

Product Information

TSSK1B(TSSK1)

 Product Number : **02-364**

Product description

Full-length human TSSK1B [1-367(end) amino acids of accession number NP_114417.1] was expressed as N-terminal GST-fusion protein using baculovirus expression system. GST-TSSK1B was purified by using glutathione sepharose chromatography. GST-TSSK1B was cleaved by PreScission protease and GST-free TSSK1B (42 kDa) was collected as flow-through fraction from glutathione sepharose chromatography.

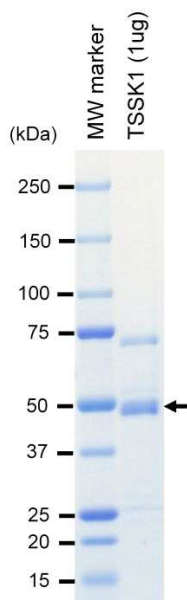
Storage buffer:

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.

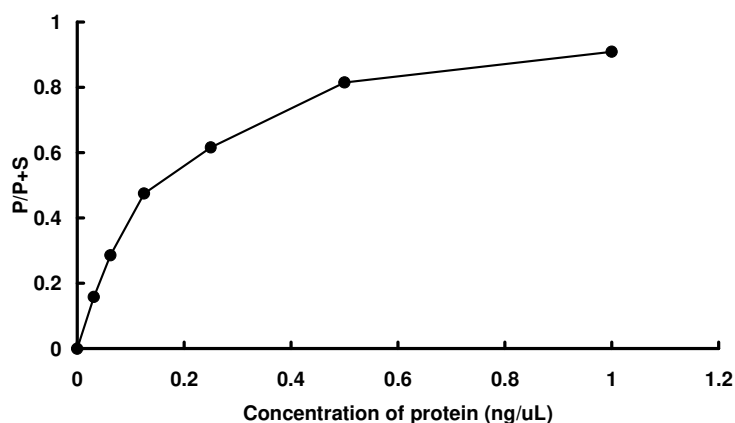
SDS-PAGE



Purity: 67 %

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 : GS peptide

 ATP : 100 μ M