

## CHKB Antibody [#C11923]

<b>Catalog Number:</b>	C11923
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q9Y259
<b>Other Names:</b>	Choline/ethanolamine kinase, Choline kinase beta (EC 2.7.1.32)(CK), Ethanolamine kinase (EC 2.7.1.82)(EK), CHKB, CHETK, CHKL
<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 peptide derived from internal of human CHKB.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	CHKB antibody detects endogenous levels of total CHKB protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Yamazaki N., J. Biol. Chem. 275:31739-31746(2000). Collins J.E., Genome Biol. 5:RESEARCH84.1-RESEARCH84.11(2004). Yu L.-R., J. Proteome Res. 6:4150-4162(2007).