

## Product Information

### GCN2(EIF2AK4)

 Product Number : **05-153**

#### Product description

Full-length human GCN2 [1-1649(end) amino acids and E556G of accession number NP\_001013725] was expressed as N-terminal GST-fusion protein (215 kDa) using baculovirus expression system. GST-GCN2 was purified by using glutathione sepharose chromatography.

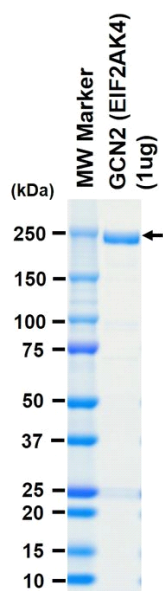
#### Storage buffer:

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.

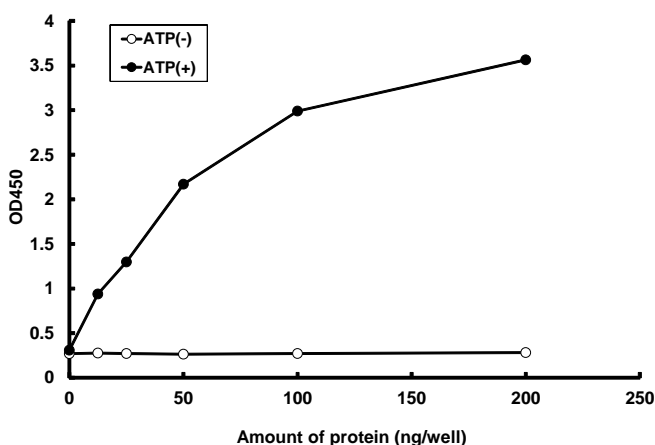
#### SDS-PAGE



Purity: 82 %

The purity was assessed by SDS-PAGE/CBB staining.

#### Activity data



The activity was determined by ELISA. The enzyme was incubated with GST-fused substrate protein and tRNA, and after stopping kinase reaction by EDTA, the reaction solution was transferred into glutathione-coated plate. Phosphorylation was detected by anti-phospho antibody and HRP-labeled anti-rabbit IgG.

Substrate : EIF2S1

 ATP : 100  $\mu$ M