

SOK1, Unactive

Full-length recombinant protein expressed in Sf9 cells

Catalog # S43-14G

Lot # M298-1

Product Description

Recombinant full-length human SOK1 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_006374](#).

Gene Aliases

STK25, YSK1, DKFZp686J1430

Formulation

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

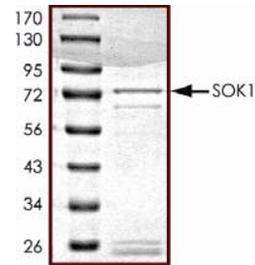
Scientific Background

SOK1 or oxidant stress-response kinase-1 is a serine/threonine kinase that is activated by autophosphorylation and by reactive oxygen intermediates, but is not activated by growth factors, alkylating agents, cytokines, or environmental stresses (1). SOK1 is expressed ubiquitously with highest expression in brain and testis. SOK1 regulates cell death after chemical anoxia, as its down-regulation by RNA interference enhances cell survival. SOK1 entry into the nucleus is important for the cell death response since SOK1 mutants unable to enter the nucleus do not induce cell death (2).

References

1. Pombo, C M. et al: Activation of a human Ste20-like kinase by oxidant stress defines a novel stress response pathway. EMBO J. 15: 4537-4546, 1996.
2. Nogueira, E. et al: SOK1 translocates from the Golgi to the nucleus upon chemical anoxia and induces apoptotic cell death. J Biol Chem. 2008 Jun 6; 283(23):16248-58.

Purity



The purity of SOK1 was determined to be >75% by densitometry. Approx. MW 74kDa.

SOK1, Unactive

Full-length recombinant protein expressed in Sf9 cells

Catalog Number S43-14G

Specific Lot Number M298-1

Purity	>75%
Concentration	0.2 µg/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.