

Anti-SGK1

Rabbit Polyclonal Antibody

Catalog # S06-63R

Lot # O2121-45

Cited Applications

WB, ELISA, IF, IHC

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the SGK1 protein

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

SGK1 antibody was raised against an 18 amino acid synthetic peptide near the carboxy terminus of human SGK1

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

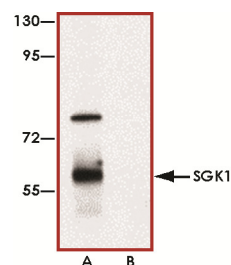
Scientific Background

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 FKHL1 (2). SGK and Akt display differences with respect to the efficacy with which they phosphorylate the three regulatory sites on FKHL1.

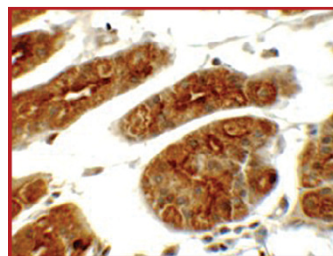
References

1. Horike N, et al: Adipose-specific expression, phosphorylation of ser794 in insulin receptor substrate-1, and activation in diabetic animals of salt-inducible kinase-2. *J. Biol. Chem.* 278: 18440-18447, 2003.
2. Dentin R, et al: Insulin modulates gluconeogenesis by inhibition of the coactivator TORC2. *Nature* 449: 366-369, 2007.

Sample Data



Western blot analysis of SGK1 in human stomach tissue lysate with SGK1 antibody at (A) 1 and (B) 2 µg/ml.



Immunohistochemistry of SGK1 in human stomach tissue with SGK1 antibody at 5 µg/ml.

Anti-SGK1

Rabbit Polyclonal Antibody

Catalog Number

S06-63R

Specific Lot Number

O2121-45

Purification

Affinity chromatography

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

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.