

p53 (Acetyl-Lys317) Antibody [#D0015]

Catalog Number:	D0015
Concentration:	1mg/ml
Swiss-Prot No.:	P04637
Other Names:	Antigen NY-CO-13, Cellular tumor antigen p53, Phosphoprotein p53, TP53, Tumor suppressor p53
All Sites:	Human: Lys317; Mouse: Lys314; Rat: Lys315
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 acetylated peptide derived from human p53 around the acetylation site of lysine 317 (Q-P-K ^A -K-K).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific acetylated peptide. The antibody against non-acetylated peptide was removed by chromatography using non-acetylated peptide corresponding to the acetylation site.
Specificity:	p53 (Acetyl-Lys317) antibody detects endogenous levels of p53 only when acetylated at lysine 317.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IF: 1:100~1:500 ELISA: 1:5000
References:	Kanashiro C.A., Proc. Natl. Acad. Sci. U.S.A. 100:15836-15841(2003). Anderson C.W., Radiat. Res. 154:473-476(2000). Shiseki M., Cancer Res. 63:2373-2378(2003).