

ADCY9 Antibody [#C20506]

Catalog Number:	C20506
Concentration:	1mg/ml
Swiss-Prot No.:	O60503
Other Names:	adenylate cyclase 9; EC 4.6.1.1; mKIAA0520
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg ²⁺ and Ca ²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized peptide derived from internal of human ADCY9.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ADCY9 antibody detects endogenous levels of total ADCY9 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:10000
References:	Hacker B.M., Genomics 50:97-104(1998). Small K.M., Pharmacogenetics 13:535-541(2003). Tantisira K.G., Hum. Mol. Genet. 14:1671-1677(2005).