

## CACNG7 Antibody [#C19525]

<b>Catalog Number:</b>	C19525
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P62955
<b>Other Names:</b>	CCG7; Neuronal voltage-gated calcium channel gamma-7 subunit; Voltage-dependent calcium channel gamma-7 subunit
<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 CACNG7.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	CACNG7 antibody detects endogenous levels of total CACNG7 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>	Burgess D.L., Genomics 71:339-350(2001). Chu P.-J., Gene 280:37-48(2001). Moss F.J., EMBO J. 21:1514-1523(2002).