

TNF-alpha

Recombinant human protein expressed in E. coli cells

Catalog # T860-40N

Lot # S346-3

Product Description

Recombinant human TNF-alpha was expressed in E. coli cells and purified by proprietary chromatographic techniques. The protein accession number is [P01375](#).

Gene Aliases

TNF, TNF α , DIF, TNFA, TNFSF2

Endotoxin Level

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

Formulation

Recombinant protein stored in 1 X PBS, 10% 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.

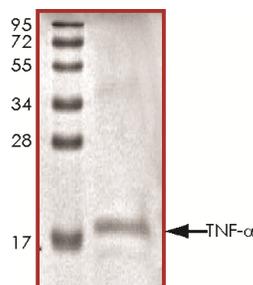
Scientific Background

TNF-alpha is a multifunctional proinflammatory cytokine that belongs to the tumor necrosis factor (TNF) superfamily and it is secreted predominantly by monocytes/macrophages that have effects on lipid metabolism, coagulation, insulin resistance, and endothelial function (1). TNF-alpha is involved in the regulation of a wide spectrum of biological processes including cell proliferation, differentiation, apoptosis, lipid metabolism, and coagulation and implicated in a variety of diseases. TNF-alpha also participates in the homeostatic activity-dependent regulation of synaptic connectivity and play roles in synaptic plasticity and modulating responses to neural injury (2).

References

1. Shirai, T. et al: Cloning and expression in Escherichia coli of the gene for human tumour necrosis factor. Nature 313: 803-806, 1985.
2. Stellwagen D. et al: Synaptic scaling mediated by glial TNF-alpha. Nature. 2006 Apr 20; 440(7087):1054-9.

Purity



The purity was determined to be **>95%** by densitometry analysis in reducing SDS-PAGE Coomassie blue gel.

Approx. MW of the major band was **18 kDa (17,484 Daltons)** by MW calculation).

TNF-alpha

Recombinant human protein expressed in E. coli cells

Catalog Number **T860-40N**

Specific Lot Number **S346-3**

Purity	>95%
Concentration	0.2 µg/µl
Stability	1 yr 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
www.signalchem.com

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