

## ADCY1 Antibody [#C12031]

<b>Catalog Number:</b>	C12031
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q08828
<b>Other Names:</b>	ATP pyrophosphate-lyase, CYA1, Ca/calmodulin activated adenylyl cyclase, EC 4.6.1.1, adenylate cyclase 1, adenylate cyclase 1 (brain), adenylate cyclase, type I
<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 ADCY1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	ADCY1 antibody detects endogenous levels of total ADCY1 protein.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	WB: 1:500~1:1000   ELISA: 1:40000
<b>References:</b>	Hillier L.W., Nature 424:157-164(2003). Villacres E.C., Genomics 16:473-478(1993).