

# Quantitative analysis of estradiol and testosterone in plasma for clinical research using the TSQ Altis triple quadrupole mass spectrometer

## Authors

Xiaolei Xie, Kristine Van Natta,  
Neloni Wijeratne, Claudia Martins

Thermo Fisher Scientific,  
San Jose, CA

## Keywords

Estradiol, LC-MS/MS, LLE,  
testosterone, TSQ Altis MS

## Goal

To develop a sensitive LC-MS/MS method for quantitative analysis of estradiol and testosterone in plasma for clinical research using liquid chromatographic separation coupled to a triple quadrupole mass spectrometer.

## Introduction

Analysis of estradiol and testosterone in plasma samples for clinical research requires a sensitive analytical method. Liquid chromatography coupled with tandem mass spectrometry (LC-MS/MS) has been widely adopted as an analytically sensitive and selective technique for estradiol and testosterone analysis in complex matrices such as human serum or plasma.

## Experimental

### Sample preparation

To prepare the samples, 10  $\mu\text{L}$  of spiking solution (final concentration range: 0.5–10 ng/mL) and 20  $\mu\text{L}$  of internal standard (2 ng/mL testosterone- $^{13}\text{C}_3$  and 20 ng/mL estradiol-d5) were added to 400  $\mu\text{L}$  of plasma. Then 2 mL of MTBE were added and the sample was vortexed. After liquid-liquid extraction (LLE), the MTBE layer was evaporated under nitrogen, and 200  $\mu\text{L}$  of 50:50 methanol/water were added to reconstitute. From this, 10  $\mu\text{L}$  were injected in triplicate for LC-MS/MS analysis.

## Liquid chromatography

Chromatographic separation was performed using a Thermo Scientific™ Vanquish™ Flex Binary HPLC system equipped with a Thermo Scientific™ Accucore™ aQ C18 Polar Endcapped LC column (100 × 2.1 mm, 2.6 μm particle size, P/N 17326-102130). Mobile phases A and B were 0.5 mM ammonium fluoride in Fisher Chemical™ Optima™ grade water and pure methanol, respectively. The column temperature was 40 °C. The total run time was 9 minutes (Table 1).

Table 1. LC gradient.

| No | Time (min) | Flow (mL/min) | %B  | Curve |
|----|------------|---------------|-----|-------|
| 1  | 0          | 0.25          | 30  | 5     |
| 2  | 1          | 0.25          | 30  | 5     |
| 3  | 1.5        | 0.25          | 55  | 5     |
| 4  | 5          | 0.25          | 85  | 5     |
| 5  | 6          | 0.25          | 100 | 5     |
| 6  | 7          | 0.25          | 100 | 5     |
| 7  | 7.01       | 0.25          | 30  | 5     |
| 8  | 9          | 0.25          | 30  | 5     |

## Mass spectrometry

MS analysis was carried out on a Thermo Scientific™ TSQ Altis™ triple quadrupole mass spectrometer equipped with heated electrospray ionization. Table 2 shows the mass spectrometer source parameters.

Two selected reaction monitoring (SRM) transitions were monitored for estradiol and testosterone and their isotope-labeled internal standards to provide ion ratio

confirmations (IRC). The scans were run in timed selected reaction monitoring (t-SRM) mode with a cycle time of 0.4 seconds. Table 3 shows SRM properties used in this analysis.

Table 2. Source parameters for the TSQ Altis mass spectrometer.

| Ion Source Parameter   | Value  |
|------------------------|--------|
| Positive Ion           | 3500 V |
| Sheath Gas             | 40 Arb |
| Aux Gas                | 12 Arb |
| Sweep Gas              | 1 Arb  |
| Ion Transfer Tube Temp | 350 °C |
| Vaporizer Temp         | 350 °C |

