



Chk1(Phospho-Ser317) Antibody

Catalog Number 11141

(For Research Use Only)

- Catalog Number:** 11141-1, 11141-2, 11141
Amount: 50µg/50µl, 100µg/100µl, 200µg/200µl
Form of Antibody: Rabbit IgG in phosphate buffered saline (without Mg²⁺ and Ca²⁺), 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 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 phosphopeptide. The antibody against non-phosphopeptide was removed by chromatography using non-phosphopeptide corresponding to the phosphorylation site.
Specificity/Sensitivity: Chk1(Phospho-Ser317) Antibody detects endogenous levels of Chk1 only when phosphorylated at Serine 317.
Reactivity: Human, Rat
Applications: WB: 1:500~1:1000 ELISA: 1:4000
Swiss-Prot No. : O14757
References: Zhang YW, et al. (2005) Mol Cell; 19(5): 607-18.
Bhounik 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.

