



Chk2(Ab-68) Antibody

Catalog Number 21087

(For Research Use Only)

- Catalog Number:** 21087-1, 21087-2, 21087
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 non-phosphopeptide derived from human Chk2 around the phosphorylation site of Threonine 68.
Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity/Sensitivity: Chk2(Ab-68) Antibody detects endogenous levels of total Chk2 protein.
Reactivity: Human, Mouse, Rat
Applications: WB: 1:500~1:1000 ELISA: 1:10000
Swiss-Prot No. : O96017
References: Bhoumik A, et al. (2005) Mol Cell; 18(5): 577-87.
Gorgoulis VG, et al. (2005) Nature; 434(7035): 907-13.
Falck J, et al. (2005) Nature; 434(7033): 605-11.
Jin ZH, et al. (2005) Oncogene; 24(12): 1973-81.
Li J, et al. (2005) J Biol Chem; 280(12): 12041-50.

