

## Product Information

### PDHK1(PDK1)

Product Number : **10-123**

#### Product description

Full-length human PDHK1[1-436(end) amino acids of accession number NP\_002601.1] was expressed as N-terminal GST-fusion protein (77 kDa) using baculovirus expression system. GST-PDHK1 was purified by using glutathione sepharose chromatography.

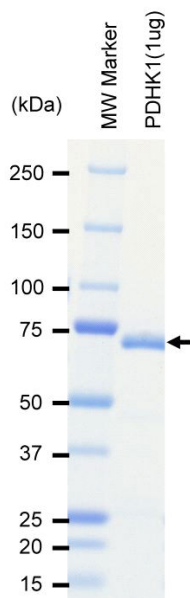
#### Storage buffer:

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

#### Storage and Handling:

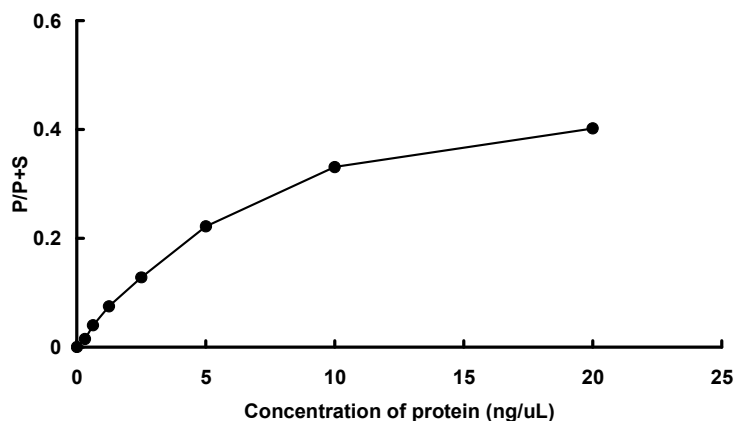
Store at -80C.  
Avoid repeating freeze-thaws.

#### SDS-PAGE



Purity: 95 %  
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 : PDHKtide

ATP : 100 µM