

## Product Information

### PBK

Product Number : **05-168**

#### Product description

Full-length human PBK [1-322(end) amino acids of accession number NP\_060962.2] was expressed as N-terminal GST-fusion protein (63 kDa) using baculovirus expression system. GST-PBK was purified by using glutathione sepharose chromatography.

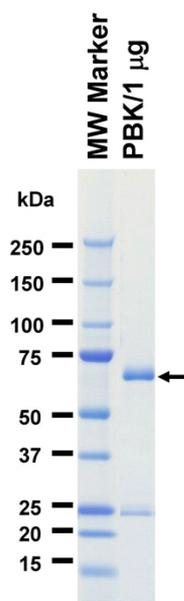
#### Storage buffer:

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

#### Storage and Handling:

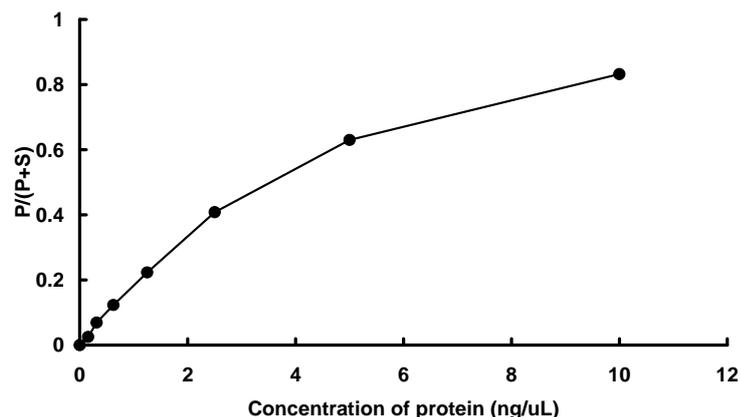
Store at -80C.  
 Avoid repeating freeze-thaws.

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

Substrate : Histone H3 peptide

ATP : 100 μM