

## Product Information

### MSK1(RPS6KA5)

Product Number : **01-147**

#### Product description

Full-length human MSK1 [1-802(end) amino acids of accession number NP\_004746.2] was co-expressed as N-terminal GST-fusion protein (117 kDa) with human His-tagged Erk2 [1-360 amino acids of accession number NP\_002736.3] using baculovirus expression system. GST-MSK1 was purified by using glutathione sepharose chromatography.

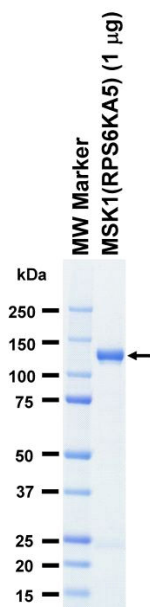
#### Storage buffer:

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.

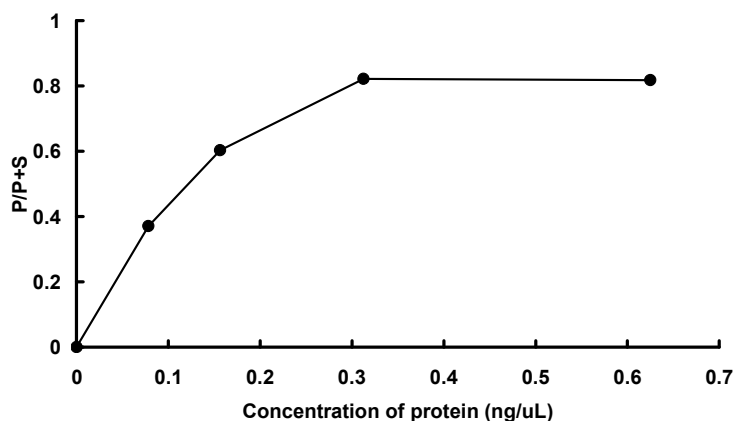
#### SDS-PAGE



Purity: 94 %

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 : Crosstide

ATP : 100 µM