

SGK1, Unactive

Full-length recombinant human protein expressed in Sf9 cells

Catalog # S06-14G

Lot # B2009-7

Product Description

Full-length recombinant human SGK1 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_005627](#).

Gene Aliases

SGK

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 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.

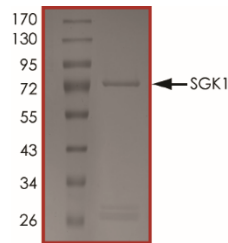
Scientific Background

SGK1 is a member of the serum- and glucocorticoid-induced protein kinase that is activated *in vitro* by 3-phosphoinositide-dependent protein kinase-1 (PDK1) and *in vivo* in response to signals that activate phosphatidylinositol (PI) 3-kinase (1). SGK1 mRNA is expressed in all tissues and the level of SGK1 mRNA is increased by stimulation with serum or dexamethasone. SGK1 promotes cell survival by phosphorylating and inactivating FKHRL1 (2). SGK and AKT display differences with respect to the efficacy with which they phosphorylate the three regulatory sites on FKHRL1.

References

1. Kobayashii, T. et al: Characterization of the structure and regulation of two novel isoforms of serum- and glucocorticoid-induced protein kinase. *Biochem J.* 1999 Nov 15;344 Pt 1:189-97.
2. Brunet, A. et al: Protein kinase SGK mediates survival signals by phosphorylating the forkhead transcription factor FKHRL1 (FOXO3a). *Mol Cell Biol.* 2001 Feb;21(3):952-65.

Purity



The purity was determined to be **>85%** by densitometry. Approx. MW **75kDa**.

SGK1, Unactive

Recombinant protein expressed in Sf9 cells

Catalog Number	S06-14G
Lot #	B2009-7
Purity	>85%
Concentration	0.05µg/µl
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.

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.