



6405 Mira Mesa Blvd., Suite 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

Human Tumor Necrosis Factor-alpha Recombinant

Catalog #: 90244-A
Lot#: 220721

Description: Recombinant human TNF-alpha is a disulfide-linked monomer protein, consisting of 158 amino acids and migrates as an approximately 17 kDa protein under reducing conditions.

Source: Optimized DNA sequence encoding human TNF- α mature chain was expressed in *E. coli*.

Formulation: Lyophilized from a 0.2 μ m filtered solution (1 mg/ml) containing phosphate buffered saline, pH 7.2.

Reconstitution: A quick mix of the vial followed by reconstitution in distilled water to concentration not less than 0.1 mg/ml. This solution can then be diluted into other buffers.

Storage: The lyophilized protein is stable for at least 2 years from date of receipt when stored at -20°C. Upon reconstitution, store in working aliquots at 2 - 8°C for up to one month, or at -20°C for up to six months, in the presence of a carrier protein. Avoid repeated freeze/thaw cycles.

Purity: >95%, as determined by SDS-PAGE and HPLC.

Endotoxin Level: Endotoxin level was found to be < 0.1 ng/ μ g (1EU/ μ g), using the LAL gel clot method.

Biological Activity: The ED50, determined by cytolysis of murine L929 cells in the presence of Actinomycin D is \leq 0.05 ng/ml, corresponding to a specific activity of > 1.9 x 10⁷ units/mg.

Protein Sequence:
MSTESMIRDVELAEEALPKKTGGPQG
SRRCLFSLFSFLIVAGATTLFCLLHFG
VIGPQREEFPRDLSLISPLAQAVRSSH
RTPSDKPVAVHVVANPQAEGQLQWLN
RRANALLANGVELRDNQLVVPSEGLY
LIYSQVLFKGQGPCSTHVLTTHTISRIA
VSYQTKVNLLSAIKSPCQRETPEGAEA
KPWYEPIYLGGVFQLEKGDRLSAEINR
PDYLDFAESGQVYFGIHAL

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.
To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: support@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com