

HCK Antibody [#B0938]

Catalog Number:	B0938
Concentration:	1mg/ml
Swiss-Prot No.:	P08631
Other Names:	B-cell/myeloid kinase, BMK, EC 2.7.10.2, Hemopoietic cell kinase, P56-HCK and P60-HCK, Tyrosine-protein kinase HCK, kinase Hck, p59-HCK/p60-HCK
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 HCK.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	HCK antibody detects endogenous levels of total HCK protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
References:	Ota T., Nat. Genet. 36:40-45(2004). Deloukas P., Nature 414:865-871(2001). Lock P., Mol. Cell. Biol. 11:4363-4370(1991).