

## p53 (Ab-379) Antibody [#D0036]

<b>Catalog Number:</b>	D0036
<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 $Mg^{2+}$ and $Ca^{2+}$ ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized non-acetylated peptide derived from human p53 around the acetylation site of lysine 379 (R-H-K <sup>A</sup> -K-L).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	p53 (Ab-379) antibody detects endogenous levels of total p53 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000
<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).