

## p70 S6 Kinase (Ab-421/424) Antibody [#B7190]

- Catalog Number:** B7190-1, B7190-2, B7190  
**Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl  
**Swiss-Prot No. :** P23443  
**All Names:** p70 S6 KINASE ALPHA, ribosomal protein S6 kinase, 70kDa, polypeptide 1  
**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 p70 S6 Kinase around the phosphorylation site of threonine 421 or serine 424(R-T-P-V-S).  
**Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.  
**Specificity/Sensitivity** p70 S6 Kinase (Ab-421/424) antibody detects endogenous levels of total p70 S6 Kinase protein.  
**Reactivity:** Human, Mouse, Rat  
**Applications:** IHC: 1:50~1:100      ELISA: 1:10000  
**References:** BA Moser, Mol. Cell. Biol., Sep 1997; 17: 5648 - 5655.  
Nicholas Lehman, FASEB J, Apr 2007; 21: 1075 - 1087.  
Bernard Masri, FASEB J, Sep 2004; 10.1096/fj.04-1930fje.  
Julian Gomez-Cambronero, J. Immunol., Dec 2003; 171: 6846.