

## Product Information

### RSK2(RPS6KA3)

Product Number : 01-150

#### Product description

Full-length human RSK2 [1-740(end) amino acids of accession number NP\_004577.1] was expressed as N-terminal GST-fusion protein (111 kDa) using baculovirus expression system. GST-RSK2 was purified by using glutathione sepharose chromatography.

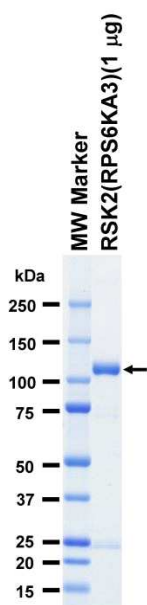
#### Storage buffer:

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.

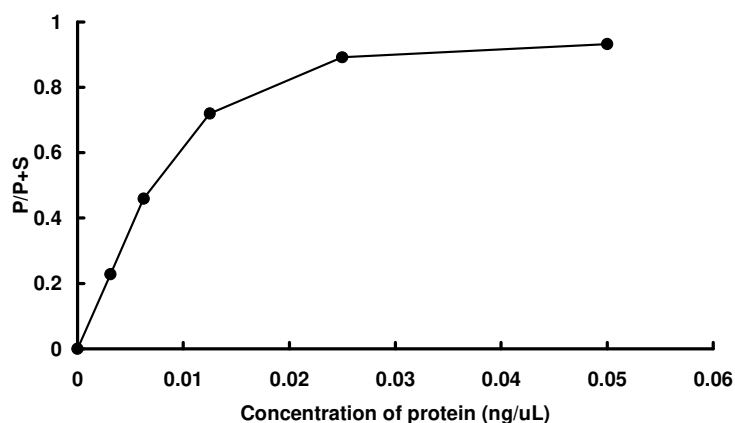
#### SDS-PAGE



Purity: 88 %

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

#### Activity data



The activity was measured by off-chip mobility shift assay(MSA). The enzyme was incubated with fluorescence-labeled substrate and Mg(or Mn)/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by MSA device.

Substrate : S6K peptide(N-FL)

ATP : 100 µM