

## p57KIP2 Antibody [#C10659]

<b>Catalog Number:</b>	C10659
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P49918
<b>Other Names:</b>	CDKN1C, CDN1C, CDNC, Cyclin-dependent kinase inhibitor 1C, Cyclin-dependent kinase inhibitor p57, KIP2
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human p57KIP2.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	p57KIP2 antibody detects endogenous levels of total p57KIP2 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100 ELISA: 1:10000
<b>References:</b>	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).