

## STAT3 Antibody [#C10589]

<b>Catalog Number:</b>	C10589
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P40763
<b>Other Names:</b>	APRF, Acute-phase response factor, Signal transducer and activator of transcription 3
<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 C-terminal of human STAT3.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	STAT3 antibody detects endogenous levels of total STAT3 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:5000
<b>References:</b>	Akira S., Cell 77:63-71(1994). Della Pietra L., Gene 213:119-124(1998). Zhang X., Science 267:1990-1994(1995). Giraud S., J. Biol. Chem. 277:8004-8011(2002).