



Anabolic steroids

Application Note

Forensic Toxicology

Authors

Agilent Technologies, Inc.

Introduction

Analyze steroids effectively with Agilent Pursuit PFP HPLC columns.



Agilent Technologies

Conditions

Column : Agilent Pursuit PFP, 4.6 mm x 150 mm 5 µm
Part No. A3050150X046

Mobile Phase : Acetonitrile: Water- 45:55

Flow Rate : 1.5 mL/min

Temperature : Ambient

Detection : 220 nm

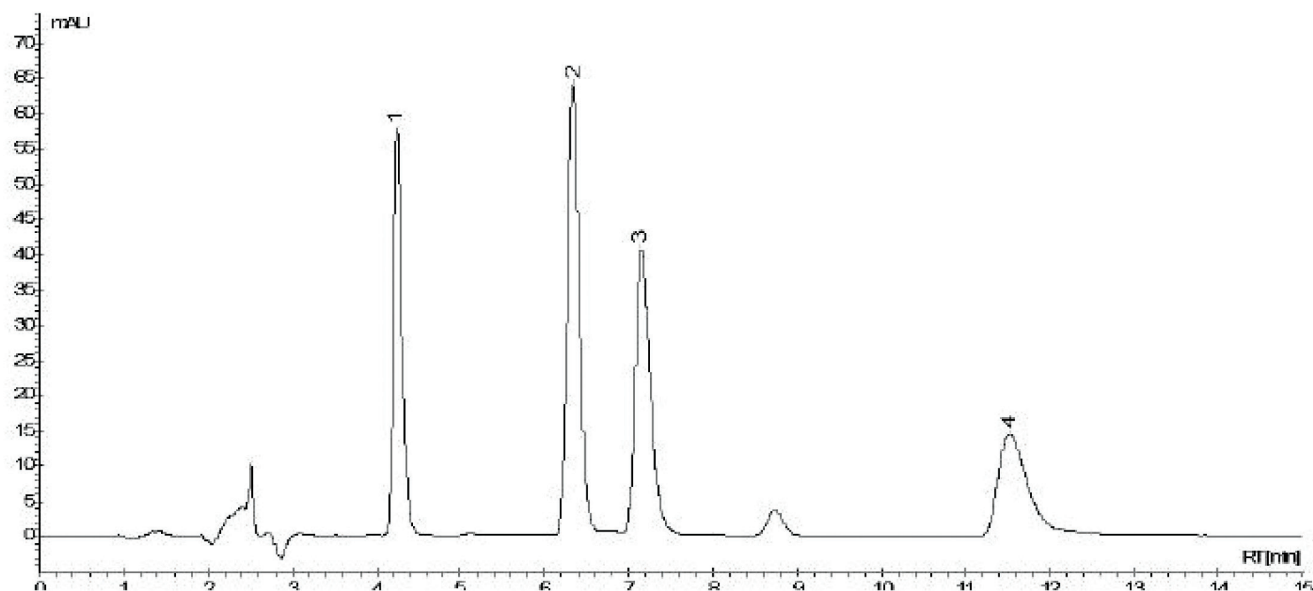
Injection Vol. : 5 µL

Sample Concentration : 1 mg/mL all components

Solvent Sample : Water

Peak identification

1. fluoxymesterone
2. methyltestosterone
3. stanozolol
4. oxymetholone



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

A02456



Agilent Technologies