

# TNF- $\alpha$ (mouse), ELISA kit

The TNF- $\alpha$  (mouse), ELISA kit is a colorimetric immunometric enzyme immunoassay kit

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

Tumor Necrosis Factor- $\alpha$  (TNF- $\alpha$ ) 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- $\alpha$  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- $\alpha$  is tightly controlled, being produced in extremely small quantities in quiescent cells, but is a major secreted factor in activated cells.

Citations: 42

[View Online »](#)

## Ordering Information

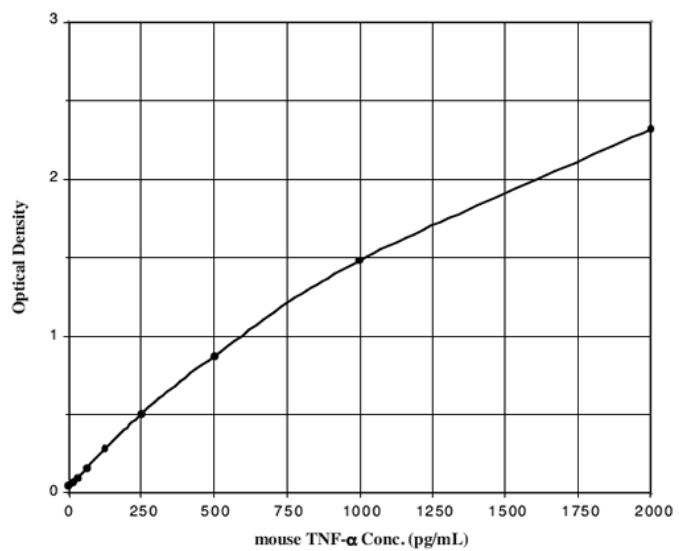
[Order Online »](#)

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** Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Tumor necrosis factor- $\alpha$ , TNFSF 2

**Application** Colorimetric detection, ELISA

**Application Notes** For the quantitative determination of mouse TNF- $\alpha$  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



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)