

Product Information

p70S6K(RPS6KB1)

Product Number : **01-154**

Product description

Human p70S6K, catalytic domain [1-421 amino acids and T412E of accession number NP_003152.1] was expressed as N-terminal GST-fusion protein (75 kDa) using baculovirus expression system. GST-p70S6K was purified by using glutathione sepharose chromatography and anion exchange chromatography.

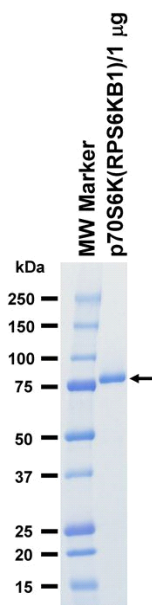
Storage buffer:

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.

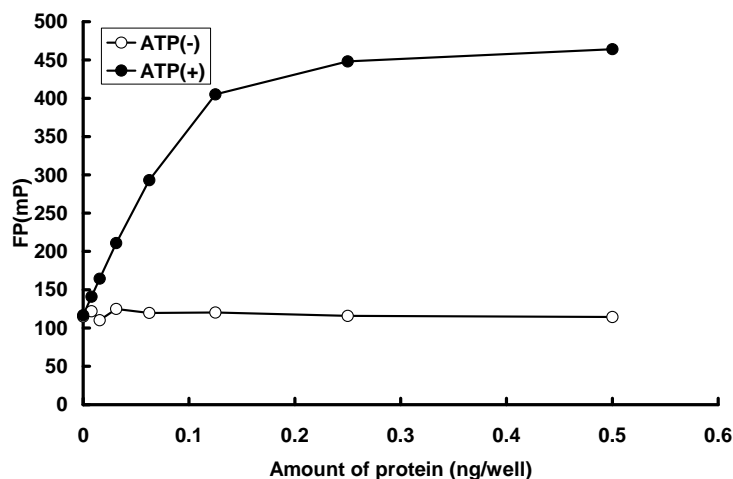
SDS-PAGE



Purity: 100 %

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 : S6K peptide

ATP : 100 µM