

TNF antibody (mAb)

Catalog Nos: 61601, 61602

RRID: AB_2793696

Isotype: IgG2a

Application(s): ELISA

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein A Chromatography

Host: Mouse

Concentration: 0.85 µg/µl

Molecular Weight: 17 kDa

Background: TNF (Tumor Necrosis Factor) is a cytokine that binds to TNFRSF1A/TNFR1 and TNFRSF1B/TNFR2. It is mainly secreted by macrophages and can induce cell death of certain tumor cell lines. It is potent pyrogen causing fever by direct action or by stimulation of interleukin-1 secretion and is implicated in the induction of cachexia. Under certain conditions it can stimulate cell proliferation and induce cell differentiation.

Immunogen: Human TNF.

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

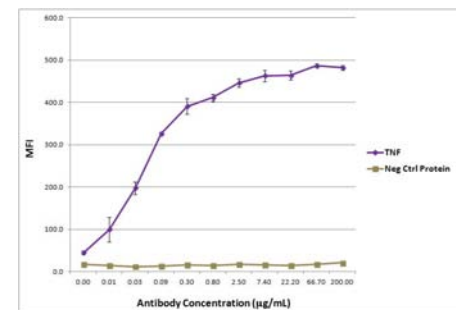
Application Notes:

Applications Validated by Active Motif:

ELISA: Individual optimization may be required.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



TNF antibody (mAb) tested by ELISA.

Recombinant TNF protein or a negative control protein was conjugated to Luminex® microspheres and then incubated in the same well with a serial dilution of TNF antibody (mAb). Biotinylated anti-mouse IgG followed by Streptavidin-PE was used to detect the bound TNF antibody. Mean Fluorescence Intensity (MFI) for each protein was measured in a Luminex® 200™.