

## HSP27 (Ab-15) Antibody [#B7110]

- Catalog Number:** B7110-1, B7110-2, B7110
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Form of Antibody:** Rabbit IgG in phosphate buffered saline (without Mg<sup>2+</sup> and Ca<sup>2+</sup>), 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 HSP27 around the phosphorylation site of serine 15.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity:** HSP27 (Ab-15) antibody detects endogenous levels of total HSP27 protein.
- Reactivity:** Human
- Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100  
ELISA: 1:20000
- Swiss-Prot No. :** P04792
- References:** Tremolada L, et al. (2005) Proteomics. 5(3): 788-795.  
McGregor E, et al. (2004) Mol Cell Proteomics. 3(2): 115-124.  
Benn SC, et al. (2002) Neuron. 36(1): 45-56.  
MacDonald JA, et al. (2002) Mol Cell Proteomics. 1(4): 314-322.