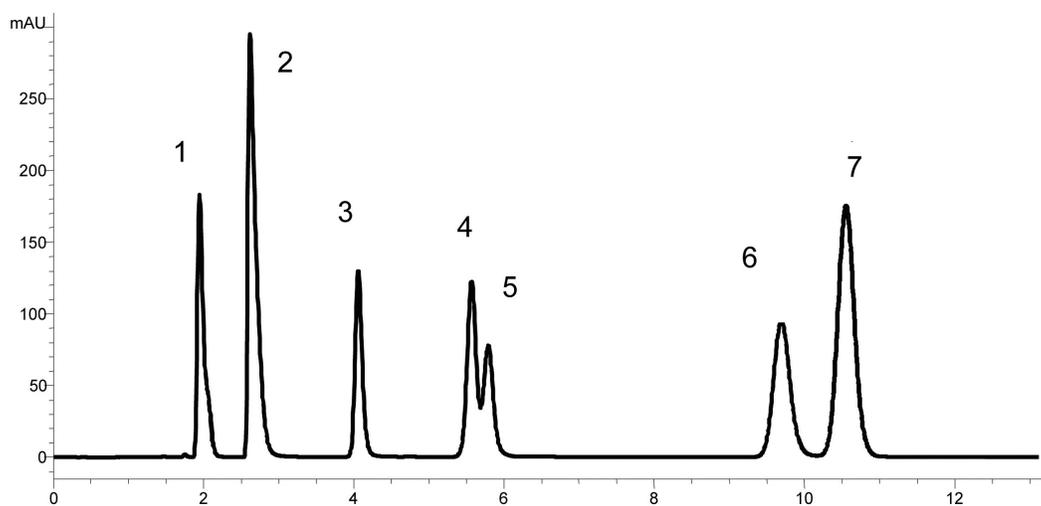
Temp : ambient

Flow : 1.0 mL/min

Detection : UV @ 220 nm

## Peak identification

1. caffeine
2. pyridine
3. aniline
4. o-toluidine
5. 4-nitro-aniline
6. 2-ethylaniline
7. 2-nitroanisole



[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

A02015



**Agilent Technologies**