

## STEAP4 Antibody [#C11207]

<b>Catalog Number:</b>	C11207
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q687X5
<b>Other Names:</b>	EC 1.16.1.-; Six-transmembrane epithelial antigen of prostate 4; SixTransMembrane protein of prostate 2; Tumor necrosis factor, alpha-induced protein 9
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human STEAP4.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	STEAP4 antibody detects endogenous levels of total STEAP4 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	Korkmaz C.G., Oncogene 24:4934-4945(2005). Ota T., Nat. Genet. 36:40-45(2004). Ohgami R.S., Nat. Genet. 37:1264-1269(2005).