

MIPT3 Antibody [#C13085]

Catalog Number:	C13085
Concentration:	1mg/ml
Swiss-Prot No.:	Q8TDR0
Other Names:	DKFZp434F124, MIP-T3, MIPT3, TNF receptor-associated factor 3 interacting protein 1
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 MIPT3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MIPT3 antibody detects endogenous levels of total MIPT3 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
References:	Ling L., J. Biol. Chem. 275:23852-23860(2000). Niu Y., FEBS Lett. 550:139-143(2003). Hillier L.W., Nature 434:724-731(2005).