

Human Thymic stromal lymphopoietin Recombinant

Catalog: 90250-B
Lot: 240409

Product Information

Description:	Recombinant human TSLP is a protein consisting of Tyr29-Gln159 amino acid residue subunits, due to glycosylation migrates as an approximately 25 kDa protein under reducing conditions in SDS-PAGE. DNA sequence encoding Human TSLP mature chain was expressed in HEK293 cells.
Species:	Human
Expression System:	HEK293
Purity:	>95% by SDS-PAGE and HPLC
Formulated In:	Lyophilized from a 0.2 µm filtered 1 mg/ml solution of 20 mM phosphate buffer, pH 7.4, 150 mM NaCl.
Endotoxin Level:	<0.1 ng/µg (1 EU/µg), using the LAL gel clot method.
MW:	Due to glycosylation migrates as an approximately 25 kDa protein on SDS-PAGE
Storage:	The lyophilized protein is stable for at least years from date of receipt at -20° C. Upon reconstitution, this cytokine can be stored in working aliquots at +4°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.
Reconstitution:	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.
Biological Activity	The activity as determined by binding Human TSLPR in an ELISA assay.