

## Product Information

### EIF4EBP1

Product Number : **00-008**

#### Product description

Full-length human EIF4EBP1 [1-118(end) amino acids of accession number NP\_004086.1] was expressed as N-terminal GST-fusion protein (39 kDa) using E.coli expression system. GST-EIF4EBP1 was purified by using glutathione sepharose chromatography.

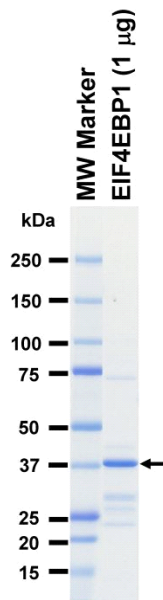
#### Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,  
1 mM DTT, 10% glycerol, pH7.5

#### Storage and Handling:

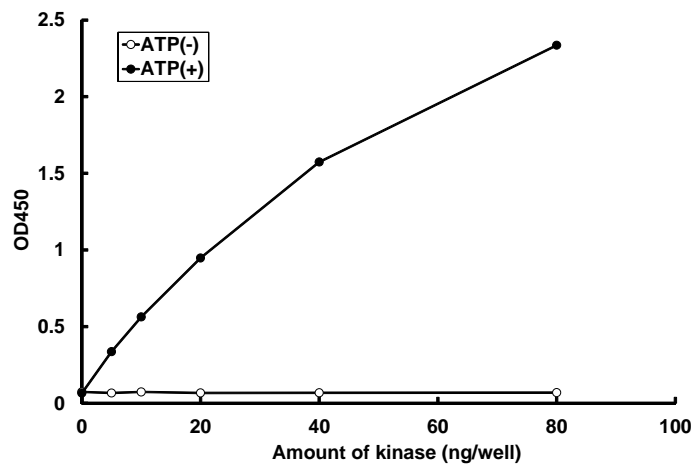
Store at -80C.  
Avoid repeating freeze-thaws.

#### SDS-PAGE



Purity: 65 %  
The purity was assessed by SDS-PAGE/CBB staining.

#### Activity data



The phosphorylational ability of the protein was determined by ELISA. The substrate protein was incubated with the indicated enzyme. Phosphorylation was detected by anti-phospho antibody and HRP-labeled anti-rabbit IgG (or HRP-labeled anti-mouse IgG).

Kinase : Erk2

ATP : 100 µM