

AKAP14 Antibody [#C11279]

Catalog Number:	C11279
Concentration:	1mg/ml
Swiss-Prot No.:	Q86UN6
Other Names:	A-kinase anchor protein 14, A-kinase anchor protein 14, A-kinase anchor protein 28 kDa, AKAP14, AKAP28
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 interinal of human AKAP14.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	AKAP14 antibody detects endogenous levels of total AKAP14 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:5000
References:	Kultgen P.L., Mol. Biol. Cell 13:4156-4166(2002). The MGC Project Team, Genome Res. 14:2121-2127(2004).