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. The larger package size offers a lower cost-per-test.

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: 17

View Online »

Ordering Information

Order Online »

ADI-901-099

5x96 wells

Manuals, SDS & CofA

View Online »

- Highly sensitive ELISA, measuring as liitle 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
- Save with bulk kit package!

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

concentrate, 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

Last modified: May 29, 2024



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