

Product Information

SRPK1

 Product Number : **04-160**

Product description

Full-length human SRPK1 [1-655(end) amino acids and V504 deletion of accession number NP_003128.3] was expressed as N-terminal GST-fusion protein (101 kDa) using E. coli expression system. GST-SRPK1 was purified by using glutathione sepharose chromatography.

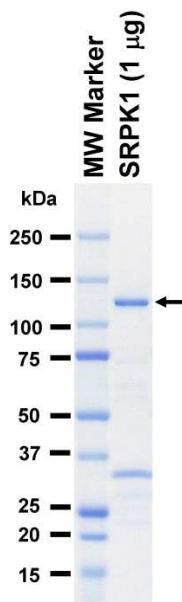
Storage buffer:

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

Storage and Handling:

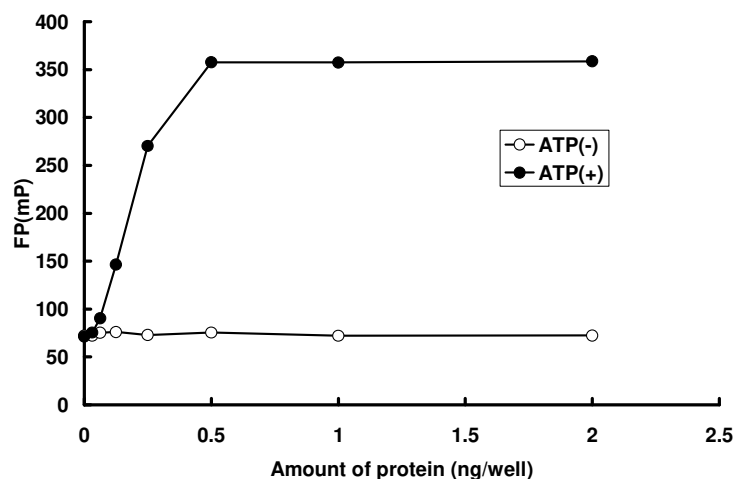
Store at -80C.
 Avoid repeating freeze-thaws.

SDS-PAGE



Purity: 56 %
 The purity was assessed by SDS-PAGE/CBB staining.

Activity data



The activity was determined by IMAP™ assay. The enzyme was incubated with fluorescein labeled peptide and phosphorylation was detected by IMAP™ technology (fluorescence polarization).

Substrate : SRPKtide

ATP : 100 µM