

## p53 (Ab-378) Antibody [#B8054]

<b>Catalog Number:</b>	B8054
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P04637
<b>Other Names:</b>	Antigen NY-CO-13; Cellular tumor antigen p53; Phosphoprotein p53; TP53; tumor protein p53; Tumor suppressor p53
<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 non-phosphopeptide derived from human p53 around the phosphorylation site of serine 378 (S-T-S <sup>P</sup> -R-H).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	p53 (Ab-378) antibody detects endogenous levels of total p53 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 92%, Positives = 92%); Rat (Identities = 92%, Positives = 92%)
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	Zakut-Houri R., EMBO J. 4:1251-1255(1985). Lamb P., Mol. Cell. Biol. 6:1379-1385(1986). Harlow E., Mol. Cell. Biol. 5:1601-1610(1985).