



Chk2 (Phospho-Ser516) Antibody

Catalog Number 11147

(For Research Use Only)

- Catalog Number:** 11147-1, 11147-2, 11147
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 Chk2 around the phosphorylation site of serine 516 (Q-P-S^P-T-S).
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: Chk2 (phospho-Ser516) antibody detects endogenous levels of Chk2 only when phosphorylated at serine 516.
Reactivity: Human
Applications: WB: 1:500~1:1000 ELISA: 1:10000
Swiss-Prot No. : O96017
References: Wu X, Chen J et al.(2003)J Biol Chem; 278(38): 36163-8.
Schwarz JK, et al.(2003)Mol Cancer Res; 1(8): 598-609.
Lou Z, et al.(2003) Nature; 421(6926): 957-61.



