

ZNF435 Antibody [#C11695]

Catalog Number:	C11695
Concentration:	1mg/ml
Swiss-Prot No.:	Q9H4T2
Other Names:	Zinc finger and SCAN domain-containing protein 16, Zinc finger protein 435, Zinc finger protein 392, ZSCAN16, ZNF392, ZNF435
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 ZNF435.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ZNF435 antibody detects endogenous levels of total ZNF435 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Ota T., Nat. Genet. 36:40-45(2004). Mungall A.J., Nature 425:805-811(2003). The MGC Project Team, Genome Res. 14:2121-2127(2004).