

p57 Kip2 (Ab-310) Antibody [#B0967]

Catalog Number:	B0967
Concentration:	1mg/ml
Swiss-Prot No.:	P49918
Other Names:	CDKN1C, CDN1C, CDNC, Cyclin-dependent kinase inhibitor 1C, Cyclin-dependent kinase inhibitor p57, KIP2
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 non-phosphopeptide derived from human p57 Kip2 around the phosphorylation site of threonine 310 (E-Q-T ^P -P-R).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	p57 Kip2 (Ab-310) antibody detects endogenous levels of total p57 Kip2 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:5000
References:	Matsuoka S., Genes Dev. 9:650-662(1995). Reid L.H., Cancer Res. 56:1214-1218(1996). Tokino T., Hum. Genet. 97:625-631(1996).