

## GLCTK Antibody [#C11307]

<b>Catalog Number:</b>	C11307
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q8IVS8
<b>Other Names:</b>	Glycerate kinase, EC 2.7.1.31, HBeAg-binding protein 4, GLYCTK, HBEBP4, LP5910
<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 GLCTK.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	GLCTK antibody detects endogenous levels of total GLCTK protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Guo J.-H., DNA Seq. 17:1-7(2006). Wan D., Proc. Natl. Acad. Sci. U.S.A. 101:15724-15729(2004). Ota T., Nat. Genet. 36:40-45(2004).