

## TSG6 Antibody [#C10812]

<b>Catalog Number:</b>	C10812
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P98066
<b>Other Names:</b>	TSG6, tumor necrosis factor, alpha-induced protein 6
<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 TSG6.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	TSG6 antibody detects endogenous levels of total TSG6 protein.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:5000
<b>References:</b>	Lee T.H., J. Cell Biol. 116:545-557(1992). Nentwich H.A., J. Biol. Chem. 277:15354-15362(2002). Wisniewski H.-G., Biochemistry 33:7423-7429(1994).