

ABHD8 Antibody [#C14220]

Catalog Number:	C14220
Concentration:	1mg/ml
Swiss-Prot No.:	Q96I13
Other Names:	abhydrolase domain-containing 8; EC 3.-.-.-
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 ABHD8.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ABHD8 antibody detects endogenous levels of total ABHD8 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Ota T., Nat. Genet. 36:40-45(2004). The MGC Project Team; Genome Res. 14:2121-2127(2004).