

SF1 (Ab-82) Antibody [#B0572]

Catalog Number:	B0572
Concentration:	1mg/ml
Swiss-Prot No.:	Q15637
Other Names:	BBP, CW17, Mammalian branch point binding protein mBBP, SF01, Splicing factor 1, Splicing factor SF1, Transcription factor ZFM1, ZFM1, ZFP162, ZNF162, Zinc finger gene in MEN1 locus, Zinc finger protein 162, Mzfm
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 non-phosphopeptide derived from human SF1 around the phosphorylation site of serine 82 (S-P-S ^P -P-E).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	SF1 (Ab-82) antibody detects endogenous levels of total SF1 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
References:	Susan Y. Park, Development, May 2005; 132: 2415 - 2423. Dujin Zhou, Mol. Endocrinol., Jul 2000; 14: 986. Dagmar Wilhelm, Genes & Dev., Jul 2002; 16: 1839. Philipp Gut, Development, Oct 2005; 132: 4611 - 4619.