

## CDK2 Antibody [#C10474]

<b>Catalog Number:</b>	C10474
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P24941
<b>Other Names:</b>	Cell division protein kinase 2, EC 2.7.11.22, kinase Cdk2, p33 protein kinase
<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 CDK2.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	CDK2 antibody detects endogenous levels of total CDK2 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Elledge S.J., EMBO J. 10:2653-2659(1991). Gray N.S., Science 281:533-538(1998). Brown N.R., Cell Cycle 6:1350-1359(2007).