

## Chk1 (Ab-317) Antibody [#B7041]

- Catalog Number:** B7041-1, B7041-2, B7041
- 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 Chk1 around the phosphorylation site of serine 317.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity:** Chk1 (Ab-317) antibody detects endogenous levels of Chk1 protein around serine 317.
- Reactivity:** Human, Rat
- Applications:** WB: 1:500~1:1000    IHC:1:50~1:100  
ELISA: 1:10000
- Swiss-Prot No. :** O14757
- References:** Zhang YW, et al. (2005) Mol Cell; 19(5): 607-18.  
Bhoumik A, et al. (2005) Mol Cell; 18(5): 577-87.  
Rocha S, et al. (2005) EMBO J.  
Clarke CA, et al. (2005) Biochem J.  
Yu X, et al. (2004) Mol Cell Biol; 24(21): 9478-86.