

# TSLP (CT) antibody (pAb)

## Rabbit Anti-Human TSLP (CT)

### Instruction Manual

Catalog Number	PK-AB718-4021
Synonyms	TSLP Antibody: Thymic stromal lymphopietin
Description	Thymic stromal lymphopietin (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.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	TSLP antibody was raised in rabbits against a 19 amino acid peptide from near the carboxy terminus of human TSLP.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human
Specificity	
Formulation	Antibody is supplied in PBS containing 0.02% sodium azide.
Reconstitution	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.
Storage & Stability	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.
Applications	E, WB, IHC Note: 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 1 - 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.
Images	Available upon request.
References	Ziegler SF and Liu Y-J. Thymic stromal lymphopietin 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 lymphopietin: 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 lymphopietin (TSLP) and signaling mechanisms leading to proliferation. <i>Leukemia</i> 2001; 15:1286-92.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4021P; TSLP Peptide Cat.No. PK-AB718-1205; Jurkat Cell Lysate

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.