

## USP48 Antibody [#C19392]

<b>Catalog Number:</b>	C19392
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q86UV5
<b>Other Names:</b>	Ubiquitin carboxyl-terminal hydrolase 48; EC 3.1.2.15; Ubiquitin thioesterase 48; Ubiquitin-specific-processing protease 48; Deubiquitinating enzyme 48; USP48; USP31
<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 USP48.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	USP48 antibody detects endogenous levels of total USP48 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 91%, Positives = 100%); Rat (Identities = 91%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100 ELISA: 1:5000
<b>References:</b>	Lockhart P.J., DNA Seq. 15:9-14(2004). Gregory S.G., Nature 441:315-321(2006). Ota T., Nat. Genet. 36:40-45(2004).