

SGK3, Unactive

Full-length recombinant protein expressed in Sf9 cells

Catalog # S08-14G

Lot #D073-1

Product Description

Recombinant full-length human SGK3 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_013257](#).

Gene Aliases

CISK, SGKL

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 0.25mM DTT, 0.1mM EGTA, 0.1mM EDTA, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

Scientific Background

SGK3 is a member of the SGK family and encodes a phosphoprotein with a PX (phox homology) domain. PKD1 can phosphorylate and activate SGK3 in vitro (1). A 10-fold increase in PKD1 increases the phosphorylation status of SGK3. When expressed in human embryonic kidney cells, IGF1 and peroxide significantly activate SGK3, and the activation could be reduced by preincubation with inhibitors of PI3 kinase. A yeast 2-hybrid screen found direct interaction between human SGK3 and GSK3 β . SGK3 phosphorylates several target proteins and has a role in neutral amino acid transport and activation of potassium and chloride channels (2).

References

1. Kobayashi, T. et al: Characterization of the structure and regulation of two novel isoforms of serum- and glucocorticoid-induced protein kinase. *Biochem. J.* 344: 189-197, 1999.
2. Gamper, N. et al: K⁺ channel activation by all three isoforms of serum- and glucocorticoid-dependent protein kinase SGK. *Europ. J. Physiol.* 445: 60-66, 2002.

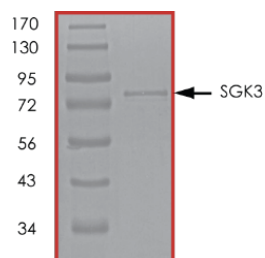
To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com

www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.

THE PRODUCT IS PROTECTED BY THE FOLLOWING PATENT/ APPLICATION: US 7, 083, 511

Purity



The purity of SGK3 was determined to be **>90%** by densitometry, approx. MW **82kDa**.

SGK3, Unactive

Full-length recombinant protein expressed in Sf9 cells

Catalog Number

S08-14G

Specific Lot Number

D073-1

Purity

>90%

Concentration

0.2 $\mu\text{g}/\mu\text{l}$

Stability

1yr at -70°C from date of shipment

Storage & Shipping

Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.