



Chk1(Ab-280) Antibody

Catalog Number 21113

(For Research Use Only)

- Catalog Number:** 21113-1, 21113-2, 21113
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: This antiserum was produced against synthesized non-phosphopeptide derived from human Chk1 around the phosphorylation site of Serine 280.
Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity/Sensitivity: Chk1(Ab-280) Antibody detects endogenous levels of total Chk1 protein.
Reactivity: Human
Applications: WB: 1:500~1:1000 ELISA: 1:10000
Swiss-Prot No. : O14757
References: Conn CW, et al. (2004)Dev Cell; 7(2): 275-81.
King FW, et al. (2004)Cell Cycle; 3(5): 634-7.
Shtivelman E, et al. (2002) Curr Biol; 12(11): 919-24.

