

## GRP75 Antibody [#C0216]

- Catalog Number:** C0216-1, C0216-2, C0216
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No.:** P38646
- All Names:** HSPA9B, MOT2, HEAT-SHOCK 70-KD PROTEIN 9B, MORTALIN 2, MORTALIN, PERINUCLEAR, heat shock 70kDa protein 9B (mortalin-2)
- 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 peptide derived from human GRP75.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** GRP75 antibody detects endogenous levels of total GRP75 protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:300~1:1000    IHC: 1:50~1:100    IF: 1:500~1:1000    ELISA: 1:5000
- References:** LA Mizzen, J. Biol. Chem., Dec 1989; 264: 20664 - 20675.  
Laura E. Lantry, J. Nucl. Med., Jul 2006; 47: 1144 - 1152.  
Xiaoyuan Chen, J. Nucl. Med., Aug 2004; 45: 1390 - 1397.  
Charles F. Gillespie, J Biol Rhythms, Jun 1996; 11: 137 - 144.