

sRANK Ligand antibody (pAb)

Rabbit Anti-Human/Mouse/Rat sRANK Ligand (TNFSF11, TRANCE, OPGL)

Instruction Manual

Catalog Number	PK-AB718-3963
Synonyms	sRANK Ligand Antibody: sRANK-L, TNFSF11, TRANCE, OPGL
Description	The receptor activator of NF- κ B ligand (RANK-L) is a recently discovered member of the TNF-ligand family involved in the regulation of the T cell-dependent immune response, lymph node organogenesis and bone formation. RANK-L exists as both a normal, transmembrane form and a truncated, soluble form (sRANK-L), both of which can stimulate the receptor. Activation of T cells, such as by treatment with interleukin-7, induces RANK-L production and leads to an increase of osteoclast formation and bone loss. Finally, sRANK-L can activate the antiapoptotic kinase Akt through a signaling complex involving Src kinase and TRAF6, suggesting sRANK-L may also play a role in regulating apoptosis.
Quantity	100 μ g
Source / Host	Rabbit
Immunogen	sRANK-L antibody was raised against a 14 amino acid peptide from near the center of human sRANK-L.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
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 Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. sRANK-L antibody can be used for detection of sRANK-L by Western blot at 0.25 - 0.5 μ g/mL. Antibody can also be used for immunohistochemistry starting at 5 μ g/mL.
Images	Available upon request.
References	Wong BR, Rho J, Arron J, et al. TRANCE is a novel ligand of the tumor necrosis factor receptor family that activates c-Jun N-terminal kinase in T cells. <i>J. Biol. Chem.</i> 1997; 272:25190-4. Kong YY, Yoshida H, Sarosi I, et al. OPGL is a key regulator of osteoclastogenesis, lymphocyte development and lymph-node organogenesis. <i>Nature</i> 1999; 397:315-23. Weitzmann MN, Cenci S, Rifas L, et al. Interleukin-7 stimulates osteoclast formation by up-regulating the T-cell production of soluble osteoclastogenic cytokines. <i>Blood</i> 2000; 96:1873-8. Bharti AC, Takada Y, Shishodia S, et al. Evidence that receptor activator of the nuclear factor (NF)-kappaB ligand can suppress cell proliferation and induce apoptosis through activation of a NF-kappaB-independent and TRAF6-dependent mechanism. <i>J. Biol. Chem.</i> 2004; 279:6065-76.
Images	Available upon request.
Related Products	sRANK-L Peptide, Cat. No. PK-AB718-3963P Liver Lysate, Cat. No. PK-AB718-1304 Liver Lysate, Cat. No. PK-AB718-1464

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