

## Product Information

### CGK2(PRKG2)

Product Number : **01-143**

#### Product description

Full-length human CGK2 [1-762(end) amino acids of accession number NP\_006250.1] was expressed as N-terminal GST-fusion protein (114 kDa) using baculovirus expression system. GST-CGK2 was purified by using glutathione sepharose chromatography.

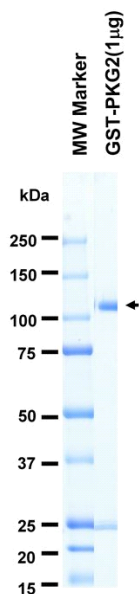
#### Storage buffer:

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.

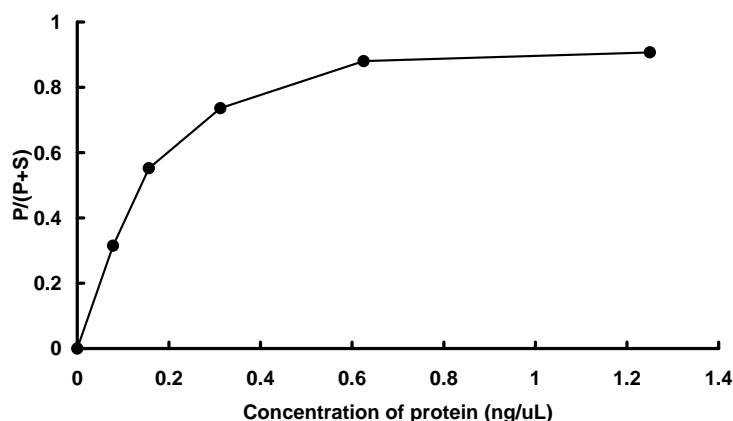
#### SDS-PAGE



Purity: 74 %

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, Mg(or Mn)/ATP, and cGMP. The phosphorylated and unphosphorylated substrates were separated and detected by LabChip™3000.

Substrate : Kemptide

ATP : 100  $\mu$ M