

TNF- α (human), ELISA kit

Highly sensitive ELISA for measuring TNF- α , the primary mediator of immune regulation.

The TNF- α (human), EIA kit is a colorimetric immunometric enzyme immunoassay kit with results in 4 hours. Absorbance is read at 450 nm. Assay up to 39 samples in duplicate per 96-well plate. The broad dynamic range makes assaying different pathologies easy.

Tumor Necrosis Factor- α (TNF- α) is a 17.5kDa, 157 amino acid protein that is a potent lymphoid factor, which exerts cytotoxic effects on a wide range of tumor cells and other target cells. TNF- α has been suggested to play a pro-inflammatory role and has been detected in synovial fluid of patients with rheumatoid arthritis. It is the primary mediator of immune regulation. The biosynthesis of TNF- α is tightly controlled, being produced in extremely small quantities in quiescent cells, but is a major secreted factor in activated cells.

Citations: 17

[View Online »](#)

Ordering Information

[Order Online »](#)

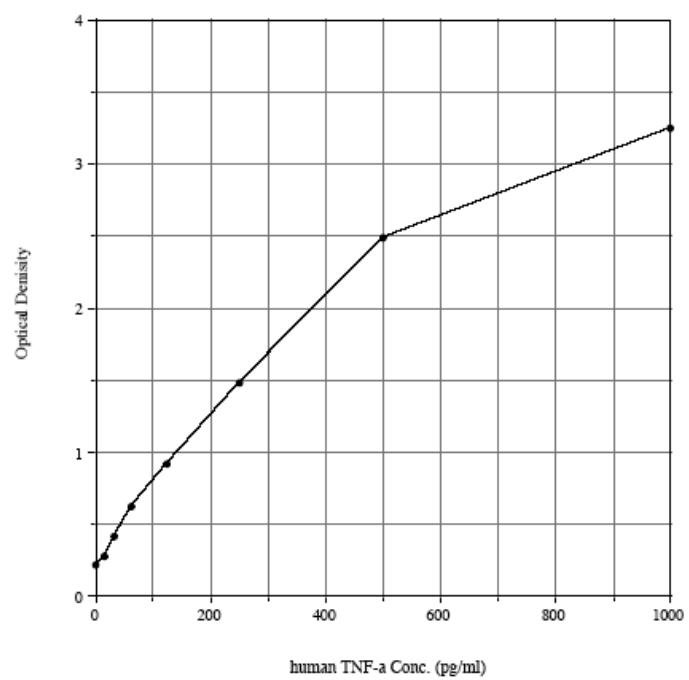
ADI-900-099

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Highly sensitive ELISA, measuring as little as 8.4 pg/ml
- High throughput format with results for up to 39 samples in duplicate in just 4 hours
- Stable liquid color coded reagents and pre-coated plates reduce errors
- Fully quantitative results that surpass semi-quantitative Western blot analysis



Handling & Storage

Use/Stability Store all components at +4°C, except standard at -20°C.

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Tumor necrosis factor- α , TNFSF 2

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of human TNF- α in culture supernatants, plasma, and serum.

Assay Time 4 hours

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

Contents Microtiter plate, Conjugate, Antibody, Assay buffer 13, Wash buffer concentrate, Standard, TMB Substrate, Stop solution 2

Sensitivity 8.43 pg/ml (range 15.63 – 1,000 pg/ml)

Species Reactivity Human

Technical Info / Product Notes **Cited samples:**
[Cytokine ELISAs Cited Samples](#)

UniProt ID P01375

Wavelength 450 nm



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)