

## Product Information

### HH498(TNNI3K)

 Product Number : **09-112**

#### Product description

Human HH498, catalytic domain [1-727 amino acid of accession number NP\_057062.1] was expressed as N-terminal GST-fusion protein (108 kDa) using baculovirus expression system. GST-HH498 was purified by using glutathione sepharose chromatography.

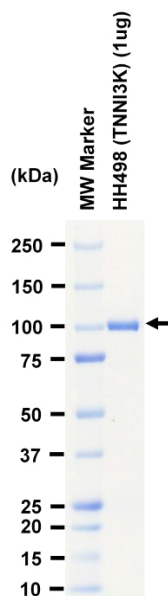
#### Storage buffer:

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

#### Storage and Handling:

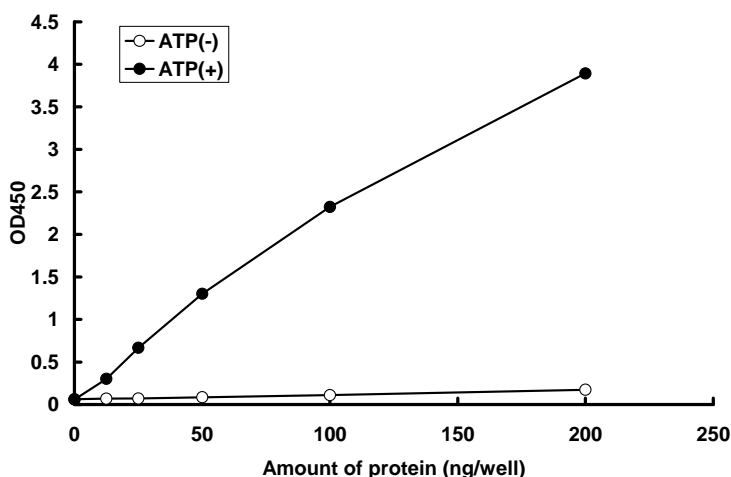
Store at -80C.  
 Avoid repeating freeze-thaws.

#### SDS-PAGE



Purity: 97 %  
 The purity was assessed by SDS-PAGE/CBB staining.

#### Activity data



The activity was determined by ELISA. The enzyme was incubated with biotinylated peptide in streptavidin-coated ELISA plate. Phosphorylation was detected by HRP-labeled anti-phospho antibody.

Substrate : Poly-(GT)

ATP : 100  $\mu$ M