

## Product Information

### PEK(EIF2AK3)

Product Number : **05-155**

#### Product description

Human PEK, cytoplasmic domain [536-1116(end) amino acids and A704S of accession number NP\_004827.4] was expressed as N-terminal GST-fusion protein (94 kDa) using E.coli expression system. GST-PEK was purified by using glutathione sepharose chromatography and anion exchange chromatography.

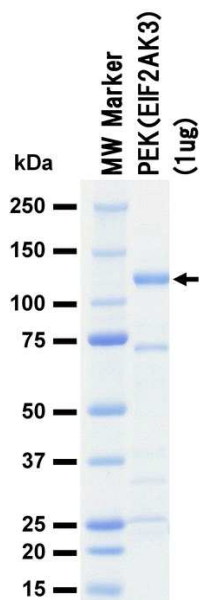
#### Storage buffer:

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.

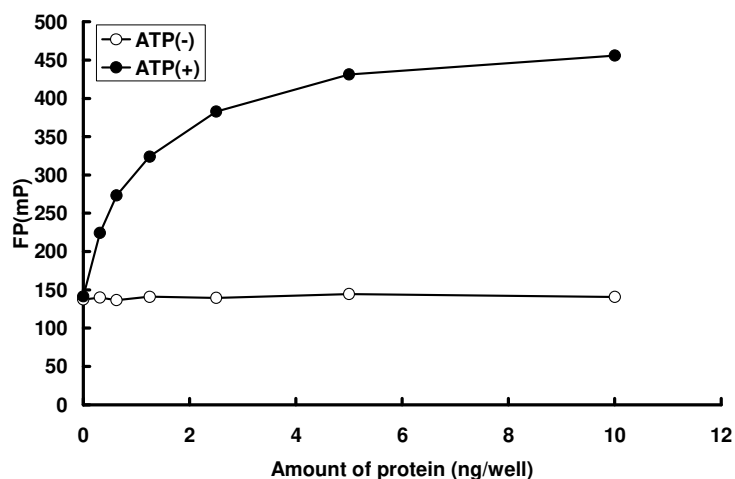
#### SDS-PAGE



Purity: 58 %

The purity was assessed  
by SDS-PAGE/CBB  
staining.

#### Activity data



The activity was determined by IMAP™ assay. The enzyme was incubated with fluorescein labeled peptide and phosphorylation was detected by IMAP™ technology (fluorescence polarization).

Substrate : SRPKtide

ATP : 100  $\mu$ M