

BAIAP2L1 Antibody [#C14787]

Catalog Number:	C14787
Concentration:	1mg/ml
Swiss-Prot No.:	Q9UHR4
Other Names:	BAIAP2L1; BI2L1
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 BAIAP2L1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	BAIAP2L1 antibody detects endogenous levels of total BAIAP2L1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%) Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Hu R.-M., Proc. Natl. Acad. Sci. U.S.A. 97:9543-9548(2000). Hillier L.W., Nature 424:157-164(2003). Scherer S.W., Science 300:767-772(2003).