

Product Information

PKR(EIF2AK2)

Product Number : **05-156**

Product description

Human PKR , catalytic domain [252-551(end) amino acids of accession number NP_002750.1] was expressed as N-terminal GST-fusion protein (62 kDa) using baculovirus expression system. GST-PKR was purified by using glutathione sepharose chromatography and anion exchange chromatography.

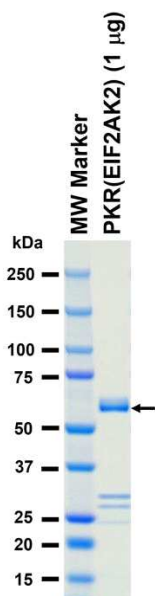
Storage buffer:

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.

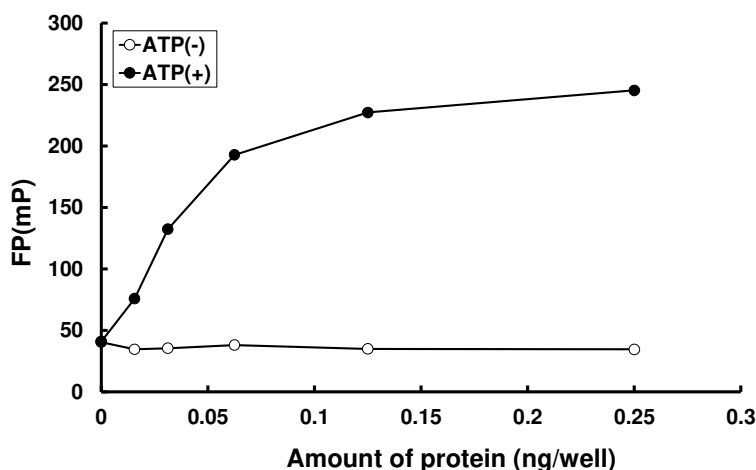
SDS-PAGE



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