

Product Information

BTN-GCN2(EIF2AK4)

Product Number : **05-453-20N**

Product description

Full-length human GCN2 [1-1649(end) amino acids and E556G of accession number NP_001013725] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (192 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

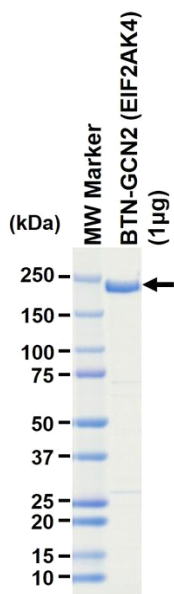
Storage buffer:

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.

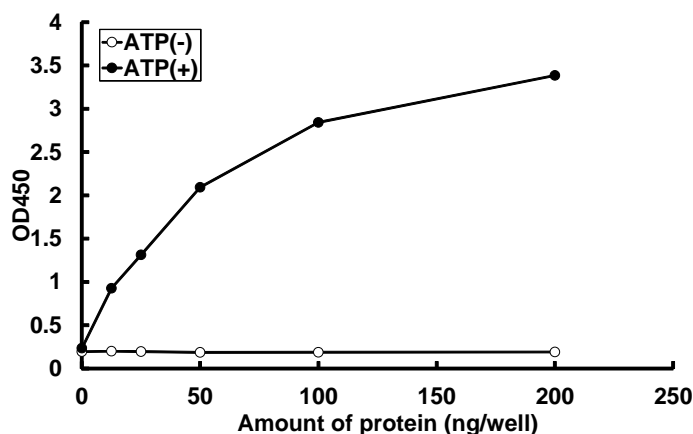
SDS-PAGE



Purity: 88 %

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 μ M