

## GRK7 Antibody [#C11400]

<b>Catalog Number:</b>	C11400
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q8WTQ7
<b>Other Names:</b>	G protein-coupled receptor kinase 7 [Precursor], EC 2.7.11.16, G protein-coupled receptor kinase GRK7, GRK7, GPRK7
<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 GRK7.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	GRK7 antibody detects endogenous levels of total GRK7 protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:5000
<b>References:</b>	Chen C.-K., Mol. Vis. 7:305-313(2001). Weiss E.R., J. Neurosci. 21:9175-9184(2001).