

## UBAP2L Antibody [#C19423]

<b>Catalog Number:</b>	C19423
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q14157
<b>Other Names:</b>	UBAP2L; ubiquitin-associated protein 2-like; UBP2L
<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 UBAP2L.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	UBAP2L antibody detects endogenous levels of total UBAP2L protein.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Nagase T., DNA Res. 2:167-174(1995). Gregory S.G., Nature 441:315-321(2006). The MGC Project Team; Genome Res. 14:2121-2127(2004).