

Product Information

PGK(PRKG1)

Product Number : **01-142**

Product description

Full-length human PGK [1-686(end) amino acids of accession number NP_006249.1] was expressed as N-terminal GST-fusion protein (105 kDa) using baculovirus expression system. GST-PGK was purified by using glutathione sepharose chromatography.

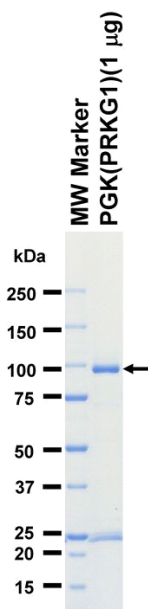
Storage buffer:

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.

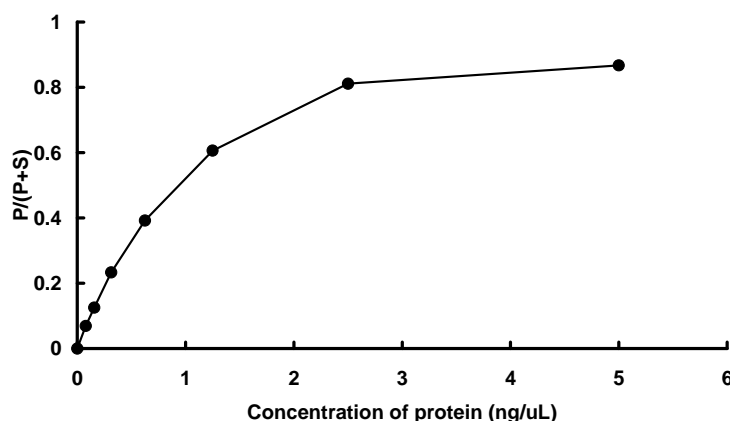
SDS-PAGE



Purity: 63 %

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 μ M