

HSP90 cyto Antibody [#C0234]

- Catalog Number:** C0234-1, C0234-2, C0234
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No.:** P07900
- All Names:** HSP89-ALPHA, HSP89A, HSPC1, HSPCA, LAP2, HEAT-SHOCK 90-KD PROTEIN 1 ALPHA, LIPOPOLYSACCHARIDE-ASSOCIATED PROTEIN 2, LPS-ASSOCIATED PROTEIN 2
- 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 peptide derived from human HSP90 cyto.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** HSP90 cyto antibody detects endogenous levels of total HSP90 cyto protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:300~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:40000
- References:** Kai Yang, Mol. Biol. Cell, Mar 2006; 17: 1461 - 1471.
Adnan Ali, Mol. Cell. Biol., Sep 1998; 18: 4949 - 4960.
Hee Chul Lee, J. Biol. Chem., Feb 2002; 277: 7430 - 7437.
Kazuya Terasawa, J. Biochem. (Tokyo), Apr 2005; 137: 443 - 447.