

## STAT1 Antibody [#C10603]

<b>Catalog Number:</b>	C10603
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P42224
<b>Other Names:</b>	Signal transducer and activator of transcription 1, Signal transducer and activator of transcription 1-alpha/beta, Transcription factor ISGF-3 components p91/p84
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), 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 STAT1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	STAT1 antibody detects endogenous levels of total STAT1 protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000   ELISA: 1:20000
<b>References:</b>	Schindler C., Proc. Natl. Acad. Sci. U.S.A. 89:7836-7839(1992). Hillier L.W., Nature 434:724-731(2005). Yan R., Nucleic Acids Res. 23:459-463(1995).