

Product Information

MSK2(RPS6KA4)

Product Number : **01-148**

Product description

Full-length human MSK2 [1-772(end) amino acids of accession number NP_003933.1] was co-expressed as N-terminal GST-fusion protein (114 kDa) with human His-tagged Erk2 [1-360 amino acids of accession number NP_002736.3] using baculovirus expression system. GST-MSK2 was purified by using glutathione sepharose chromatography.

Storage buffer:

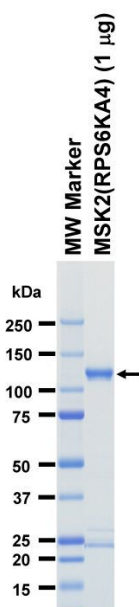
50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,
1 mM DTT, 10% glycerol, pH7.5

Storage and Handling:

Store at -80C.

Avoid repeating freeze-thaws.

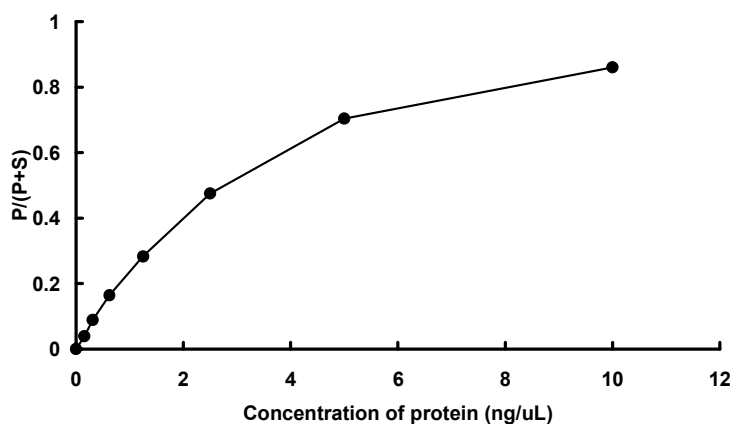
SDS-PAGE



Purity: 76 %

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 μ M