

Product Information

p70S6K(RPS6KB1)

Product Number : **01-156**

Product description

Human p70S6K, catalytic domain [1-421 amino acids and T412E of accession number NP_003152.1] was co-expressed as N-terminal GST-fusion protein (75 kDa) with His-tagged PDK1 [1-556(end) amino acids of accession number NP_002604.1] using baculovirus expression system. GST-p70S6K was purified by using glutathione sepharose chromatography.

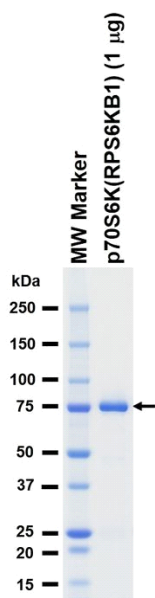
Storage buffer:

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.

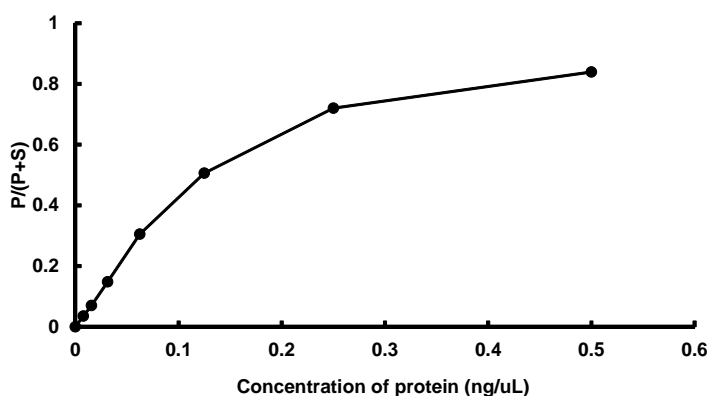
SDS-PAGE



Purity: 97 %

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