

STEA2 Antibody [#C11375]

Catalog Number:	C11375
Concentration:	1mg/ml
Swiss-Prot No.:	Q8NFT2
Other Names:	EC 1.16.1.-; Six-transmembrane epithelial antigen of prostate 2; SixTransMembrane protein of prostate 1; Prostate cancer-associated protein 1
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg ²⁺ and Ca ²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized peptide derived from C-terminal of human STEA2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	STEA2 antibody detects endogenous levels of total STEA2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 92%, Positives = 100%)
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
References:	Korkmaz K.S., J. Biol. Chem. 277:36689-36696(2002). Porkka K.P., Lab. Invest. 82:1573-1582(2002). Ohgami R.S., Blood 108:1388-1394(2006).