



**Signosis**  
BioSignal Capture

---

## CDK2(Ab-160) Antibody

Catalog Number 21111

(For Research Use Only)

---

- Catalog Number:** 21111-1, 21111-2, 21111  
**Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl  
**Form of Antibody:** 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.  
**Storage/Stability:** Store at -20°C/1 year  
**Immunogen:** The antiserum was produced against synthesized non-phosphopeptide derived from human CDK2 around the phosphorylation site of Threonine160.  
**Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.  
**Specificity/Sensitivity:** CDK2(Ab-160) Antibody detects endogenous levels of total CDK2 protein.  
**Reactivity:** Human, Mouse, Rat  
**Applications:** WB: 1:500~1:1000      ELISA: 1:10000  
**Swiss-Prot No. :** P24941  
**References:** Ukomadu C, et al.(2003) J Biol Chem; 278(7): 4840-6.  
Morris MC, et al.(2002)J Biol Chem; 277(26): 23847-53.  
Brown NR, et al.(1999)J Biol Chem; 274(13): 8746-56.  
Liu Y, et al.(2004) J Biol Chem; 279(6): 4507-14.