

TGF- α ELISA kit

Highly sensitive ELISA kit enabling detection of TGF- α in 3.5 hours.

The TGF- α ELISA kit is a colorimetric immunometric immunoassay kit with results in 3.5 hours.

TGF- α is a ligand for the epidermal growth factor receptor, which activates a signaling pathway for cell proliferation, differentiation and development. TGF- α act as either a transmembrane-bound ligand or a soluble ligand. TGF- α is thought to be involved in oncogenesis and wound healing.

Ordering Information

[Order Online »](#)

ENZ-KIT132-0001	96 wells
-----------------	----------

Manuals, SDS & CofA

[View Online »](#)

- Sensitive measurement of TGF- α , detecting as little as 9.375 pg/ml
- Negligible reactivity from similar growth factors
- High throughput format with results in 3.5 hours
- Fully quantitative results that surpass semi-quantitative Western blot analysis

Handling & Storage

Short Term Storage +4°C

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Transforming growth factor α

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of TGF- α in serum, plasma, cell culture supernatant, tissue homogenates in human samples.

Assay Time 3.5 hrs

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

Contents Microtiter Plate, Standard, Detection Antibody, Detection Antibody Diluent, Sample Diluent, Wash Buffer Concentrate, Conjugate, Conjugate Diluent, Substrate, and Stop Solution.

Sensitivity 9.375 pg/ml

Species Reactivity Human

UniProt ID P01135

Wavelength 450 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-be@enzolifesciences.com)

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