

PPP1R1C Antibody [#C18013]

Catalog Number:	C18013
Concentration:	1mg/ml
Swiss-Prot No.:	Q8WVI7
Other Names:	Protein phosphatase 1 regulatory subunit 1C
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 PPP1R1C.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PPP1R1C antibody detects endogenous levels of total PPP1R1C protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:10000
References:	The MGC Project Team; Genome Res. 14:2121-2127(2004). Li N., Submitted (MAR-2002) to the EMBL/GenBank/DDBJ databases. Bechtel S., BMC Genomics 8:399-399(2007).