

53BP1 (Ab-6) Antibody [#B0760]

Catalog Number:	B0760
Concentration:	1mg/ml
Swiss-Prot No.:	Q12888
Other Names:	P531, TP53B, TP53BP1, TRP53BP1, Tumor suppressor p53-binding protein 1, p53-binding protein 1
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 53BP1 around the phosphorylation site of serine 6 (T-G-S ^P -Q-L).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	53BP1 (Ab-6) antibody detects endogenous levels of total 53BP1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:10000
References:	Iwabuchi K., J. Biol. Chem. 273:26061-26068(1998). Iwabuchi K., Proc. Natl. Acad. Sci. U.S.A. 91:6098-6102(1994). Beausoleil S.A., Proc. Natl. Acad. Sci. U.S.A. 101:12130-12135(2004).