

4E-BP1 (Ab-69) Antibody [#B0402]

- Catalog Number:** B0402-1, B0402-2, B0402
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** Q13541
- All Names:** 4EBP1, EIF4EBP1, Eukaryotic translation initiation factor 4E binding protein 1, Insulin-stimulated EIF-4E binding protein PHAS-I, P/OKCL.6, PHAS-1, PHAS-I
- 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 non-phosphopeptide derived from human 4E-BP1 around the phosphorylation site of threonine 70 (A-K-T-P-P).
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** 4E-BP1 (Ab-69) antibody detects endogenous levels of total 4E-BP1 protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100
ELISA: 1:5000
- References:** Anne-Claude Gingras, Genes & Dev., Jun 1999; 13: 1422.
Guangming Liu, J. Biol. Chem., Mar 2002; 277: 8810 - 8816.
Ya-Ping Jiang, J. Biol. Chem., Mar 2001; 276: 10943 - 10951.
Anne-Claude Gingras, Genes & Dev., Feb 1998; 12: 502.