

## ST5 Antibody [#C10797]

<b>Catalog Number:</b>	C10797
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P78524
<b>Other Names:</b>	Suppression of tumorigenicity 5, HeLa tumor suppression 1, DENN domain-containing protein 2B, ST5, DENND2B, HTS1
<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 ST5.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	ST5 antibody detects endogenous levels of total ST5 protein.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Lichy J.H., Nucleic Acids Res. 24:4700-4708(1996). Amid C., Cytogenet. Cell Genet. 93:284-290(2001). Lichy J.H., Cell Growth Differ. 3:541-548(1992).