

## RGS7 Antibody [#C18326]

<b>Catalog Number:</b>	C18326
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P49802
<b>Other Names:</b>	RGS7; Regulator of G-protein signaling 7
<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 RGS7.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	RGS7 antibody detects endogenous levels of total RGS7 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 ELISA: 1:20000
<b>References:</b>	Kim E., Proc. Natl. Acad. Sci. U.S.A. 96:6371-6376(1999). Koelle M.R., Cell 84:115-125(1996). Puhl H.L. III, Submitted (MAR-2002) to the EMBL/GenBank/DDBJ databases.