

Application Report 07

Steroids on Discovery C18

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

Key Words

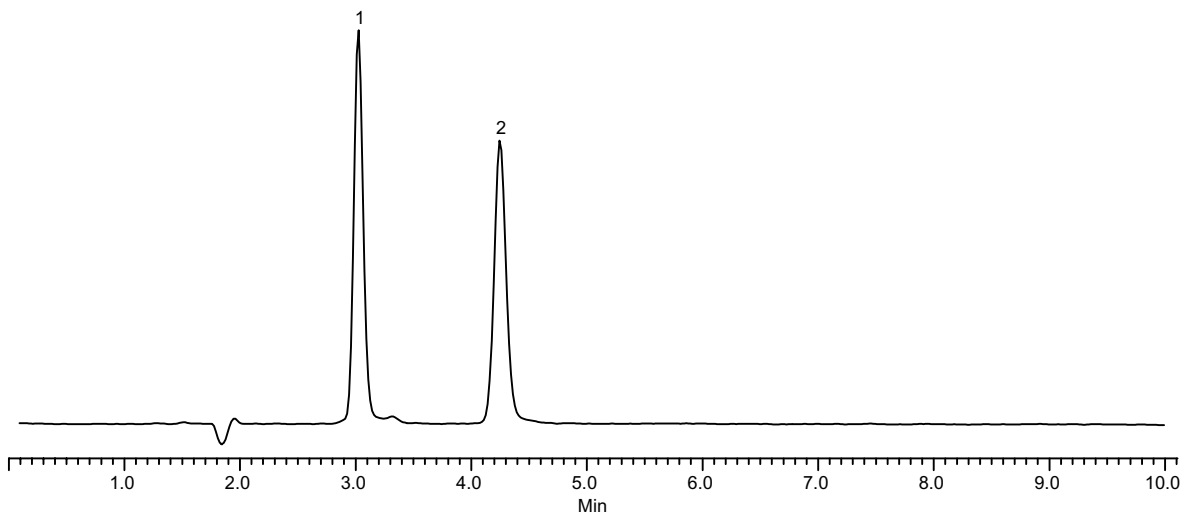
Discovery C18, oral contraceptives, norethindrone, norethindrone acetate, 504955

Author: Carmen T. Santasania

Raw Data File Name: mx19c18

Acquisition System: W-2690

Notebook Reference:1458-45



G002042

Conditions

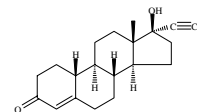
Column: Discovery C18 column, 15cm x 4.6mm ID, 5µm
Cat. No.: 504955
Mobile Phase: A:Water/B:Methanol 30/70
Flow Rate: 1.0mL/min
Temperature: 30°C
Detection: UV at 254nm
Injection Volume: 10µL
Sample: 25µg/mL of each in 50/50 water/methanol

Peak IDs

1. Norethindrone
2. Norethindrone acetate

Structures

Norethindrone - G002040



Norethindrone acetate - G002041

