

KANK2 Antibody [#C14496]

Catalog Number:	C14496
Concentration:	1mg/ml
Swiss-Prot No.:	Q63ZY3
Other Names:	Ankyrin repeat domain-containing protein 25; SRC-1-interacting protein; SRC1-interacting protein; Matrix-remodeling-associated protein 3
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 KANK2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	KANK2 antibody detects endogenous levels of total KANK2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:40000
References:	Nagase T., DNA Res. 7:143-150(2000). Ota T., Nat. Genet. 36:40-45(2004). Olsen J.V., Cell 127:635-648(2006).