

## Casein Kinase I $\alpha$ (Ab-321) Antibody [#B1194]

<b>Catalog Number:</b>	B1194
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P48729
<b>Other Names:</b>	CK1-alpha, CKI-alpha, CSNK1A1, Casein kinase I, alpha isoform, EC 2.7.11.1, KC1A, kinase
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized non-phosphopeptide derived from human Casein Kinase I $\alpha$ around the phosphorylation site of threonine 321 (A-Q-T <sup>P</sup> -P-T).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	Casein Kinase I $\alpha$ (Ab-321) antibody detects endogenous levels of total Casein Kinase I $\alpha$ protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:5000
<b>References:</b>	Tapia C., FEBS Lett. 349:307-312(1994). Fish K.J., J. Biol. Chem. 270:14875-14883(1995). Liu C., I 108:837-847(2002).