

AKAP10 Antibody [#C10110]

Catalog Number:	C10110
Concentration:	1mg/ml
Swiss-Prot No.:	O43572
Other Names:	A kinase anchor protein 10, mitochondrial [Precursor], Protein kinase A-anchoring protein 10, PRKA10, Dual specificity A kinase-anchoring protein 2, D-AKAP-2, AKAP10
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 AKAP10.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	AKAP10 antibody detects endogenous levels of total AKAP10 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Wang L., Proc. Natl. Acad. Sci. U.S.A. 98:3220-3225(2001). Chatterjee T.K., Submitted (DEC-1997) to the EMBL/GenBank/DDBJ databases. The MGC Project Team, Genome Res. 14:2121-2127(2004).