

## ADCY5/6 Antibody [#C12036]

<b>Catalog Number:</b>	C12036
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	O43306/O95622
<b>Other Names:</b>	ATP pyrophosphate-lyase, CYA6, Ca(2)-inhibitable adenylyl cyclase, EC 4.6.1.1, KIAA0422, adenylate cyclase, type VI
<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 ADCY5/6.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	ADCY5/6 antibody detects endogenous levels of total ADCY5/6 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Wicker R., Biochim. Biophys. Acta 1493:279-283(2000). Ishikawa K., DNA Res. 4:307-313(1997). Nakajima D., DNA Res. 9:99-106(2002).