

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: 30

[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



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)