

AKAP1 Antibody [#C11427]

Catalog Number:	C11427
Concentration:	1mg/ml
Swiss-Prot No.:	Q92667
Other Names:	A kinase anchor protein 1, mitochondrial precursor, A-kinase anchor protein 149 kDa, AKAP 149, AKAP149, AKP1, D-AKAP-1, Dual specificity A-kinase anchoring protein 1, PRKA1, Protein kinase A anchoring protein 1, S-AKAP84, Spermatid A-kinase anchor protein 84
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 AKAP1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	AKAP1 antibody detects endogenous levels of total AKAP1 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000
References:	Trendelenburg G., Biochem. Biophys. Res. Commun. 225:313-319(1996). Lin R.-Y., J. Biol. Chem. 270:27804-27811(1995). Hillman R.T., Genome Biol. 5:RESEARCH008.1-RESEARCH008.16(2004).