

SGOL1 Antibody [#C11176]

Catalog Number:	C11176
Concentration:	1mg/ml
Swiss-Prot No.:	Q5FBB7
Other Names:	Shugoshin-like 1, hSgo1, Serologically defined breast cancer antigen NY-BR-85, SGO1
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 SGOL1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	SGOL1 antibody detects endogenous levels of total SGOL1 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	McGuinness B.E., PLoS Biol. 3:433-449(2005). Kitajima T.S., Curr. Biol. 15:353-359(2005). Tang Z., Proc. Natl. Acad. Sci. U.S.A. 101:18012-18017(2004).