

Product Information

TTK

Product Number : **05-169**

Product description

Full-length human TTK [1-857(end) amino acids of accession number NP_003309.2] was expressed as N-terminal GST-fusion protein (124 kDa) using baculovirus expression system. GST-TTK was purified by using glutathione sepharose chromatography.

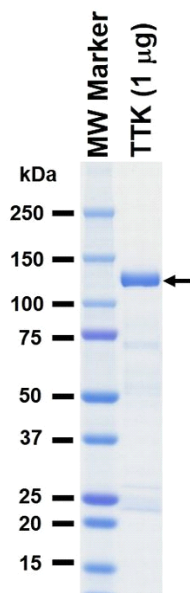
Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.1% CHAPS,
 1 mM DTT, 10% glycerol, pH 7.5

Storage and Handling:

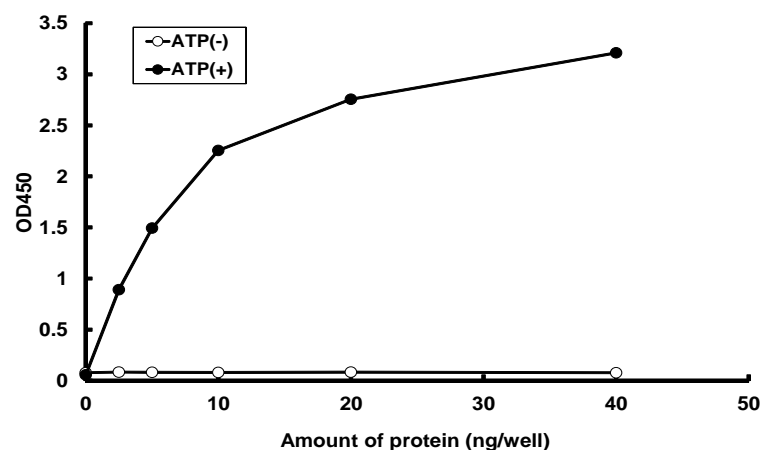
Store at -80C.
 Avoid repeating freeze-thaws.

SDS-PAGE



Purity: 78 %
 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 : Lyn substrate peptide

ATP : 20 μ M