

Recombinant Human CNTF

Recombinant human protein expressed in E. coli cells

Catalog # C851-40N

Lot # G3211-17

Product Description

Recombinant Human CNTF was expressed in E. coli cells. The protein accession number is [P26441](#)

Gene Aliases

Ciliary neurotrophic factor

Endotoxin Level

<1.0 EU/µg of recombinant protein as determined by the LAL method

Formulation

Lyophilized from a 0.2 µm filtered PBS pH 7.5

Reconstitution Protocol

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

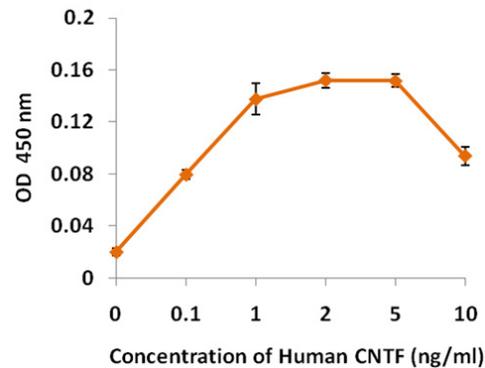
Storage and Stability

The lyophilized protein is stable for at least one year from date of receipt at -70°C. Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

Scientific Background

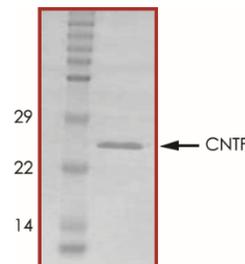
CNTF is found predominantly in peripheral nerve tissues. The main source appears to be myelin-associated Schwann cells in peripheral nerves and astrocytes in the central nervous system. CNTF is present at very high concentrations within intraocular tissues that contain the same muscle cells innervated by ciliary ganglionic neurons in vivo. CNTF receptors are expressed exclusively in the nervous system and skeletal muscle. The alpha subunit of the receptor is anchored to the cell membrane by a glycosyl-phosphatidylinositol linkage. The other receptor subunit, known as gp130, is also a component of receptors for LIF, Oncostatin M, IL-6 and IL-11, and seems to be involved also in signaling for CT-1. CNTF induces the expression of genes encoding hepatic acute phase proteins such as Haptoglobin, Alpha-1-antichymotrypsin, Alpha-2-Macroglobulin and Beta-fibrinogen in human hepatoma cells (HepG2) and in primary rat hepatocytes with a time course and dose-response comparable with that of IL-6.

Specific Activity



The ED50 as determined by the dose-dependent proliferation of human erythroleukemic cell line was found to be <2 ng/mL

Purity



Recombinant Human CNTF resolved on a 15% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **22.0 kDa**

Recombinant Human CNTF

Recombinant human protein expressed in E. coli cells

Catalog Number	C851-40N
Specific Lot Number	G3211-17
Purity	95%
Stability	6 mos. at -20°C from date of shipment
Storage & Shipping	The lyophilized protein is stable for at least one year from date of receipt at -70°C. Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

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.