

Recombinant Human TWEAK Receptor (TNFRSF12A)

Recombinant human protein expressed in E. coli cells

Catalog # T883-40N

Lot # G3223-23

Product Description

Recombinant Human TWEAK Receptor (TNFRSF12A) was expressed in E. coli cells. The protein accession number is [Q9NP84](#)

Gene Aliases

Fibroblast growth factor-inducible immediate-early response protein 14, Tweak-receptor, CD-antigen: CD266

Endotoxin Level

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

Formulation

Lyophilized from a 0.2 µm filtered solution in PBS

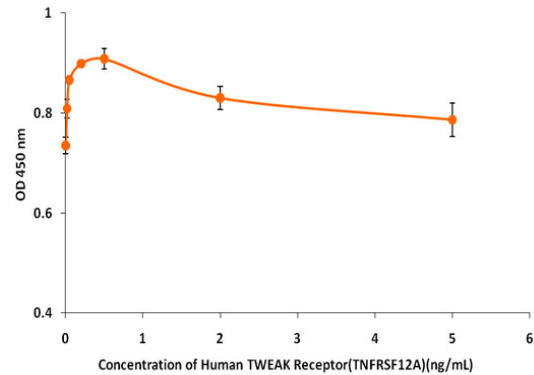
Storage and Stability

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

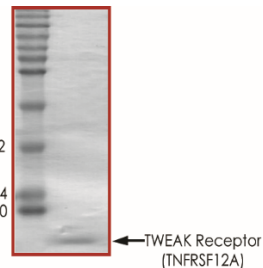
Tumor necrosis factor receptor superfamily member 12A (TNFRSF12A) is the receptor for TNFSD12/TWEAK and is the smallest member of the TNF receptor superfamily. It is composed of a signal peptide, an extracellular domain, a membrane-anchoring domain, and a cytoplasmic "death" domain, which can activate a cell's apoptotic mechanisms. It also promotes endothelial cell proliferation and angiogenesis, and modulates cellular adhesion to matrix proteins. TNFRSF12A expression can be found in the spleen, thymus, small intestine, and colon, and can be activated by membrane-anchored or soluble versions of TWEAK.

Specific Activity



The ED50 is 2.0-12.0 µg/mL as determined by its ability to inhibit TWEAK-dependent proliferation in human umbilical vein endothelial cells (with use of 1 µg/mL rhTWEAK)

Purity



Recombinant Human TWEAK Receptor (TNFRSF12A) resolved on a 10% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **5.6 kDa**

Recombinant Human TWEAK Receptor (TNFRSF12A)

Recombinant human protein expressed in E. coli cells

Catalog Number	T883-40N
Specific Lot Number	G3223-23
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.