

# STEAP2 (CT) antibody (pAb)

Rabbit Anti-Human/Mouse STEAP2 (CT)

## Instruction Manual

Catalog Number	PK-AB718-4307
Synonyms	STEAP2 Antibody: Six transmembrane epithelial antigen of prostate 2
Description	The six-transmembrane epithelial antigen of prostate 2 (STEAP2) is a member of a family of metalloredutases identified as cell-surface antigens in prostate tissue. Similar to two other members of the STEAP family (STEAP 3 and STEAP4), STEAP2 promotes both iron and copper reduction. STEAP2 expression in transfected cells also correlated with iron or copper uptake, suggesting that the STEAP family of proteins may function to stimulate iron and copper uptake. STEAP2 is widely expressed in many tissues in the plasma membrane, but is most highly expressed in prostate. At least three isoforms of STEAP2 are known to exist. This STEAP2 antibody does not cross-react with other STEAP proteins.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	STEAP2 antibody was raised in rabbits against a 16 amino acid peptide from near the carboxy terminus of human STEAP2.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse
Specificity	This STEAP2 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, IHC, IF    INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. STEAP2 antibody can be used for detection of STEAP2 by Western blot at 1 - 2 µg/mL. Antibody can also be used for immunohistochemistry starting at 2.5 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Porkka KP, Helenius MA and Visakorpi T. Cloning and characterization of a novel six-transmembrane protein STEAP2, expressed in normal and malignant prostate. Lab Invest. 2002; 82:1573-82. Ohgami RS, Campagna DR, McDonald A, et al. The Steap proteins are metalloredutases. Blood 2006; 108:1388-94. Ohgami RS, Campagna DR, Greer EL, et al. Identification of a ferrireductase required for efficient transferrin-dependent iron uptake in erythroid cells. Nat. Genet. 2005; 37:1264-9.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4307P; STEAP2 Peptide Cat.No. PK-AB718-1312; Human Prostate Tissue Lysate

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