

Product Information

SGK

Product Number : **01-158**

Product description

Truncated human SGK [61-431(end) amino acids and S422D of accession number NP_005618.2] was co-expressed as N-terminal GST-fusion protein (69 kDa) with His-tagged PDK1[1-556(end) amino acids of accession number NP_002604.1] using baculovirus expression system. GST-SGK was purified by using glutathione sepharose chromatography.

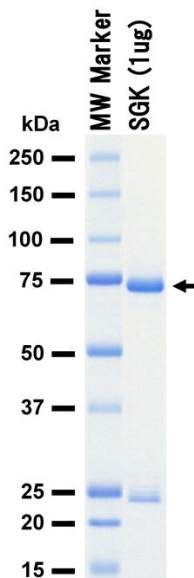
Storage buffer:

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

Storage and Handling:

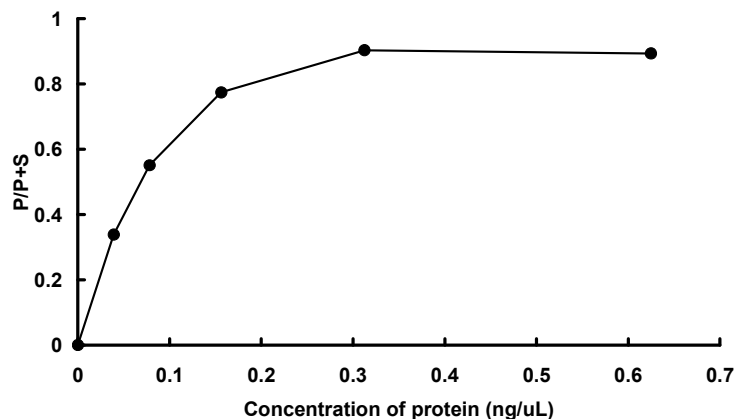
Store at -80C.
 Avoid repeating freeze-thaws.

SDS-PAGE



Purity: 73 %
 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 : SGKtide

ATP : 100 μ M