

## Product Information

### STK33

Product Number : **02-301**

#### Product description

Full-length human STK33 [1-514(end) amino acids of accession number NP\_112168.1] was expressed as N-terminal GST-fusion protein using baculovirus expression system. GST-STK33 was purified by using glutathione sepharose chromatography. GST-STK33 was cleaved by PreScission protease and GST-free STK33 (59 kDa) was collected as flow-through fraction from glutathione sepharose chromatography.

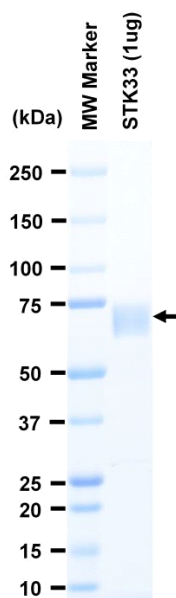
#### Storage buffer:

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.

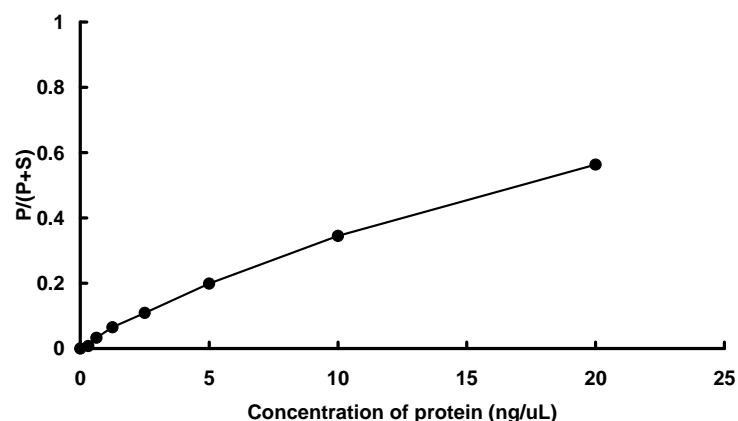
#### SDS-PAGE



Purity: 96 %

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  $\mu$ M