

ALPK2 Antibody [#C11273]

Catalog Number:	C11273
Concentration:	1mg/ml
Swiss-Prot No.:	Q86TB3
Other Names:	Alpha-protein kinase 2, EC 2.7.11.-, Heart alpha-protein kinase, ALPK2, HAK
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 ALPK2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ALPK2 antibody detects endogenous levels of total ALPK2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:40000
References:	Ota T., Nat. Genet. 36:40-45(2004). The German cDNA consortium, Submitted (JUN-2003) to the EMBL/GenBank/DDBJ databases. Ryazanov A.G., Curr. Biol. 9:R43-R45(1999).