

TNF- α (mouse), ELISA kit

The TNF- α (mouse), ELISA kit is a colorimetric immunometric enzyme immunoassay kit

The TNF- α (mouse), EIA kit is a colorimetric immunometric enzyme immunoassay kit with results in 5 hours. Absorbance is read at 450 nm.

Tumor Necrosis Factor- α (TNF- α) is a 17.5 kDa, 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: 43

[View Online »](#)

Ordering Information

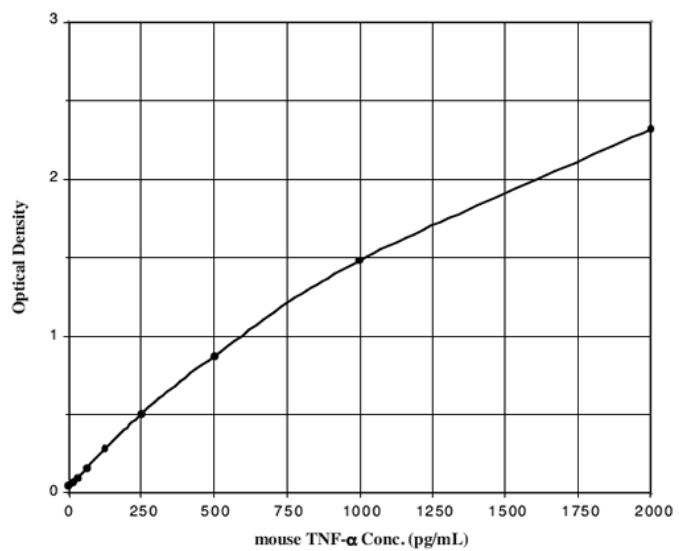
[Order Online »](#)

ADI-900-047	96 wells
-------------	----------

Manuals, SDS & CofA

[View Online »](#)

- Stable color coded reagents and pre-coated plates reduce errors
- Assay up to 38 samples in just 5 hours
- Save time, money, and precious sample compared to 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 mouse TNF- α in culture supernatants and serum. Cited sample type includes tissue (ref. 1) and plasma.

Assay Time 5 hours

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

Contents Microtiter plate, Conjugate, Antibody, Assay buffer 16, Wash buffer concentrate, Standard, TMB Substrate, Stop solution

Sensitivity 3.9 pg/ml (range 15.63 – 2,000 pg/ml)

Species Reactivity Mouse

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

UniProt ID P06804

Wavelength 450 nm

Last modified: November 5, 2025



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)