## Data analysis

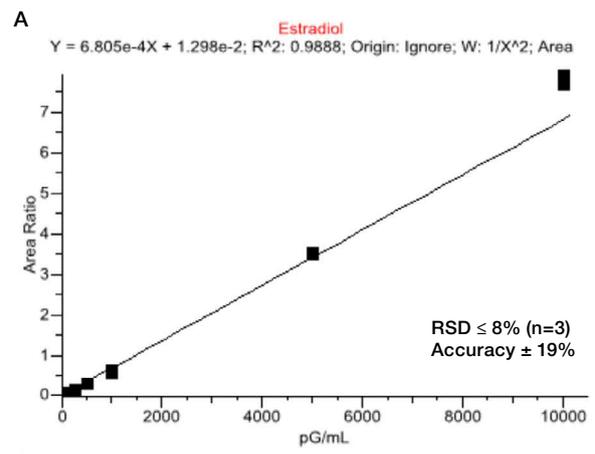
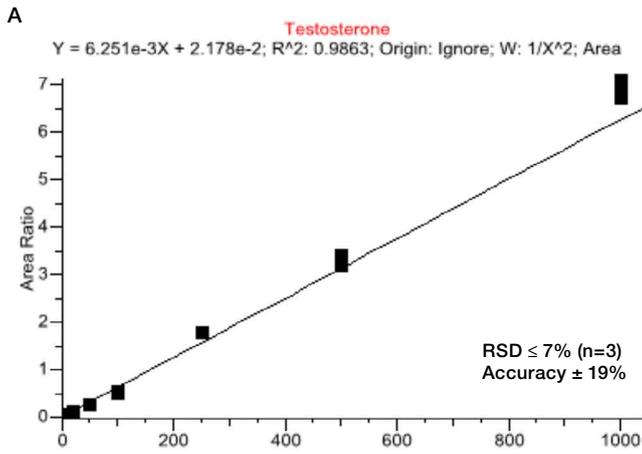
Data was acquired and processed using Thermo Scientific™ TraceFinder™ software.

## Results and discussion

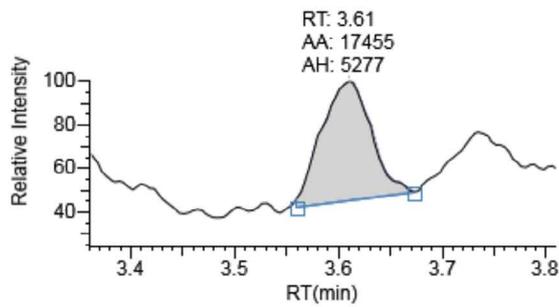
Lower limits of quantitation (LLOQ) for estradiol and testosterone were defined as the lowest concentration at which the back-calculated calibrator concentration on the linear calibration curve was within 20% of theoretical, the ion ratio was within 20% of the target, and replicate injections had a %RSD of less than 20%. The LLOQ of testosterone in plasma was 2 pg/mL (linearity range: 2–1000 pg/mL - Figure 1). For estradiol in plasma, the LLOQ was 20 pg/mL (linearity range: 20–10,000 pg/mL - Figure 2). The precisions were less than 8% and 7% for testosterone and estradiol, respectively, for all replicates at all concentrations.

Table 3. SRM properties for analysis of estradiol and testosterone.

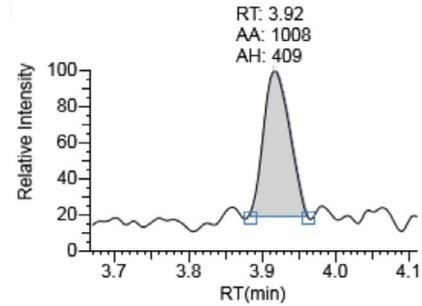
| Compound                       | Retention Time (min) | RT Window (min) | Polarity | Precursor (m/z) | Product (m/z) | Collision Energy (V) | RF Lens (V) | Quantifying or Confirming Ion |
|--------------------------------|----------------------|-----------------|----------|-----------------|---------------|----------------------|-------------|-------------------------------|
| Estradiol                      | 4.8                  | 1               | Negative | 271.1           | 183.1         | 40                   | 100         | Quantifying                   |
| Estradiol                      | 4.8                  | 1               | Negative | 271.1           | 145.1         | 40                   | 100         | Confirming                    |
| Estradiol-d5                   | 4.8                  | 1               | Negative | 276.1           | 187.1         | 41                   | 100         | Quantifying                   |
| Estradiol-d5                   | 4.8                  | 1               | Negative | 276.1           | 147.1         | 38                   | 100         | Confirming                    |
| Testosterone                   | 4.9                  | 1               | Positive | 289.1           | 109.1         | 25                   | 51          | Quantifying                   |
| Testosterone                   | 4.9                  | 1               | Positive | 289.1           | 97.1          | 22                   | 51          | Confirming                    |
| Testosterone- <sup>13</sup> C3 | 4.9                  | 1               | Positive | 292.3           | 112.1         | 25                   | 51          | Quantifying                   |
| Testosterone- <sup>13</sup> C3 | 4.9                  | 1               | Positive | 292.3           | 100.1         | 21                   | 51          | Confirming                    |



