

## ZNF436 Antibody [#C11660]

<b>Catalog Number:</b>	C11660
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q9C0F3
<b>Other Names:</b>	Zinc finger protein 436, ZNF436, KIAA1710
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human ZNF436.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	ZNF436 antibody detects endogenous levels of total ZNF436 protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Nagase T., DNA Res. 7:347-355(2000). Gregory S.G., Nature 441:315-321(2006). Sjoebloom T., Science 314:268-274(2006).