

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: 234

[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

Shipping Blue Ice

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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)