

Purified Mouse Anti-human TNF α Antibody
MAB1, monoclonalCatalog number: V1032355
Unit size: 0.1 mg**Product Details**

| | |
|--------------------|---|
| Storage Conditions | 2-8°C with minimized light exposure. Do not freeze. |
| Expiration Date | 12 months upon receiving |
| Concentration | Lot specific (please consult certificate of analysis for given lot) |
| Formulation | Phosphate-buffered saline (PBS, pH 7.2), 15 mM sodium azide, 0.2% (w/v) BSA |

Antibody Properties

| | |
|--------------------|--------------|
| Species Reactivity | Human |
| Class | Primary |
| Clonality | Monoclonal |
| Host | Mouse |
| Immunogen | TNF α |
| Clone | MAB1 |

Biological Properties

| | |
|-------------|--|
| Preparation | Antibody purified by affinity chromatography and then conjugated with under optimal conditions |
| Application | WB, ELISA |

Applications

Tumor necrosis factor (also known as TNF- α , TNF-a, Tumor necrosis factor ligand superfamily member 2 or Cachectin) is a transmembrane protein with a molecular weight of 26 kDa, found in the phagocytic cup, integral component of plasma membrane and recycling endosome of cells. In humans, TNF- α has been thought to be involved with important functions like cytokine activity. It is the subject of extensive research because of the fact that it acts in the lipopolysaccharide-mediated signaling pathway, intrinsic apoptotic signaling pathway in response to DNA damage and extrinsic apoptotic signaling pathway via death domain receptors. Sequencing of TNF- α has exemplified it contains 2 types of conserved structural units: extracellular and cytoplasmic domain. TNF- α is a suppressor of fat cell differentiation, myosin-light-chain-phosphatase activity and interleukin-6 production while also is a positive regulator of interleukin-8 production, ERK1 and ERK2 cascade and protein catabolic process. It is an integral part of organismal processes, in particular, microglial cell activation, JNK cascade and cortical actin cytoskeleton organization. TNF- α binds to protease, transcription regulatory region sequence-specific DNA and identical protein. It has been found to be involved in establishment of endothelial barrier, endothelial cell apoptotic process and transcription by RNA polymerase II. TNF- α is clinically significant because abnormalities in its function have been closely linked to diseases such as psoriatic arthritis.