

p53(Phospho-Ser15) Antibody [#8A7180]

Catalog Number:	8A7180
Concentration:	1mg/ml
Origin of Antibody:	Rabbit
Swiss-Prot No.:	P04637
Other Names:	Antigen NY-CO-13; Cellular tumor antigen p53; Phosphoprotein p53; TP53; tumor protein p53; Tumor suppressor p53
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 phosphopeptide derived from human p53 around the phosphorylation site of Serine 15 (P-L-S ^P -Q-E).
Purification:	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.
Specificity:	p53(Phospho-Ser15) Antibody detects endogenous levels of p53 only when phosphorylated at Serine15.
Reactivity:	Human, Rat.
Applications:	IHC: 1:50~1:100 ELISA: 1:10000 WB: 1:500~1:1000 IP: Various Dilution
References:	Shieh, S. Y. et al. (1999) <i>EMBO J.</i> 18, 1815-1823. Honda, R. et al. (1997) <i>FEBS Lett.</i> 420, 25-27. Hirao, A. et al. (2000) <i>Science</i> 287, 1824-1827.