



---

## 4E-BP1 (Phospho-Thr36) Antibody

Catalog Number 11222

(For Research Use Only)

---

**Catalog Number:** 11222-1, 11222-2, 11222

**Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl

**Swiss-Prot No. :** Q13541

**All Names:** 4E-BP1, eIF4E-binding protein 1, Phosphorylated heat- and acid-stable protein regulated by insulin 1, PHAS-I

**All Sites:** Human: Thr36 or Thr37; Mouse: Thr35; Rat: Thr35

**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 phosphopeptide derived from human 4E-BP1 around the phosphorylation site of threonine 36 (S-T-T<sup>P</sup>-P-G).

**Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific phosphopeptide. The antibody against non-phosphopeptide was removed by chromatography using non-phosphopeptide corresponding to the phosphorylation site.

**Specificity/Sensitivity:** 4E-BP1 (phospho-Thr36) antibody detects endogenous levels of 4E-BP1 only when phosphorylated at threonine 36. This antibody may cross-react with 4E-BP2 and 4E-BP3 when phosphorylated at equivalent sites.

**Reactivity:** Human, Mouse, Rat

**Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100  
ELISA: 1:10000

**References:** Gingras AC, et al. Genes Dev Nov. 2005 Feb 15.  
Mothe-Satney I, et al.(2000) J Biol Chem:15(21): 2852-64.  
Gingras AC, et al.(1999) Genes Dev:13(11): 1422-37.  
Lal L, et al.(2005) Blood;105(4): 1669-7.

