

## 53BP1(Phospho-Ser25) Antibody [#8A0886]

<b>Catalog Number:</b>	8A0886
<b>Concentration:</b>	1mg/ml
<b>Origin of Antibody:</b>	Rabbit
<b>Swiss-Prot No.:</b>	Q12888
<b>Other Names:</b>	P531, TP53B, TP53BP1, TRP53BP1, Tumor suppressor p53-binding protein 1, p53-binding protein 1.
<b>All Sites:</b>	Human: Ser25; Mouse: Ser25; Rat: Ser30
<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 phosphopeptide derived from human 53BP1 around the phosphorylation site of serine 25 (E-D-SP-Q-P).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific phosphopeptide. The antibody against non-phosphopeptide was removed by chromatography using non-phosphopeptide corresponding to the phosphorylation site.
<b>Specificity:</b>	53BP1 (Phospho-Ser25) antibody detects endogenous levels of 53BP1 only when phosphorylated at serine 25.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100 IF: 1:100~1:500 ELISA: 1:40000
<b>References:</b>	Irene M. Ward. Mol. Cell. Biol., Nov 2005; 25: 10079 - 10086. Fiona Pryde. J. Cell Sci., May 2005; 118: 2043 - 2055. Julio C. Morales. J. Biol. Chem., Apr 2003; 278: 14971 - 14977.