

MPRIP Antibody [#C16762]

Catalog Number:	C16762
Concentration:	1mg/ml
Swiss-Prot No.:	Q6WCQ1
Other Names:	Myosin phosphatase Rho-interacting protein, Rho-interacting protein 3; M-RIP; RIP3; p116Rip
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 MPRIP.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MPRIP antibody detects endogenous levels of total MPRIP protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:40000
References:	Surks H.K., J. Biol. Chem. 278:51484-51493(2003). Koga Y., J. Biol. Chem. 280:4983-4991(2005). Inazawa J., Submitted (DEC-2002) to the EMBL/GenBank/DDBJ databases.