

ZNF541Antibody [#C11671]

Catalog Number:	C11671
Concentration:	1mg/ml
Swiss-Prot No.:	Q9H0D2
Other Names:	Zinc finger protein 541, ZNF541, ZN541
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 ZNF541.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ZNF541 antibody detects endogenous levels of total ZNF541 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Wiemann S., Genome Res. 11:422-435(2001). Yu L.-R., J. Proteome Res. 6:4150-4162(2007). Sjoeblom T., Science 314:268-274(2006).