

## PPP1R7 Antibody [#C18018]

<b>Catalog Number:</b>	C18018
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q15435
<b>Other Names:</b>	PP1R7; Protein phosphatase 1, regulatory (Inhibitor) subunit 7; Protein phosphatase 1, regulatory subunit 7; Protein phosphatase-1 regulatory subunit 7; Sds22; Yeast SDS22 homolog
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human PPP1R7.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	PPP1R7 antibody detects endogenous levels of total PPP1R7 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%)
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:10000
<b>References:</b>	Renouf S., FEBS Lett. 375:75-78(1995). Ceulemans H., Eur. J. Biochem. 262:36-42(1999). Kalnine N., Submitted (OCT-2004) to the EMBL/GenBank/DDBJ databases.