

Instruction Manual

Catalog Number	PK-AB718-4313
Synonyms	STEAP4 Antibody: Six transmembrane epithelial antigen of prostate 4, STAMP2
Description	The six-transmembrane epithelial antigen of prostate 4 (STEAP4) is a member of a family of metalloreductases identified as cell-surface antigens in prostate tissue. Similar to two other members of the STEAP family (STEAP 2 and STEAP3), STEAP4 promotes both iron and copper reduction. STEAP4 is highly expressed in placenta, lung, heart and prostate tissues, and overexpressed in prostate cancer cells compared to normal prostate cells. Overexpression of STEAP4 in prostate cells significantly increases cell growth and colony formation, suggesting STEAP4 may have a role in cell proliferation and prostate cancer progression. At least two isoforms are known to exist for this protein. This STEAP4 antibody does not cross-react with other STEAP proteins.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	STEAP4 antibody was raised in rabbits against a 14 amino acid peptide from near the center of human STEAP4.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	At least two isoforms are known to exist for this protein. This STEAP4 antibody does not cross-react with other STEAP proteins.
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 Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. STEAP4 antibody can be used for detection of STEAP4 by Western blot at 0.5 - 1 µg/mL.
Images	Available upon request.
References	Korkmaz CG, Korkmaz KS, Kurys P, et al. Molecular cloning and characterization of STAMP2, an androgen-regulated six transmembrane protein that is overexpressed in prostate cancer. <i>Oncogene</i> 2005; 24:4934-45. Ohgami RS, Campagna DR, McDonald A, et al. The Steap proteins are metalloreductases. <i>Blood</i> 2006; 108:1388-94.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4313P; STEAP4 Peptide Cat.No. PK-AB718-1464 - Rat Liver Tissue Lysate

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