## Testosterone high sensitivity ELISA kit

A highly sensitive competitive immunoassay for the measurement of Testosterone in human serum and plasma samples.

A highly sensitive competitive immunoassay for the measurement of Testosterone in human serum and plasma samples. Results are read on a colorimetric plate reader at 405 nm.

Testosterone is one of the most important androgens secreted into the bloodstream and is synthesized from pregnenolone which is itself formed from cholesterol. Testosterone is the main androgen secreted by the Leydig cells of the testes and effects both primary and secondary sexual development such as muscle mass and sex drive. In adult humans, testosterone circulates in plasma predominately bound to proteins, including specific sex hormone binding globulin (SHBG) and nonspecific proteins such as albumin.

Citations: 27

View Online »

**Ordering Information** 

**Order Online** »

ADI-901-176

5x96 wells

Manuals, SDS & CofA

View Online »

- Excellent sensitivity
- Validated for use with human samples
- Stable color-coded reagents reduce errors
- Save with bulk kit package!

## **Handling & Storage**

Short Term Storage +4°C

Long Term Storage +4°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

**Application** Colorimetric detection, ELISA

**Application Notes**For the quantitative determination of human testosterone

in plasma, serum, and urine samples.

Assay Time 3 hours

Compatibility This product is compatible with the Absorbance 96 Plate

Reader.

**Contents** GxM Microtiter plate, Assay buffer, Standard, Antibody,

Conjugate, Wash buffer concentrate, pNpp Substrate,

Stop solution

**Sensitivity** 2.6 pg/ml (range 3.9 – 1000 pg/ml)

Species Reactivity Human

Wavelength 405 nm

Last modified: May 29, 2024

