

## GRP78 Antibody [#C0217]

- Catalog Number:** C0217-1, C0217-2, C0217  
**Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl  
**Swiss-Prot No.:** P11021  
**All Names:** BIP, GLUCOSE-REGULATED PROTEIN 78-KD, HEAT-SHOCK 70-KD PROTEIN 5, IMMUNOGLOBULIN HEAVY CHAIN-BINDING PROTEIN, heat shock 70kDa protein 5 (glucose-regulated protein, 78kDa)  
**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 GRP78.  
**Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.  
**Specificity/Sensitivity:** GRP78 antibody detects endogenous levels of total GRP78 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:** Uma K. Misra, J. Biol. Chem., Oct 2002; 277: 42082 - 42087.  
CA Clairmont, J. Biol. Chem., Feb 1992; 267: 3983 - 3990.  
J Tooze, J. Cell Biol., Jul 1989; 109: 35.  
Jong K. Yun, PNAS, Dec 1997; 94: 13903.