



Barbiturates

Application Note

Forensic Toxicology

Authors

Agilent Technologies, Inc.

Introduction

HPLC with Agilent Pursuit C18 5 μm material in a ChromSep cartridge analyzes eight barbiturates in 18 minutes.



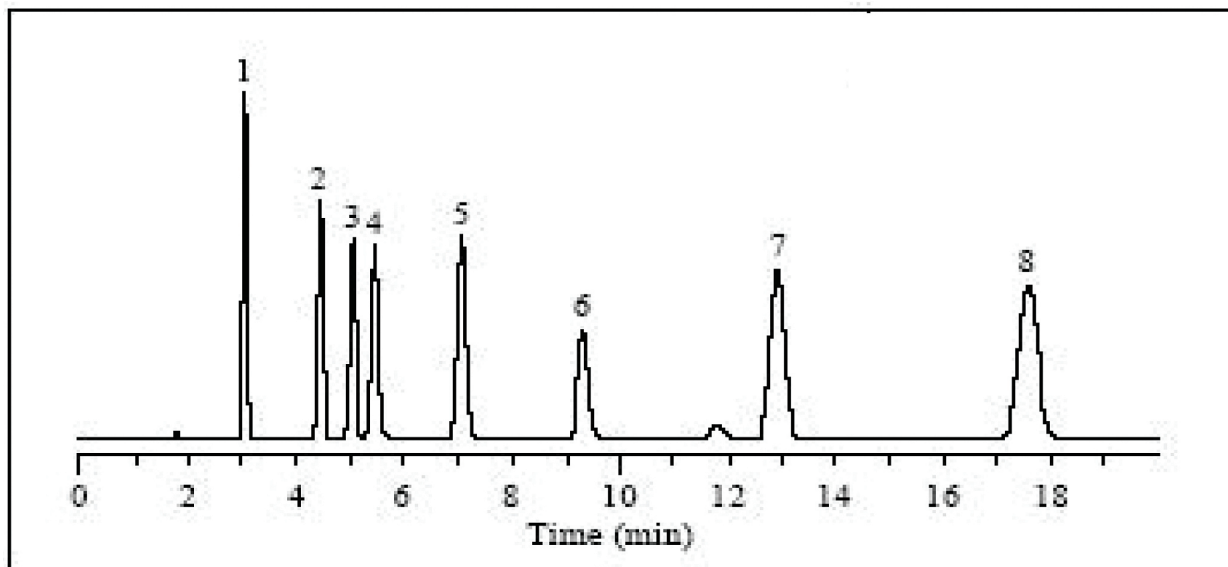
Agilent Technologies

Conditions

Technique : HPLC
Column : Agilent Pursuit C18, 4.6 x 150 mm, 5 µm
Part no. : Conventional A3000150X046
ChromSep Cartridge System A3000150C046
Temperature : Ambient
Eluent : A: MeOH
B: Water
A:B = 45:55
Flow rate : 1.0 mL/min
Detection : UV @ 214 nm

Peak identification

1. barbital
2. allobarbitol
3. phenobarbital
4. aprobarbital
5. butabarbital
6. mephobarbital
7. pentobarbital
8. secobarbital



www.agilent.com/chem

For Forensic Use only.

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

A02313



Agilent Technologies