

KAD6 Antibody [#C11948]

Catalog Number:	C11948
Concentration:	1mg/ml
Swiss-Prot No.:	Q9Y3D8
Other Names:	Adenylate kinase isoenzyme 6, EC 2.7.4.3, ATP-AMP transphosphorylase 6, AK6, AD-004, CGI-137
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 KAD6.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	KAD6 antibody detects endogenous levels of total KAD6 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Lai C.-H., Genome Res. 10:703-713(2000). Hu R.-M., Proc. Natl. Acad. Sci. U.S.A. 97:9543-9548(2000). Ren H., Proc. Natl. Acad. Sci. U.S.A. 102:303-308(2005).