

## Product Information

### CLK1

Product Number : **04-126**

#### Product description

Human CLK1, catalytic domain [129-484(end) amino acids of accession number NP\_004062.2] was expressed as N-terminal GST-fusion protein (69 kDa) using baculovirus expression system. GST-CLK1 was purified by using glutathione sepharose chromatography.

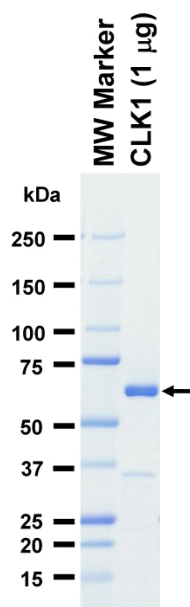
#### Storage buffer:

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

#### Storage and Handling:

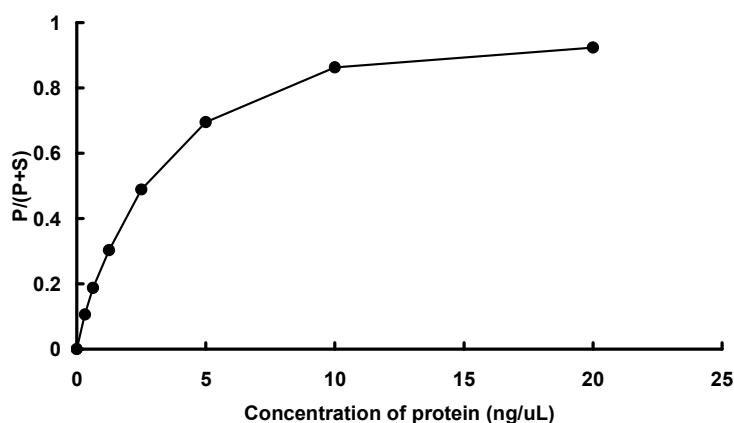
Store at -80C.  
Avoid repeating freeze-thaws.

#### SDS-PAGE



Purity: 92 %  
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 : DYRKtide

ATP : 100 µM