

## ADCY4 Antibody [#C12034]

<b>Catalog Number:</b>	C12034
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q8NFM4
<b>Other Names:</b>	Adenylate cyclase, type IV; Adenylyl cyclase 4; ATP pyrophosphate-lyase 4; EC 4.6.1.1
<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 ADCY4.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	ADCY4 antibody detects endogenous levels of total ADCY4 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000   ELISA: 1:5000
<b>References:</b>	Ludwig M.G., J. Recept. Signal Transduct. Res. 22:79-110(2002). The MGC Project Team; Genome Res. 14:2121-2127(2004). Ota T., Nat. Genet. 36:40-45(2004).