

Steroids used in hormone replacement therapies

Application Note

Clinical Research

Authors

Agilent Technologies, Inc.

Introduction

Agilent Pursuit PFP columns are ideal for the HPLC analysis of steroids when developing drugs.



Agilent Technologies

Conditions

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

Mobile Phase : Acetonitrile: Water + 0.1% HCOOH, 40:60

Flow Rate : 0.8 mL/min

Temperature : Ambient

Detection : 220 nm

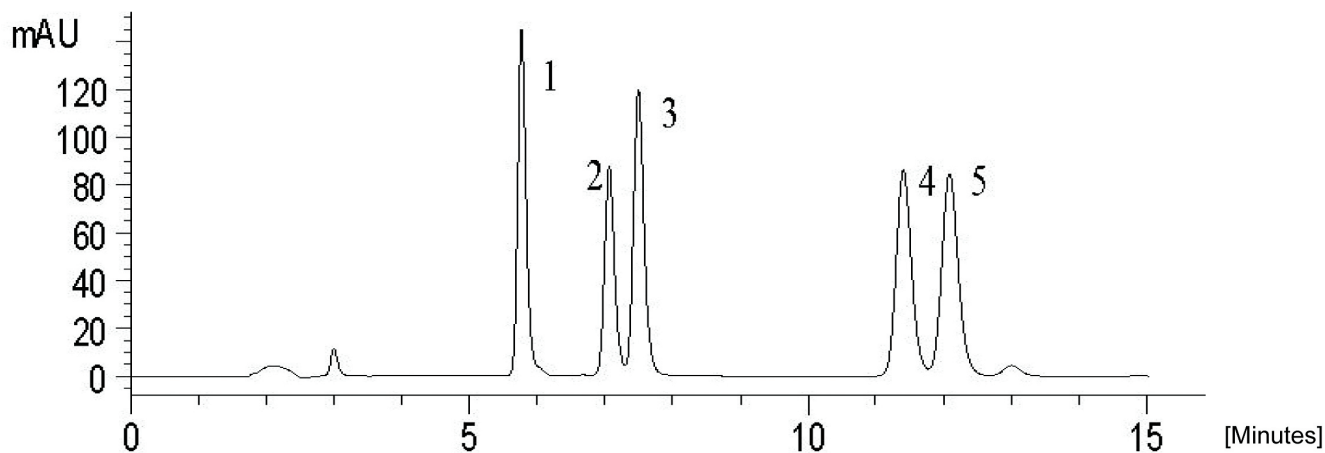
Injection Vol. : 5 μ L

Sample Concentration : 1 mg/mL all components

Solvent Sample : Water

Peak identification

1. norethindrone
2. 17- α -ethynylestradiol
3. d (-)- norgestrel
4. progesterone
5. 19-norethindrone acetate



www.agilent.com/chem

For Research Use Only. Not for use in diagnostic procedures.

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

A02453



Agilent Technologies