

ABHD12 Antibody [#C14209]

Catalog Number:	C14209
Concentration:	1mg/ml
Swiss-Prot No.:	Q8N2K0
Other Names:	ABD12; ABHD12; abhydrolase domain containing 12
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 ABHD12.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ABHD12 antibody detects endogenous levels of total ABHD12 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Ota T., Nat. Genet. 36:40-45(2004). Deloukas P., Nature 414:865-871(2001). Mural R.J., Submitted (JUL-2005) to the EMBL/GenBank/DDBJ databases.