TL1A, His-Tag, Avi-Tag Recombinant

Product Information

Description:	Recombinant human TL1A (TNF-like ligand 1A (TL1A, also known as Vascular endothelial growth inhibitor, VEGI or TNFSF15), encompassing amino acids 72 to 251(end). This construct contains an N-terminal His-tag (6xHis) followed by an Avi-Tag [™] . This protein was affinity purified. The protein runs as two bands in SDS-PAGE due to variable glycosylation, as confirmed by PNGase treatment.
Background:	TNF-like ligand 1A (TL1A, also known as Vascular endothelial growth inhibitor, VEGI or TNFSF15) is an anti-angiogenic cytokine. It is an important mediator of inflammation, participates in innate and adaptive immune homeostasis through binding to its receptor, DR3, activating downstream signaling. Numerous studies showed that soluble TL1A can be detected in the serum of patients with T-cell mediated autoimmune diseases like rheumatoid arthritis, psoriatic arthritis, and inflammatory bowel disease. In addition, recent clinical studies suggested that anti-TL1A antibody treatment is a promising therapeutic approach in inflammatory and auto-immune disorder.
Species:	Human
Construct:	TL1A (His-Avi-72-251(end))
Concentration:	4.32 mg/ml
Expression System:	HEK293
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
MW:	23 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Genbank Accession:	NM_005118
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

