

ZC3H7B Antibody [#C11837]

Catalog Number:	C11837
Concentration:	1mg/ml
Swiss-Prot No.:	Q9UGR2
Other Names:	KIAA1031, ROXN, Rotavirus 'X' associated non-structural protein, Z3H7B, ZC3H7B, zinc finger CCCH-type containing 7B
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 ZC3H7B.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ZC3H7B antibody detects endogenous levels of total ZC3H7B protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Ota T., Nat. Genet. 36:40-45(2004). Rush J., Nat. Biotechnol. 23:94-101(2005). Kikuno R., DNA Res. 6:197-205(1999).