

## ZNRF2 Antibody [#C15596]

<b>Catalog Number:</b>	C15596
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q8NHG8
<b>Other Names:</b>	E3 ubiquitin-protein ligase ZNRF2, EC 6.3.2.-; Zinc/RING finger protein 2; RING finger protein 202; Protein Ells2
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), 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 ZNRF2.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	ZNRF2 antibody detects endogenous levels of total ZNRF2 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 91%, Positives = 91%)
<b>Applications:</b>	IF: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	Araki T., J. Neurosci. 23:9385-9394(2003). Guo J.H., Submitted (JUL-2002) to the EMBL/GenBank/DDBJ databases. Olsen J.V., Cell 127:635-648(2006).