

## Product Information

### MRCK $\alpha$ (CDC42BPA)

Product Number : **01-107**

#### Product description

Truncated human MRCK $\alpha$  [1-574 amino acids of accession number NP\_003598.2] was expressed as N-terminal GST-fusion protein (93 kDa) using baculovirus expression system. GST-MRCK $\alpha$  was purified by using glutathione sepharose chromatography.

#### Storage buffer:

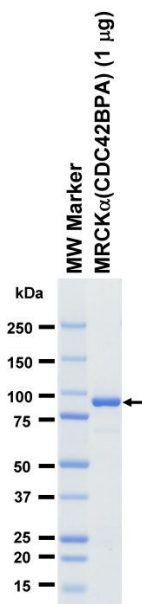
50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,  
1 mM DTT, 10% glycerol, pH7.5

#### Storage and Handling:

Store at -80C.

Avoid repeating freeze-thaws.

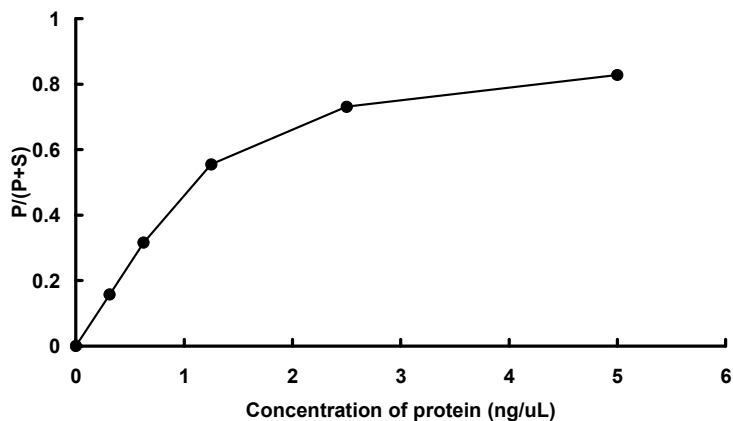
#### SDS-PAGE



Purity: 97 %

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  $\mu$ M