

SLPI antibody (pAb)

Rabbit Anti-Human/Mouse SLPI

Instruction Manual

Catalog Number	PK-AB718-4249
Synonyms	SLPI Antibody; Secretory leukocyte protease inhibitor, HUSI
Description	Secretory leukocyte protease inhibitor (SLPI) is produced at mucosal surfaces, primarily the upper respiratory tract and is thought to play an important role in the antiprotease defense mechanism of the lung. SLPI forms inhibitory complexes with numerous proteolytic enzymes such as neutrophil elastase, and has been shown to possess anti-inflammatory, anti-viral, and antibacterial activities. Its expression in oral epithelial cells is stimulated by HIV-1 gp120, suggesting that SLPI is a component of the oral mucosal response to HIV-1. In peripheral blood monocytes, SLPI can inhibit NF- κ B activation by inhibiting I κ B degradation in the cytoplasm and competing for NF- κ B binding sites in the nucleus. This attenuation of the inflammatory response may also act to suppress liver metastases and other cancer cell invasions, but promote blood-borne metastasis via an invasion-independent pathway.
Quantity	100 μ g
Source / Host	Rabbit
Immunogen	SLPI antibody was raised in rabbits against a 17 amino acid peptide from near the center of human SLPI.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse
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 μ l 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 INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. SLPI antibody can be used for detection of SLPI by Western blot at 1 - 2 μ g/mL. Antibody can also be used for immunohistochemistry starting at 10 μ g/mL.
Images	Available upon request.
References	Abe T, Kobayashi N, Yoshimura K, et al. Expression of the secretory leukoprotease inhibitor gene in epithelial cells. <i>J. Clin. Invest.</i> 1991; 87:2207-15. Hiemstra PS, Fernie-King BA, McMichael J, et al. Antimicrobial peptides: mediators of innate immunity as templates for the development of novel anti-infective and immune therapies. <i>Curr. Pharm. Des.</i> 2004; 10:2891-905. Jana NK, Gray LR and Shugars DC. Human immunodeficiency virus type 1 stimulates the expression and production of secretory leukocyte protease inhibitor (SPLI) in oral epithelial cells: a role for SLPI in innate mucosal immunity. <i>J. Virol.</i> 2005; 79:6432-40 Taggart CC, Cryan S-A, Weldon S, et al. Secretory leucoprotease inhibitor binds to NF- κ B binding sites in monocytes and inhibits p65 binding. <i>J. Exp. Med.</i> 2005; 202:1659-68.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4249P; SLPI Peptide Cat.No. PK-AB718-1224; Daudi Cell Lysate

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