Testosterone ELISA kit

Easy-to-use, highly sensitive ELISA kit for measuring testosterone with results in 3h

The Testosterone ELISA kit is a colorimetric competitive enzyme immunoassay kit with results in 3 hours. Absorbance is read at 405 nm. Ready-to-use liquid color-coded reagents are provided to reduce error. The reproducible results lot-after-lot provide excellent consistency, ideal for longitudinal studies.

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: 235

View Online »

Ordering Information

Order Online »

ADI-901-065

5x96 wells

Manuals, SDS & CofA

View Online »

- Sensitive measurement of testosterone, detecting as little as 5.67 pg/ml
- Simple and ready-to-use liquid color-coded reagents reduces error
- Steroid displacement reagent provided to assay free or total testosterone
- High throughput format with results in 3 hours
- Excellent consistency lot-after-lot ideal for longitudinal studies
- Fully quantitative results that surpass semi-quantitative
 Western blot analysis
- · Save with bulk kit package!

Handling & Storage

Long Term Storage +4°C

Blue Ice Shipping

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of testosterone in

> culture supernatants, plasma, serum, and saliva from any species. Not suitable for human serum and plasma. For those samples use product ADI-901-176. Cited sample

type includes tissue and feces.

Assay Time 3 hours

Compatibility This product is compatible with the **Absorbance 96 Plate**

Reader.

Contents GxM Microtiter plate, Conjugate, Antibody, Assay buffer 3,

Wash buffer concentrate, Standard, pNpp Substrate, Stop

solution, Steroid displacement reagent

Crossreactivity Testosterone (100%), 19-hydroxytestosterone (14.6%),

Androstendione (7.20%), Dehydroepiandrosterone

(0.72%), Estradiol (0.40%) and <0.001%:

Dihydrotestosterone, Estriol, Aldosterone, Corticosterone, Cortisol, Cortisone, Estrone, Progesterone, Pregenolone

Sensitivity 5.67 pg/ml (range 7.81 - 2,000 pg/ml)

Species Reactivity Species independent

Technical Info / Product Notes Cited samples:

Hormone ELISAs: Cited Samples

Wildlife Endocrinology: Cited Sample Types

Application Note:

Validation of Two Steroid Hormone ELISA Kits with Ovine

Blood Samples.

Wavelength 405 nm

Last modified: May 29, 2024

