

Product Information

SGK2

Product Number : **01-159**

Product description

Full-length human SGK2 [1-367(end) amino acids and S356D of accession number NP_733794.1] was co-expressed as N-terminal GST-fusion protein (68 kDa) with His-tagged PDK1 [1-556(end) amino acids of accession number NP_002604.1] using baculovirus expression system. GST-SGK2 was purified by using glutathione sepharose chromatography.

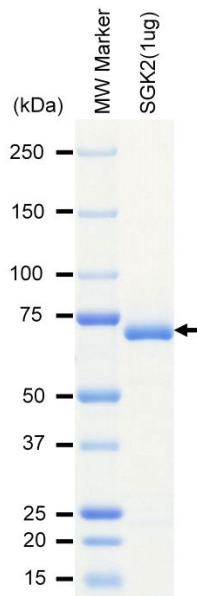
Storage buffer:

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

Storage and Handling:

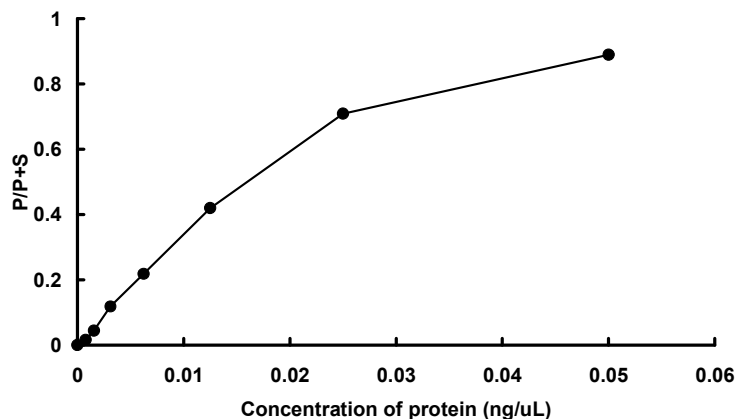
Store at -80C.
 Avoid repeating freeze-thaws.

SDS-PAGE



Purity: 97 %
 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