

## SENP3 Antibody [#C0367]

<b>Catalog Number:</b>	C0367
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q9H4L4
<b>Other Names:</b>	Sentrin-specific protease 3; EC=3.4.22.-; Sentrin/SUMO-specific protease SENP3; SUMO-1-specific protease 3; SSP3; SUSP3
<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 SENP3.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	SENP3 antibody detects endogenous levels of total SENP3 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 91%, Positives = 100%)
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:5000
<b>References:</b>	Yeh E.T.H., Gene 248:1-14(2000). Choi S.J., Submitted (OCT-1999) to the EMBL/GenBank/DDBJ databases. Wiemann S., Genome Res. 11:422-435(2001).