

## SSBP2 Antibody [#C10803]

<b>Catalog Number:</b>	C10803
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P81877
<b>Other Names:</b>	Sequence-specific single- stranded-DNA-binding protein 2; Single-stranded DNA-binding protein 2; SSB2; SSDP2
<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 SSBP2.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	SSBP2 antibody detects endogenous levels of total SSBP2 protein.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:10000
<b>References:</b>	Castro P.D., Genomics 80:78-85(2002). Hu R.-M., Proc. Natl. Acad. Sci. U.S.A. 97:9543-9548(2000). Wiemann S., Genome Res. 11:422-435(2001).