

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.

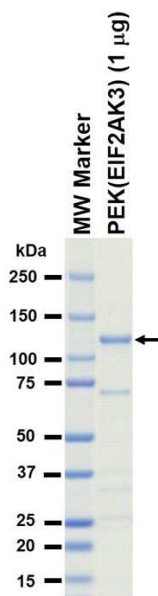
Storage buffer:

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.

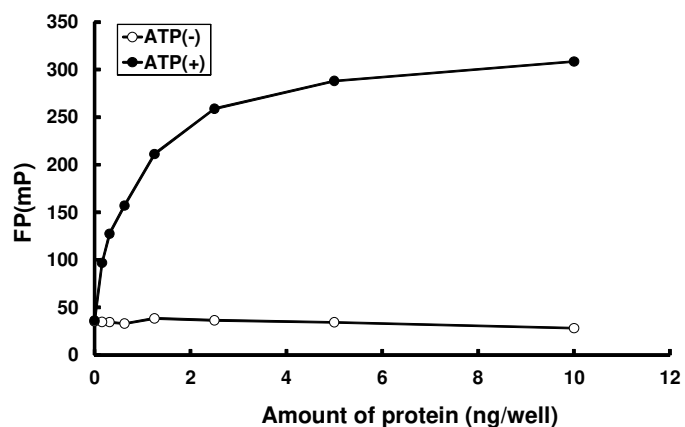
SDS-PAGE



Purity: 70 %

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 μ M