

Cyclosome 1 Antibody [#C12817]

Catalog Number:	C12817
Concentration:	1mg/ml
Swiss-Prot No.:	Q9H1A4
Other Names:	ANAPC1, ANC1, Anaphase promoting complex subunit 1, Cyclosome subunit 1, Mitotic checkpoint regulator, Protein Tsg24, TSG24
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 Cyclosome 1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Cyclosome 1 antibody detects endogenous levels of total Cyclosome 1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:10000
References:	Joergensen P.M., Gene 262:51-59(2001). Ota T., Nat. Genet. 36:40-45(2004). Kraft C., EMBO J. 22:6598-6609(2003).