

# Human Thymic stromal lymphopoietin Recombinant

Catalog: 90250-A  
Lot: 240409

## Product Information

<b>Description:</b>	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.
<b>Species:</b>	Human
<b>Expression System:</b>	HEK293
<b>Purity:</b>	>95% by SDS-PAGE and HPLC
<b>Formulated In:</b>	Lyophilized from a 0.2 µm filtered 1 mg/ml solution of 20 mM phosphate buffer, pH 7.4, 150 mM NaCl.
<b>Endotoxin Level:</b>	<0.1 ng/µg (1 EU/µg), using the LAL gel clot method.
<b>MW:</b>	Due to glycosylation migrates as an approximately 25 kDa protein on SDS-PAGE
<b>Storage:</b>	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.
<b>Reconstitution:</b>	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.
<b>Biological Activity</b>	The activity as determined by binding Human TSLPR in an ELISA assay.