

## RNF138 Antibody [#C15572]

<b>Catalog Number:</b>	C15572
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q8WVD3
<b>Other Names:</b>	E3 ubiquitin-protein ligase RNF138; EC 6.3.2.-; RING finger protein 138; Nemo-like kinase-associated RING finger protein; NLK-associated RING finger protein
<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 RNF138.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	RNF138 antibody detects endogenous levels of total RNF138 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 92%, Positives = 100%); Rat (Identities = 92%, Positives = 100%)
<b>Applications:</b>	WB: 1:500~1:1000   ELISA: 1:40000
<b>References:</b>	Wang L., Submitted (APR-2000) to the EMBL/GenBank/DDBJ databases. Ota T., Nat. Genet. 36:40-45(2004). Kalline N., Submitted (MAY-2003) to the EMBL/GenBank/DDBJ databases.