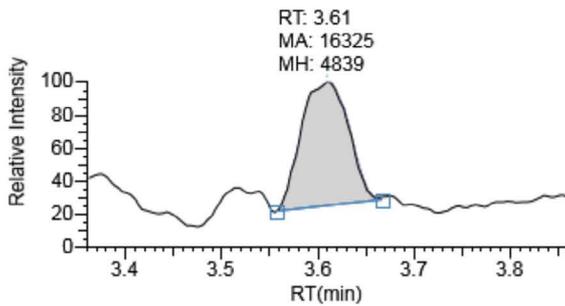
**B** Plasma\_Testosterone\_2pGmL\_01 Testosterone m/z: 109.100



**B** HESI\_E2\_Cal-06\_001 Estradiol m/z: 183.0710



**C** Plasma\_Testosterone\_2pGmL\_01 Testosterone m/z: 97.000



**C** HESI\_E2\_Cal-06\_001 Estradiol m/z: 145.0400

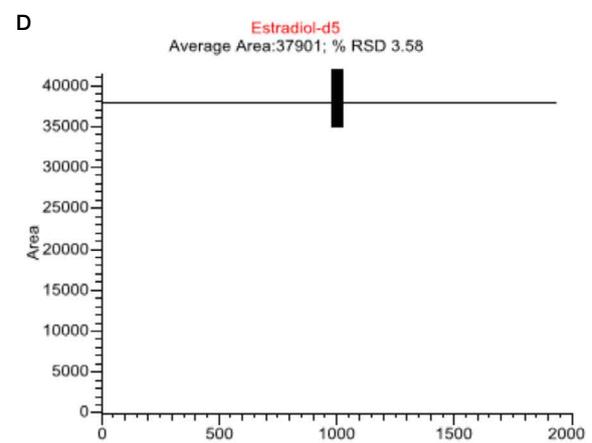
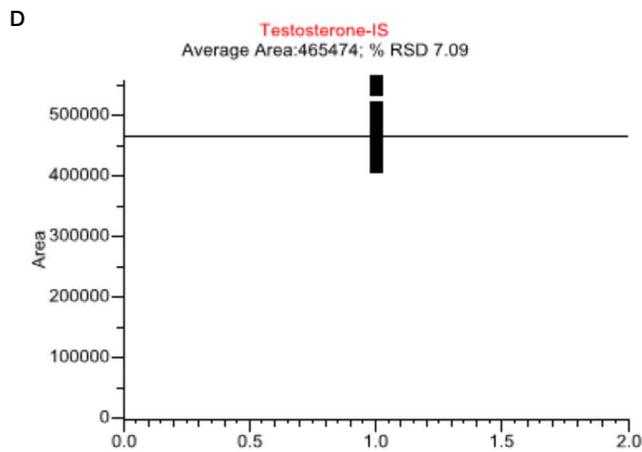
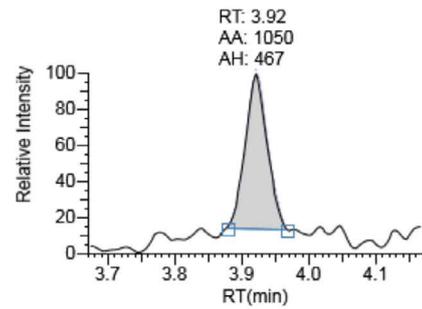


Figure 1. Testosterone data: (A) calibration curve, (B) LLOQ chromatogram for quantifying ion, (C) LLOQ chromatogram for confirming ion, and (D) internal standard.

Figure 2. Estradiol data: (A) calibration curve, (B) LLOQ chromatogram for quantifying ion, (C) LLOQ chromatogram for confirming ion, and (D) internal standard.

## Conclusion

- The TSQ Altis triple quadrupole mass spectrometer provides the superior sensitivity required for the analysis of estradiol and testosterone in plasma for clinical research.
- Limits of quantitation in plasma of 2 pg/mL for testosterone and 20 pg/mL for estradiol were obtained with the methodology described in this technical note.

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

Find out more at [thermofisher.com/Altis-Quantis](https://thermofisher.com/Altis-Quantis)

**ThermoFisher**  
S C I E N T I F I C