

## Product Information

### VRK1

 Product Number : **03-110**

#### Product description

Full-length human VRK1 [1-396(end) amino acids of accession number NP\_003375.1] was expressed as N-terminal GST-fusion protein (72 kDa) using E. coli expression system. GST-VRK1 was purified by using glutathione sepharose chromatography.

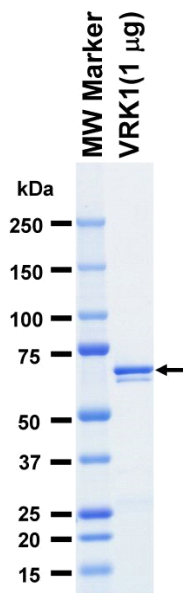
#### Storage buffer:

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.

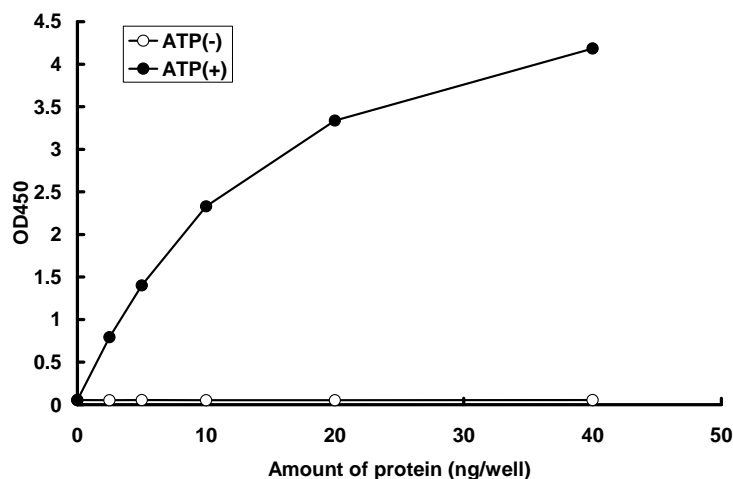
#### SDS-PAGE



Purity: 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 anti-phospho antibody and HRP-labeled anti-rabbit IgG (or HRP-labeled anti-mouse IgG).

Substrate : ATF2

 ATP : 100  $\mu$ M