

KPB1/2 Antibody [#C10626]

Catalog Number:	C10626
Concentration:	1mg/ml
Swiss-Prot No.:	P46020
Other Names:	KPB1, PHKA, Phosphorylase B kinase alpha regulatory chain, skeletal muscle isoform, Phosphorylase kinase alpha M subunit, kinase PHKA1
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 peptide derived from internal of human KPB1/2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	KPB1/2 antibody detects endogenous levels of total KPB1/2 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:20000
References:	Wuellrich A., J. Biol. Chem. 268:23208-23214(1993). The MGC Project Team, Genome Res. 14:2121-2127(2004). Burwinkel B., Eur. J. Hum. Genet. 11:516-526(2003).