

ZNF95 Antibody [#C11936]

Catalog Number:	C11936
Concentration:	1mg/ml
Swiss-Prot No.:	Q9Y2L8
Other Names:	Zinc finger protein with KRAB and SCAN domains 5, Zinc finger protein 95 homolog, Zfp-95, ZKSCAN5, KIAA1015, ZFP95
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 ZNF95.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ZNF95 antibody detects endogenous levels of total ZNF95 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IF: 1:500~1:1000 ELISA: 1:20000
References:	Dreyer S.D., Genomics 62:119-122(1999). Nagase T., DNA Res. 6:63-70(1999).