

## PPP2R1B Antibody [#C18649]

<b>Catalog Number:</b>	C18649
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P30154
<b>Other Names:</b>	2AAB; beta isoform of regulatory subunit A, protein phosphatase 2; PP2A, A, PR65-beta isoform; PP2A, A, R1-beta isoform; PR65B; protein phosphatase 2 (formerly 2A), regulatory A (PR 65), beta isoform; protein phosphatase 2 (formerly 2A), regulatory A, beta isoform; protein phosphatase 2, structural/regulatory A, beta; serine/threonine protein phosphatase 2A, 65 kDa regulatory A, beta isoform; serine/threonine-protein phosphatase 2A 65 kDa regulatory A beta isoform
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), 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 PPP2R1B.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	PPP2R1B antibody detects endogenous levels of total PPP2R1B protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:40000
<b>References:</b>	Baysal B.E., Gene 217:107-116(1998). Wang S.S., Science 282:284-287(1998). Baysal B.E., Eur. J. Hum. Genet. 9:121-129(2001).