

## Product Information

### MRCK $\beta$ (CDC42BPB)

Product Number : **01-108**

#### Product description

Truncated human MRCK $\beta$  [1-473 amino acids of accession number NP\_006026.3] was expressed as N-terminal GST-fusion protein (82 kDa) using baculovirus expression system. GST-MRCK $\beta$  was purified by using glutathione sepharose chromatography.

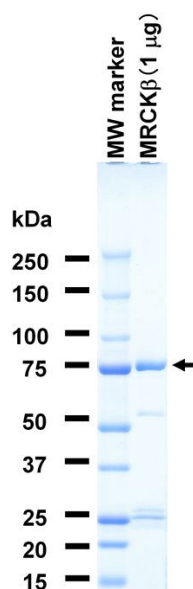
#### Storage buffer:

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

#### Storage and Handling:

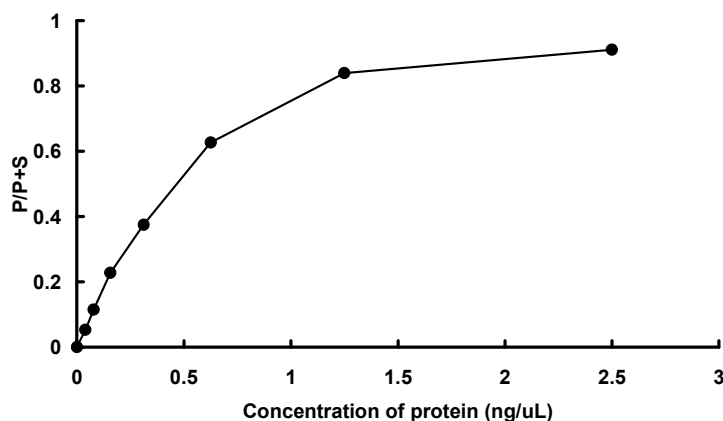
Store at -80C.  
Avoid repeating freeze-thaws.

#### SDS-PAGE



Purity: 61 %  
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<sup>TM</sup>3000.

Substrate : DAPK1tide

ATP : 100  $\mu$ M