

Product Information

BUBR1(BUB1B)

 Product Number : **05-105**

Product description

Full-length human BUBR1 [1-1050(end) amino acids of accession number NP_001202.4] was expressed as N-terminal GST-fusion protein (147 kDa) using baculovirus expression system. GST-BUBR1 was purified by using glutathione sepharose chromatography.

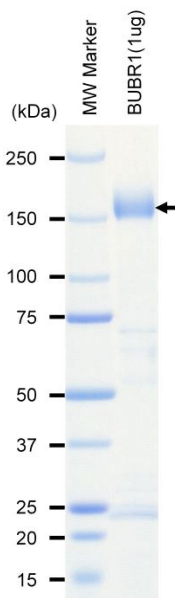
Storage buffer:

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.

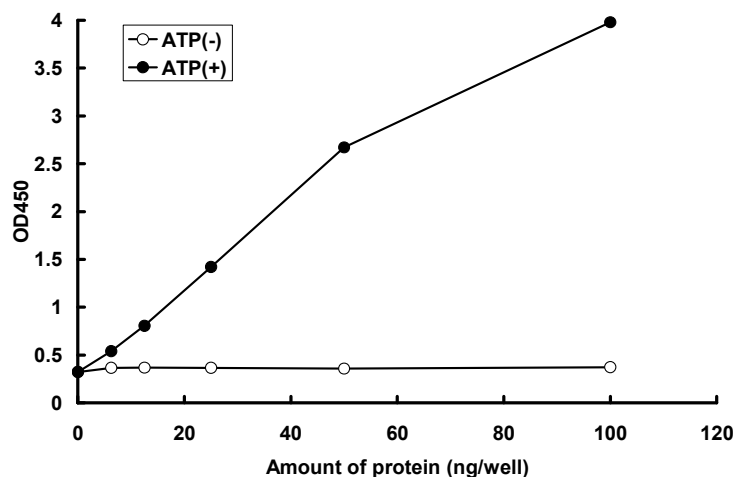
SDS-PAGE



Purity: 74 %

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

Activity data



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

Substrate : Histone H1

 ATP : 100 μ M