

# Application Report 321

## Analysis of the Anabolic Androgenic Steroid Methandienone and Hydroxy Metabolite Using Ascentis™ C8

This application demonstrates the suitability of the Ascentis C8 for the analysis of the anabolic androgenic steroid methandienone and its metabolite 6- $\beta$ -hydroxymethandienone. Structures along with the optimized chromatogram obtained on the Ascentis C8 are presented.

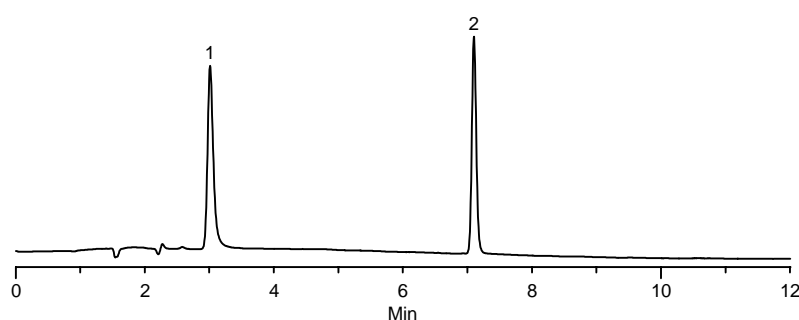
### Key Words

methandienone, 55609, 72-63-9, 6- $\beta$ -hydroxymethandienone, Ascentis C8, 581424-U

Author: Hugh M. Cramer

Acquisition System: Hitachi Mt Fuji

Notebook Reference: 1558



G003241

### Conditions

column: Ascentis C8, 15 cm x 4.6 mm I.D., 5  $\mu$ m particles (581424-U)  
mobile phase A: water  
mobile phase B: acetonitrile  
flow rate: 1.0 mL/min.  
temp.: 35 °C  
det.: UV at 254nm  
injection: 10  $\mu$ L  
sample: as indicated 60:40, water:acetonitrile  
gradient:

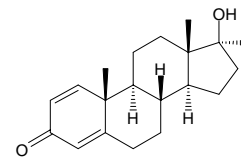
Min	%A	%B
0	60	40
2	60	40
10	0	100
12	0	100

### Peak IDs

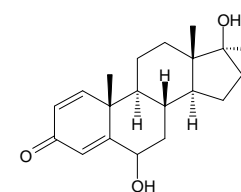
1. 6- $\beta$ -hydroxymethandienon (50  $\mu$ g/mL)
2. Methandienone (50  $\mu$ g/mL)

### Structures

6- $\beta$ -hydroxymethandienon - G003234



Methandienone - G003240



Ascentis is a trademark of Sigma-Aldrich Co.