

Application Report 06

Steroids on Discovery HS PEG

Norethindrone and Norethindrone Acetate are oral contraceptives and are used to treat a variety of reproductive illnesses in female patients.

Key Words

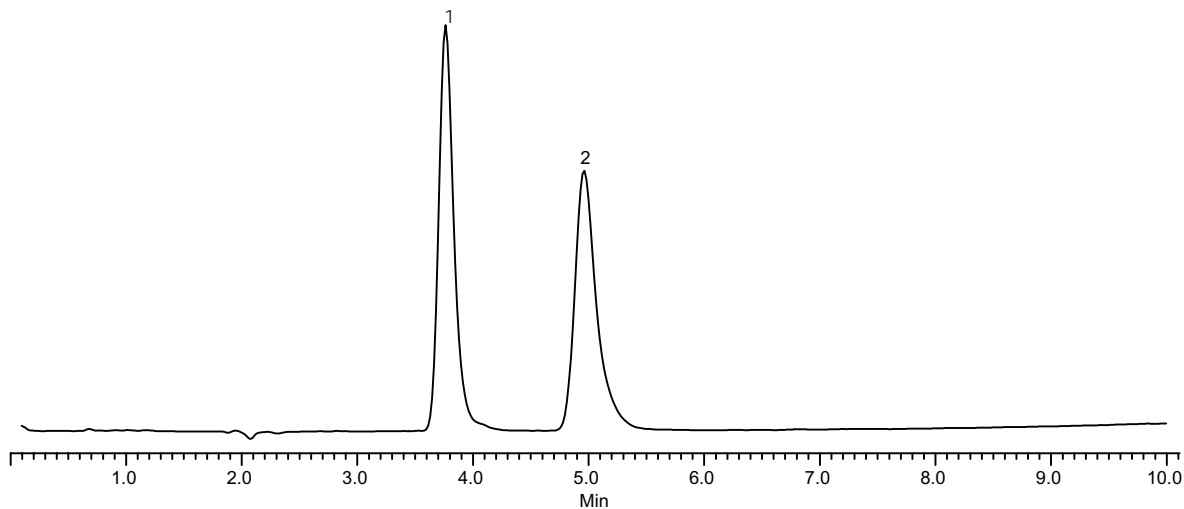
HS PEG, oral contraceptives, norethindrone, norethindrone acetate, 567416-U

Author: Carmen T. Santasania

Raw Data File Name: mx4

Acquisition System: W-2690

Notebook Reference:1458-43



G002039

Conditions

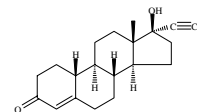
Column: Discovery HS PEG, 15cm x 4.6mm ID, 5µm
Cat. No.: 567416-U
Mobile Phase: A:Water/B:Methanol 60/40
Flow Rate: 1.0mL/min
Temperature: 30°C
Detection: UV at 254nm
Injection Volume: 10µL
Sample: 25µg/mL of each in 50/50 water/methanol

Peak IDs

1. Norethindrone
2. Norethindrone acetate

Structures

Norethindrone - G002040



Norethindrone acetate - G002041

