

p53 Antibody [#C10276]

Catalog Number:	C10276
Concentration:	1mg/ml
Swiss-Prot No.:	P04637
Other Names:	Antigen NY-CO-13, Cellular tumor antigen p53, Phosphoprotein p53, TP53, 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 peptide derived from internal of human p53.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	p53 antibody detects endogenous levels of total p53 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	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).