

Recombinant Human TNFRSF10B

Recombinant human protein expressed in E. coli cells

Catalog # T881-40BN

Lot # G3223-19

Product Description

Recombinant Human TNFRSF10B was expressed in E. coli cells. The protein accession number is [O14763](#)

Gene Aliases

Tumor necrosis factor receptor superfamily member 10B, Death receptor 5, TRAIL receptor 2, TRAIL-R2, CD262, TNF-related apoptosis-inducing ligand receptor 2

Endotoxin Level

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

Formulation

Lyophilized from a solution in Tris and NaCl

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.

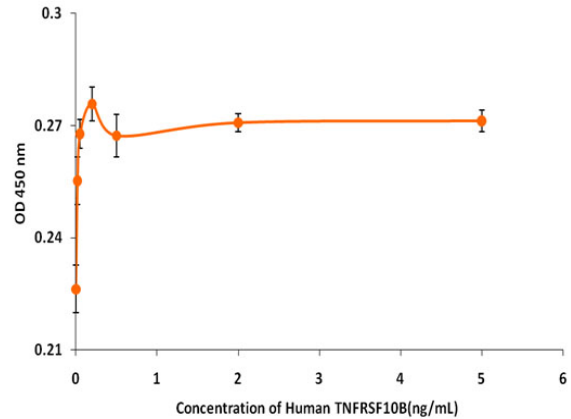
Storage and Stability

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

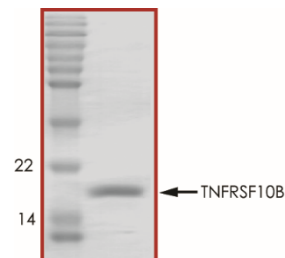
TNFRSF10B is a member of the TNFR superfamily and is a cell surface receptor that is involved in apoptosis. It contains an intracellular death domain and can be activated by binding to either membrane-anchored or soluble TRAIL/Apo2L. Diseases associated with TNFRSF10B include squamous cell carcinoma, head and neck and oral cavity cancer. Among its related pathways are Akt Signaling and Akt Signaling. It is closely related to TRAIL R1/DR4, sharing 55% amino acid sequence identity.

Specific Activity



The ED(50) was determined by the dose-dependent proliferation of human umbilical vein endothelial cells and was found to be <0.2ng/mL.

Purity



Recombinant Human TNFRSF10B resolved on a 15% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **15.0 kDa**

Recombinant Human TNFRSF10B

Recombinant human protein expressed in E. coli cells

Catalog Number	T881-40BN
Specific Lot Number	G3223-19
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.