

## MORF4L1 Antibody [#C11822]

<b>Catalog Number:</b>	C11822
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q9UBU8
<b>Other Names:</b>	Mortality factor 4-like protein 1, MORF-related gene 15 protein, MORF-related gene 15 protein, MSL3-1 protein, MORF4L1, MRG15, FWP006, HSPC008, HSPC061, PP368
<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 MORF4L1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	MORF4L1 antibody detects endogenous levels of total MORF4L1 protein.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Wan D., Proc. Natl. Acad. Sci. U.S.A. 101:15724-15729(2004). Zhang Q.-H., Genome Res. 10:1546-1560(2000). Cai Y., J. Biol. Chem. 278:42733-42736(2003).