

4E-BP1 (Ab-45) Antibody [#B7002]

- Catalog Number:** B7002-1, B7002-2, B7002
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/OKCL6, 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 45 (S-T-T^P-P-G).
Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity/Sensitivity: 4E-BP1 (Ab-45) antibody detects endogenous levels of total 4E-BP1 protein. This antibody may cross-react with 4E-BP2 and 4E-BP3.
Reactivity: Human, Mouse, Rat
Applications: WB: 1:500~1:1000 IHC: 1:50~1:100
ELISA: 1:20000
References: Gingras AC, et al. (1998) Genes Dev 12(4): 502-513.
Brugarolas J, et al. (2004) Genes Dev 18(23): 2893-2904.
Kumar V, et al. (2000) EMBO J 19(5): 1087-1097.
Moody CA, et al. (2005) J Virol 79(9): 5499-5506.
Burnett PE, et al. (1998) Proc Natl Acad Sci U S A 95(4): 1432-1437.