

## Product Information

### SLK

 Product Number : **07-129**

#### Product description

Full-length human SLK [1-1152(end) amino acids and S5N of accession number NP\_055535.1] was expressed as N-terminal GST-fusion protein (160 kDa) using baculovirus expression system. GST-SLK was purified by using glutathione sepharose chromatography.

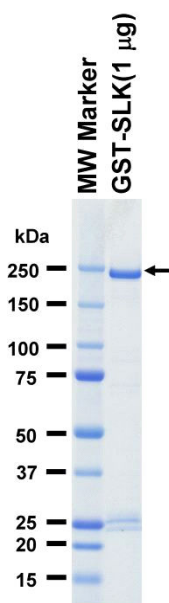
#### Storage buffer:

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

#### Storage and Handling:

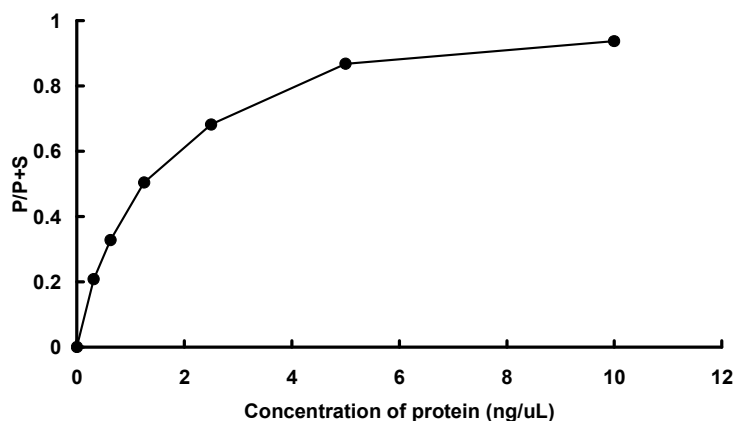
Store at -80C.  
 Avoid repeating freeze-thaws.

#### SDS-PAGE



Purity: 68 %  
 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  $\mu$ M