

KCIP1 Antibody [#C11788]

Catalog Number:	C11788
Concentration:	1mg/ml
Swiss-Prot No.:	Q9NZI2
Other Names:	Kv channel-interacting protein 1, KCHIP1, A-type potassium channel modulatory protein 1, Potassium channel-interacting protein 1, Vesicle APC-binding protein, KCNIP1, KCHIP1, VABP
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 N-terminal of human KCIP1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	KCIP1 antibody detects endogenous levels of total KCIP1 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: 1:20000
References:	An W.F., ure 403:553-556(2000). Van Hoorick D., Mol. Cell. Neurosci. 24:357-366(2003). Nakamura T.Y., FEBS Lett. 499:205-209(2001).