

## Product Information

### GPRK4(GRK4)

Product Number : **01-313**

#### Product description

Full-length human GPRK4 [1-578(end) amino acids of accession number NP\_892027.2] was expressed as N-terminal GST-fusion protein using baculovirus expression system. GST-GPRK4 was purified by using glutathione sepharose chromatography. GST-GPRK4 was cleaved by PreScission protease and GST-free GPRK4(67 kDa) was collected as flow-through fraction from glutathione sepharose chromatography.

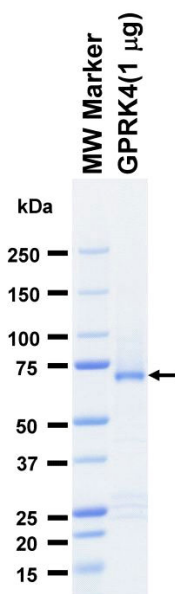
#### Storage buffer:

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

#### Storage and Handling:

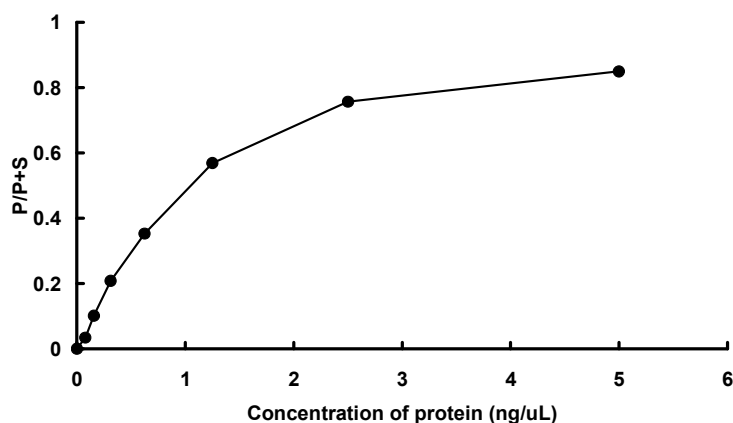
Store at -80C.  
Avoid repeating freeze-thaws.

#### SDS-PAGE



Purity: 78 %  
The purity was assessed by SDS-PAGE/CBB staining.

#### Activity data



The activity was measured by off-chip mobility shift assay. The enzyme was incubated with fluorescence-labeled substrate and Mg(or Mn)/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by LabChip™3000.

Substrate : Modified PLKtide

ATP : 100 µM