

AIBP Antibody [#C14522]

Catalog Number:	C14522
Concentration:	1mg/ml
Swiss-Prot No.:	Q8NCW5
Other Names:	AIBP; apoA-I binding; APOA1BP; apolipoprotein A-I binding
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 AIBP.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	AIBP antibody detects endogenous levels of total AIBP protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:40000
References:	Ritter M., Genomics 79:693-702(2002). Gregory S.G., Nature 441:315-321(2006). The MGC Project Team; Genome Res. 14:2121-2127(2004).