

KITH Antibody [#C10270]

Catalog Number:	C10270
Concentration:	1mg/ml
Swiss-Prot No.:	P04183
Other Names:	Thymidine kinase, EC 2.7.1.21, TK
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 C-terminal of human KITH.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	KITH antibody detects endogenous levels of total KITH protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Bradshaw H.D. Jr., Mol. Cell. Biol. 4:2316-2320(1984). Flemington E., Gene 52:267-277(1987). Kreidberg J.A., Mol. Cell. Biol. 6:2903-2909(1986).