

## Product Information

### HPK1(MAP4K1)

 Product Number : **07-410**

#### Product description

Human HPK1, catalytic domain [1-346 amino acids of accession number NP\_009112.1] was expressed as N-terminal DYKDDDDK tagged protein (41 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

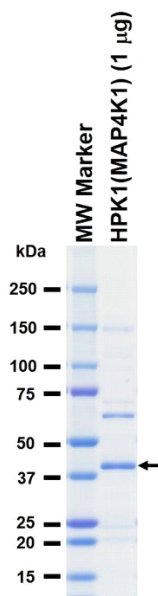
#### Storage buffer:

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.

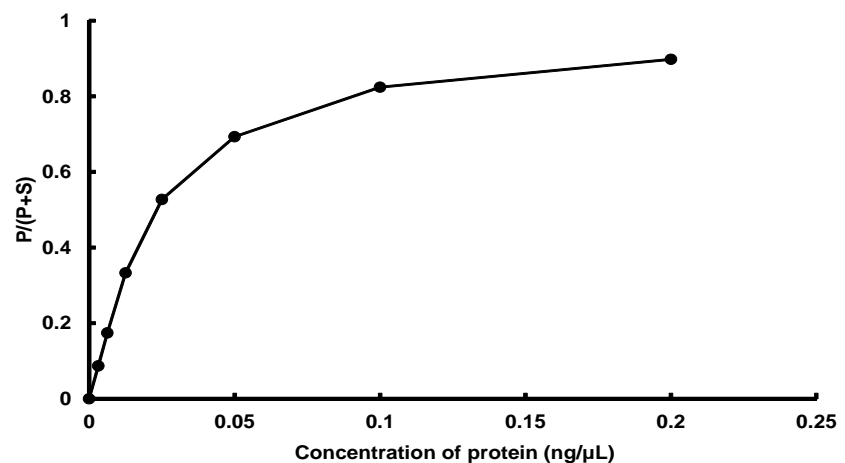
#### SDS-PAGE



Purity: 54 %

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 : S6K2 peptide

ATP : 100 µM