

MtSSB Antibody [#C10880]

Catalog Number:	C10880
Concentration:	1mg/ml
Swiss-Prot No.:	Q04837
Other Names:	Mt-SSB, MtSSB, PWP1-interacting protein 17, SSB, SSBP, single-stranded DNA-binding protein, mitochondrial
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg^{2+} and Ca^{2+}), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized peptide derived from internal of human MtSSB.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MtSSB antibody detects endogenous levels of total MtSSB protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000
References:	Tiranti V., Gene 126:219-225(1993). Matsuoka S., Science 316:1160-1166(2007). Yang C., Nat. Struct. Biol. 4:153-157(1997).