

## UBA5 Antibody [#C19417]

<b>Catalog Number:</b>	C19417
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q9GZZ9
<b>Other Names:</b>	ThiFP1; UBA5; ubiquitin-activating enzyme 5; UFM1-activating enzyme
<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 UBA5.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	UBA5 antibody detects endogenous levels of total UBA5 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:5000
<b>References:</b>	Komatsu M., EMBO J. 23:1977-1986(2004). Dou T., Submitted (OCT-2003) to the EMBL/GenBank/DDBJ databases. Wiemann S., Genome Res. 11:422-435(2001).