

FO XK1 Antibody [#C10811]

Catalog Number:	C10811
Concentration:	1mg/ml
Swiss-Prot No.:	P85037
Other Names:	FHX, FOXJ2 forkhead factor, fork head homologous X
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 FOXK1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	FOXK1 antibody detects endogenous levels of total FOXK1 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Ota T., Nat. Genet. 36:40-45(2004). Hillier L.W., Nature 424:157-164(2003). Kato M., Int. J. Mol. Med. 14:127-132(2004).