

## Product Information

### DMPK2(CDC42BPG)

 Product Number : **01-106**

#### Product description

Human DMPK2, catalytic domain [1-473 amino acids of accession number NP\_059995.2] was expressed as N-terminal GST-fusion protein (81 kDa) using baculovirus expression system. GST-DMPK2 was purified by using glutathione sepharose chromatography.

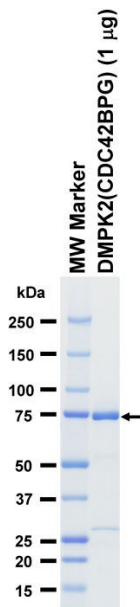
#### Storage buffer:

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.

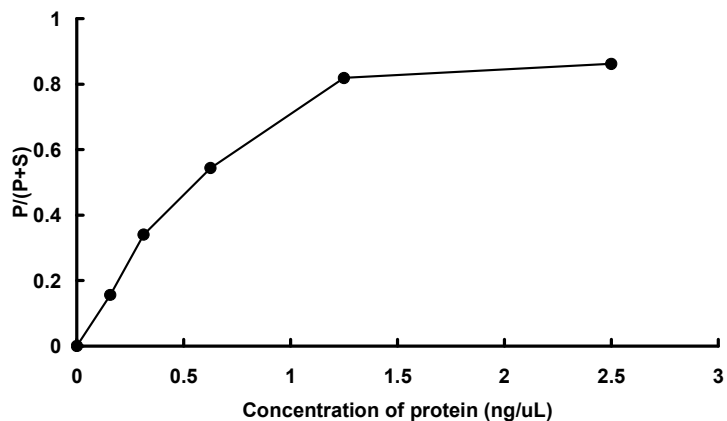
#### SDS-PAGE



Purity: 87 %

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 : DAPK1tide

ATP : 100 µM