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: 42

View Online »

Ordering Information

Order Online »

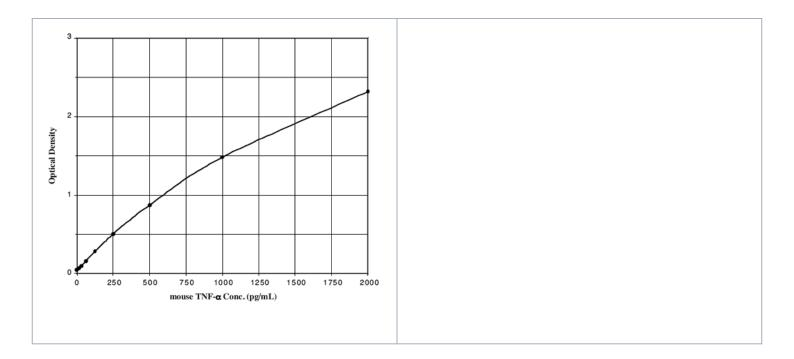
ADI-900-047

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 Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative NameTumor 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 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com