

## Product Information

### YSK1(STK25)

Product Number : **07-136**

#### Product description

Full-length human YSK1 [1-426(end) amino acids of accession number NP\_006365.2] was expressed as N-terminal GST-fusion protein (75 kDa) using baculovirus expression system. GST-YSK1 was purified by using glutathione sepharose chromatography.

#### Storage buffer:

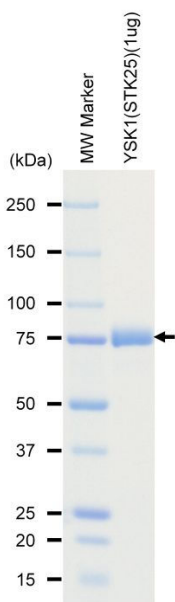
50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,  
1 mM DTT, 10% glycerol, pH7.5

#### Storage and Handling:

Store at -80C.

Avoid repeating freeze-thaws.

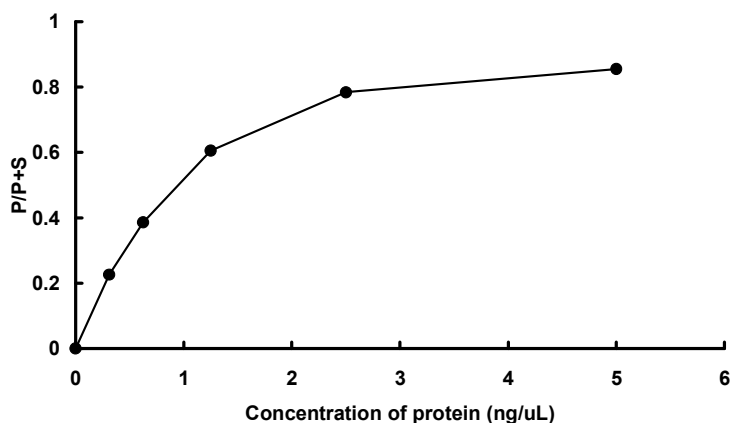
#### SDS-PAGE



Purity: 98 %

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 : Moesin-derived peptide

ATP : 100 µM