

p53 (Ab-317) Antibody [#D0035]

Catalog Number:	D0035
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 non-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 immunogen.
Specificity:	p53 (Ab-317) antibody detects endogenous levels of total p53 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
References:	Lin Liu, Mol. Cell. Biol., Feb 1999; 19: 1202 - 1209. James J. Lee, AACR Meeting Abstracts, Apr 2005; 2005: 511. Kazuyasu Sakaguchi, Genes & Dev., Sep 1998; 12: 2831.