

SERC1 Antibody [#C11754]

Catalog Number:	C11754
Concentration:	1mg/ml
Swiss-Prot No.:	Q9NRX5
Other Names:	Serine incorporator 1, Tumor differentially expressed protein 2, Tumor differentially expressed 1 protein-like, SERINC1, KIAA1253, TDE1L, TDE2, UNQ396/PRO732
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 internal of human SERC1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	SERC1 antibody detects endogenous levels of total SERC1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Nagase T., DNA Res. 6:337-345(1999). Hu R.-M., Proc. Natl. Acad. Sci. U.S.A. 97:9543-9548(2000). Clark H.F., Genome Res. 13:2265-2270(2003).