

# Application Report 319

## Analysis of Anabolic Androgenic Steroids Using Ascentis™ C8

This application demonstrates the suitability of the Ascentis C8 for the analysis of anabolic androgenic steroids. Structures along with the optimized chromatogram obtained on the Ascentis C8 are presented.

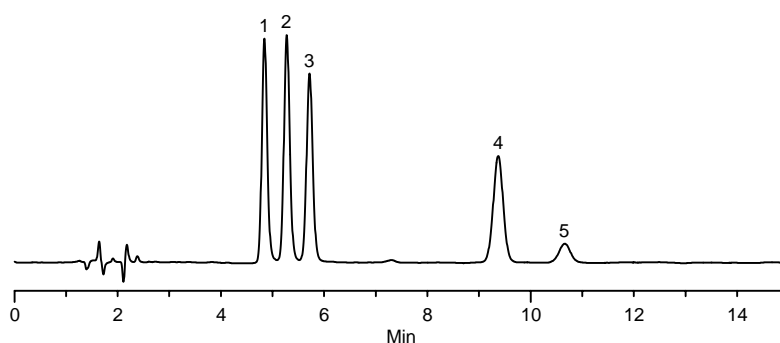
### Key Words

nandrolone, 46476, 434-22-0, methandienone, 55609, 72-63-9, testosterone, 46923, 58-22-0, norethandrolone, 52-78-8, oxymetholone, O0380, 434-07-1, Ascentis C8, 581424-U

Author: Hugh M. Cramer

Acquisition System: Hitachi Mt Fuji

Notebook Reference: 1558



G003238

### Conditions

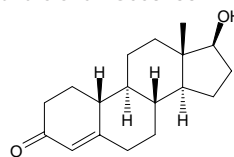
column: Ascentis C8, 15 cm x 4.6 mm I.D., 5 µm particles (581424-U)  
mobile phase: 50:50, water:acetonitrile  
flow rate: 1.0 mL/min.  
temp.: 35 °C  
det.: UV at 254 nm  
injection: 10 µL  
sample: as indicated in mobile phase

### Peak IDs

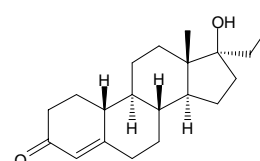
1. Nandrolone (50 µg/mL)
2. Methandienone (50 µg/mL)
3. Testosterone (50 µg/mL)
4. Norethandrolone (50 µg/mL)
5. Oxymetholone (50 µg/mL)

### Structures

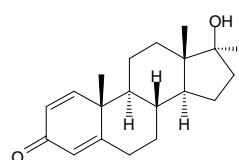
Nandrolone - G003233



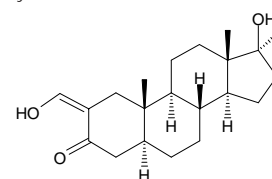
Norethandrolone - G003236



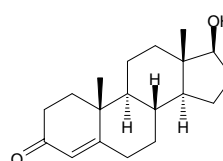
Methandienone - G003234



Oxymetholone - G003237



Testosterone - G003235



Ascentis is a trademark of Sigma-Aldrich Co.