

## TNF-beta

Recombinant mouse protein expressed in Sf9 cells

**Catalog # T861-40BG**

Lot # C1996-5

### Product Description

Mouse active form TNF-beta (34-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The protein accession number is [P09225](#).

### Gene Aliases

h1b382; LT; LT-alpha; LT-[a]; LTalpha; Ltx; LT[a]; LTA; Tnfb; TNFSF1; Tnfsf1b

### Endotoxin Level

Less than 0.1 ng/µg (IEU/µg) of TNF-beta.

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

### Storage and Stability

Stored at desiccated below -20°C. It is stable at 4°C between 2-7 days. For long term storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

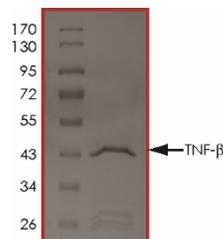
### Scientific Background

TNF-beta is a member of the tumor necrosis factor cytokine family that are primarily expressed on lymphocytes that control lipid metabolism. TNF-beta is highly inducible, secreted, and forms heterotrimers with lymphotoxin-beta which anchor lymphotoxin-alpha to the cell surface. TNF-beta acts as a novel therapeutic route for the treatment of dyslipidemia (1). The presence of TNF-beta accelerates the initiation of immune responses and protects the host from larger viral challenges (2).

### References

1. Lo, J. C. et.al: Lymphotoxin beta receptor-dependent control of lipid homeostasis. *Science* 316: 285-288, 2007.
2. Lund, F. E. et.al: Lymphotoxin-alpha-deficient mice make delayed, but effective, T and B cell responses to influenza. *J. Immun.* 169: 5236-5243, 2002.

### Purity



The purity was determined to be **>90%** by densitometry analysis in reducing SDS-PAGE Coomassie blue gel.

Approx. MW of the major band was **43 kDa**.

## TNF-beta

Recombinant mouse protein expressed in Sf9 cells

Catalog #

T861-40BG

Lot #

C1996-5

Purity

>90%

Concentration

0.1 µg/µl

Stability

2yr at -20°C from date of shipment

Storage & Shipping

Stored at desiccated below -20°C. It is stable at 4°C between 2-7 days. For long term storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**