

## ZNF187 Antibody [#C11138]

<b>Catalog Number:</b>	C11138
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q16670
<b>Other Names:</b>	SRE-ZBP; zinc finger and SCAN domain-containing 26; zinc finger protein 187; ZN187; ZSCAN26
<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 ZNF187.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	ZNF187 antibody detects endogenous levels of total ZNF187 protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:10000
<b>References:</b>	Mungall A.J., Nature 425:805-811(2003). The MGC Project Team, Genome Res. 14:2121-2127(2004). Attar R.M., Mol. Cell. Biol. 12:2432-2443(1992).