

PPP1R3C Antibody [#C18015]

Catalog Number:	C18015
Concentration:	1mg/ml
Swiss-Prot No.:	Q9UQK1
Other Names:	PPP1R3C; PPP1R5; PPR3C; protein phosphatase 1 regulatory subunit 3C; protein phosphatase 1 regulatory subunit 5; protein targeting to glycogen; PTG
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 PPP1R3C.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PPP1R3C antibody detects endogenous levels of total PPP1R3C protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Doherty M.J., FEBS Lett. 399:339-343(1996). Permana P.A., Biochem. Biophys. Res. Commun. 258:184-186(1999). Deloukas P., Nature 429:375-381(2004).