

# TSLP Receptor antibody (pAb)

## Rabbit Anti-Human TSLP Receptor

### Instruction Manual

Catalog Number	PK-AB718-4207
Synonyms	TSLP Receptor Antibody: TSLP Receptor, Thymic stromal lymphopoietin receptor
Description	Thymic stromal lymphopoietin (TSLP) has recently been identified as an important factor capable of driving dendritic cell maturation and activation. It 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. Its receptor TSLP-R will bind TSLP in a low-affinity fashion in transfected cells; co-transfection with IL-7R-alpha cDNA results in high-affinity binding and a functional heteromeric complex. This heteromeric receptor requires stat5 for TSLP-mediated signal transduction and is inhibited by SOCS-1. Despite its predicted molecular weight, TSLP-R often migrates at a higher molecular weight in SDS-PAGE.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	TSLP Receptor antibody was raised in rabbits against a 19 amino acid peptide from near the center of human TSLP receptor.
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, IF INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually.  TSLP-R antibody can be used for detection of TSLP-R by Western blot at 0.5 - 1 µg/mL. Despite its predicted molecular weight, TSLP-R often migrates at a higher molecular weight in SDS-PAGE. Antibody can also be used for immunohistochemistry starting at 5 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	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. Park LS, Martin U, Garka K, et al. Cloning of the murine thymic stromal lymphopoietin (TSLP) receptor: Formation of a functional heteromeric complex requires interleukin 7 receptor. <i>J. Exp. Med.</i> 2000; 192:659-70.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4207P; TSLP Receptor Peptide Cat.No. PK-AB718-1304; Human Liver Tissue Lysate

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