

### Instruction Manual

<b>Catalog Number</b>	PK-AB718-4025
<b>Synonyms</b>	TSLP Antibody: mTSLP, Thymic stromal lymphopoietin
<b>Description</b>	Thymic stromal lymphopoietin (TSLP) has recently been identified as an important factor capable of driving dendritic cell maturation and activation. TSLP is a four-helix-bundle cytokine that is expressed mainly by barrier epithelial cells and is a potent activator of several cell types such as myeloid dendritic cells. TSLP is involved in the positive selection of regulatory T cells, maintenance of peripheral CD4+ T cell homeostasis and the induction of CD4+ T cell-mediated allergic reaction. TSLP is also capable of supporting the growth of fetal liver and adult B cell progenitors and their differentiation to the IgM-positive stage of B cell development. Amino acid sequence analysis has shown poor homology between human and mouse TSLP although they exhibit similar biological functions and are expressed in similar tissues. Despite its predicted molecular weight, TSLP often migrates at a higher molecular weight in SDS-PAGE. At least two differentially spliced isoforms of TSLP are known to exist.
<b>Quantity</b>	100 µg
<b>Source / Host</b>	Rabbit
<b>Immunogen</b>	TSLP antibody was raised in rabbits against a 20 amino acid peptide from near the center of mouse TSLP.
<b>Purification Method</b>	Affinity chromatography purified via peptide column.
<b>Clone / IgG Subtype</b>	Polyclonal antibody
<b>Species Reactivity</b>	Mouse, Rat
<b>Specificity</b>	
<b>Formulation</b>	Antibody is supplied in PBS containing 0.02% sodium azide.
<b>Reconstitution</b>	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 µl or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
<b>Storage &amp; Stability</b>	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
<b>Applications</b>	E, WB, IHC, IF    INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually.  TSLP antibody can be used for detection of TSLP by Western blot at 0.5 - 2 µg/mL. Despite its predicted molecular weight, TSLP often migrates at a higher molecular weight in SDS-PAGE. Antibody can also be used for immunohistochemistry starting at 2.5 µg/mL. For immunofluorescence start at 20 µg/mL.
<b>Images</b>	Available upon request.
<b>References</b>	Ziegler SF and Liu Y-J. Thymic stromal lymphopoietin in normal and pathogenic T cell development and function. <i>Nature Immunol.</i> 2006; 7:709-14.  Sims JE, Williams DE, Morrissey PJ, et al. Molecular cloning and biological characterization of a novel murine lymphoid growth factor. <i>J. Exp. Med.</i> 2000; 192:671-80.  Levin SD, Koelling RM, Friend SL, et al. Thymic stromal lymphopoietin: a cytokine that promotes the development of IgM+ cells in vitro and signals via a novel mechanism. <i>J. Immunol.</i> 1999; 162:677-83.  Quentmeier H, Drexler HG, Fleckenstein D, et al. Cloning of human thymic stromal lymphopoietin (TSLP) and signaling mechanisms leading to proliferation. <i>Leukemia</i> 2001; 15:1286-92.
<b>Images</b>	Available upon request.
<b>Related Products</b>	Cat.No. PK-AB718-4025P; TSLP Peptide  Cat.No. PK-AB718-1288; A20 Cell Lysate
FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.	