

p53 (Ab-6) Antibody [#B7185]

- Catalog Number:** B7185-1, B7185-2, B7185
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 6.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity:** p53 (Ab-6) antibody detects endogenous levels of total p53 protein.
- Reactivity:** Human, Mouse
- Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100
ELISA: 1:20000
- Swiss-Prot No. :** P04637
- References:** Thompson T, et al. (2004) J Biol Chem; 279(51): 53015-22.
Jackson MW, et al. (2004) Oncogene.
Hao, M. et al. (1996) J. Biol. Chem. 271, 29380-29385.
Lu, H. et al. (1997) Mol. Cell. Biol. 17, 5923-5934.
Lohrum, M. and Scheidtmann, K.H. (1996) Oncogene 13, 2527-2539.