

Chk1(Phospho-Ser301)Antibody [#8A0862]

Catalog Number:	8A0862
Concentration:	1mg/ml
Origin of Antibody:	Rabbit
Swiss-Prot No.:	O14757
Other Names:	CHEK1, Chk1, EC 2.7.11.1, kinase Chk1
All Sites:	Human: Ser
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg^{2+} and Ca^{2+}), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized phosphopeptide derived from human Chk1 around the phosphorylation site of serine 301 (D-F-S ^P -P-V).
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:	Chk1 (Phospho-Ser301) antibody detects endogenous levels of Chk1 only when phosphorylated at serine 301.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IF: 1:100~1:500 ELISA: 1:40000
References:	Sanchez Y., Science 277:1497-1501(1997). Shieh S.-Y., Genes Dev. 14:289-300(2000). Uensal-Kacmaz K., Mol. Cell. Biol. 25:3109-3116(2005). Matsuoka S., Science 316:1160-1166(2007).