

SGOL2 Antibody [#C18741]

Catalog Number:	C18741
Concentration:	1mg/ml
Swiss-Prot No.:	Q562F6
Other Names:	Shugoshin-like 2; Tripin; SGOL2
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 SGOL2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	SGOL2 antibody detects endogenous levels of total SGOL2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:10000
References:	Oliveira V., Submitted (APR-2002) to the EMBL/GenBank/DDBJ databases. Ota T., Nat. Genet. 36:40-45(2004). Hillier L.W., Nature 434:724-731(2005).