

Product Information

STLK3(STK39)

 Product Number : **07-130**

Product description

Full-length human STLK3 [1-545(end) amino acids of accession number NP_037365.2] was co-expressed as N-terminal GST-fusion protein (87 kDa) with His-tagged WNK1 [1-491 amino acids of accession number NP_061852.1] using baculovirus expression system. GST-STLK3 was purified by using glutathione sepharose chromatography and Ni-NTA chromatography.

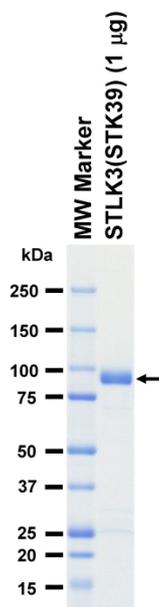
Storage buffer:

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.

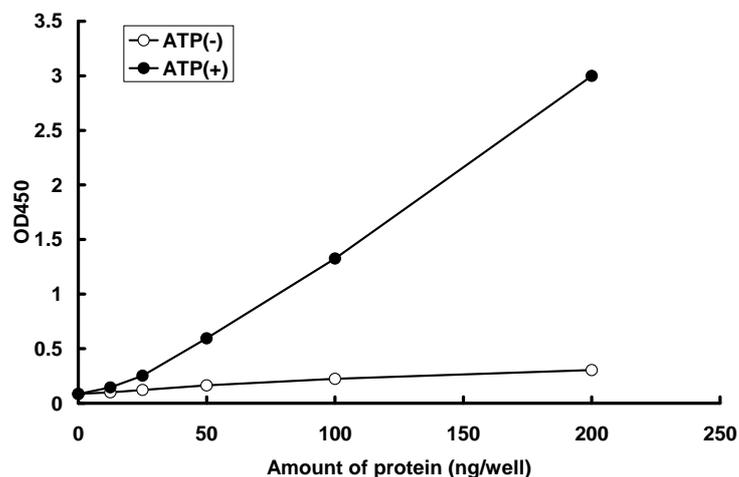
SDS-PAGE



Purity: 93 %

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/peptide, 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 (or HRP-labeled anti-mouse IgG).

Substrate : CATCHtide

 ATP : 100 μ M