



Aniline homologs

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Agilent Pursuit C18 demonstrates absence of silanol activity. Aniline homologs are separated, based on hydrophobic substitution.



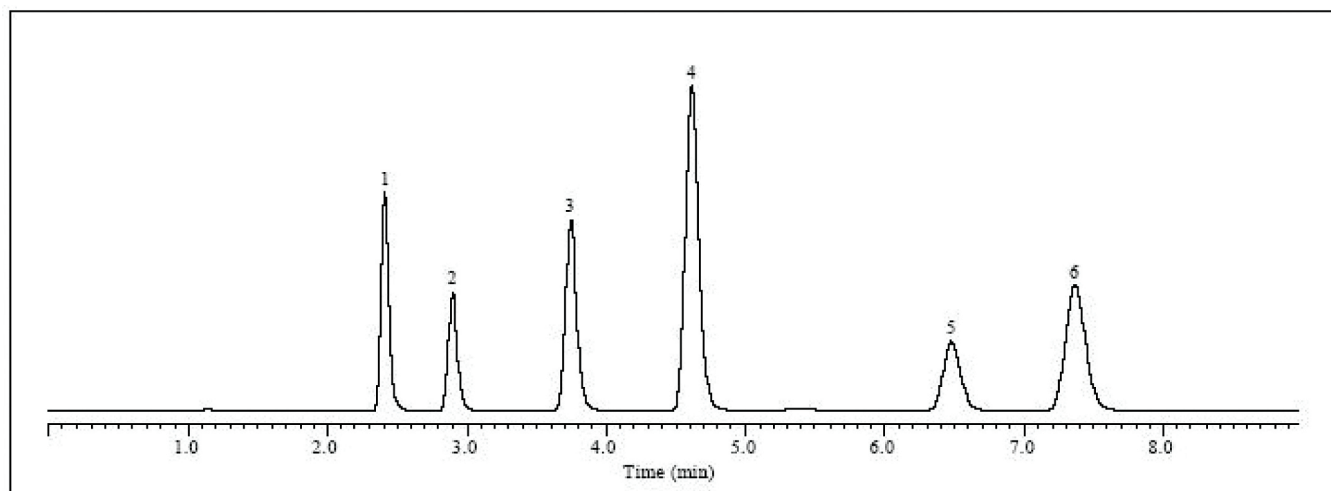
Agilent Technologies

Conditions

Technique : HPLC
Column : Agilent Pursuit C18, 4.6 x 150 mm, 5 μ m
Part no. : Conventional A3000150X046
ChromSep Cartridge System A3000150C046
Temperature : Ambient
Eluent : A: MeOH
B: Water
A:B = 60:40
Flow rate : 1.0 mL/min
Detection : UV @ 254 nm

Peak identification

1. aniline
2. o-toluidine
3. 2-ethylaniline
4. n-ethylaniline
5. n,n-dimethylaniline
6. n-propylaniline



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

A02310



Agilent Technologies