

## Product Information

### VRK2

 Product Number : **03-111**

#### Product description

Human VRK2, catalytic domain [1-483 amino acids of accession number NP\_006287.2] was expressed as N-terminal GST-fusion protein (82 kDa) using baculovirus expression system. GST-VRK2 was purified by using glutathione sepharose chromatography.

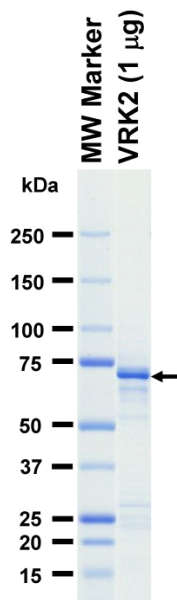
#### Storage buffer:

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.

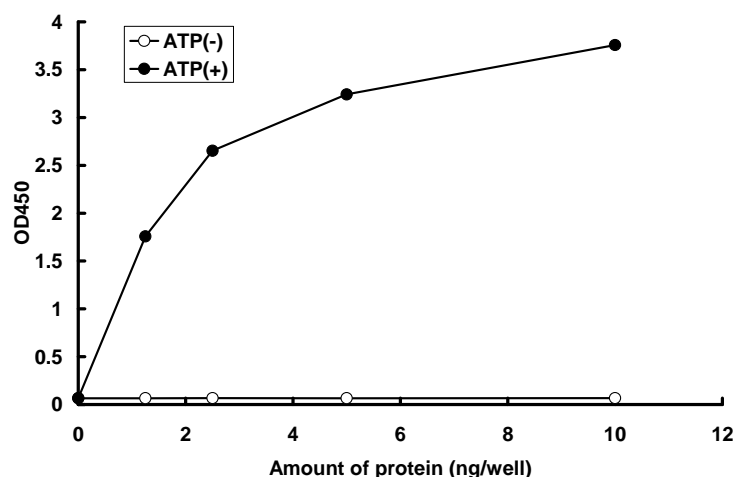
#### SDS-PAGE



Purity: 72 %

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