

p53 (Ab-37) Antibody [#B7183]

- Catalog Number:** B7183-1, B7183-2, B7183
Amount: 50µg/50µl, 100µg/100µl, 200µg/200µl
Form of Antibody: Rabbit IgG in phosphate buffered saline (without Mg²⁺ and Ca²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Storage/Stability: Store at -20°C/1 year
Immunogen: The antiserum was produced against synthesized non-phosphopeptide derived from human p53 around the phosphorylation site of Serine 37.
Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity/Sensitivity: p53(Ab-37) Antibody detects endogenous levels of p53 protein around Serine 37.
Reactivity: Human
Applications: WB: 1:500~1:1000 ELISA: 1:20000
Swiss-Prot No. : P04637
References: Dhavan, R. and Tsai, L.H. (2001) Nat Rev Mol Cell Biol. 2: 749-759.
Patrick, G. N. et al. (1998) J Biol Chem. 273: 24057-24064.
Mayo LD, et al.(2005) J Biol Chem. 280(28):25953-25959.
Thompson T, et al.(2004) J Biol Chem. 279(51): 53015-53022.
Jackson MW, et al.(2004) Oncogene. 23(25):4477-4487.