

Anti-SIK1

Rabbit Polyclonal Antibody

Catalog # S14-63R

Lot # O2121-43

Cited Applications

WB, ELISA, IHC

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

Specificity

Recognizes the SIK1 protein

Cross Reactivity

Human

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

SIK1 antibody was raised against an 18 amino acid synthetic peptide near the center of human SIK1

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

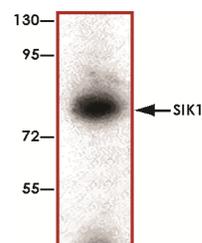
Scientific Background

SIK is a protein kinase that is involved in regulating AMPK-related kinases (1). SIK may mediate the physiological effects of LKB1, including its tumour suppressor function. SIK is also involved in signaling by various proteins like STRAD, NUA1, NUA2, BRSK1, BRSK2, QIK, QSK, SIK, MARK1, MARK2, MARK3, MARK4 and MELK that are related to AMPK. Activation of SIK1 by phosphorylation on thr322 can lead to an increase in the catalytic activity of sodium/potassium ATPase alpha subunit at the plasma membrane (2). This results in an increase in intracellular sodium in intact mammalian cells.

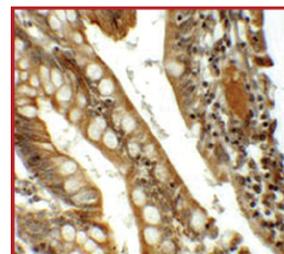
References

1. Takemori, H. Et al: TORC-SIK cascade regulates CREB activity through the basic leucine zipper domain. FEBS J. 2007; 274 (13): 3202-9.
2. Sjoström, M.; SIK1 is part of a cell sodium-sensing network that regulates active sodium transport through a calcium-dependent process. Proc. Nat. Acad. Sci. 104: 16922-16927, 2007.

Sample Data



Western blot analysis of SIK1 in human small intestine tissue lysate with SIK1 antibody at 1 µg/ml.



Immunohistochemistry of SIK1 in human small intestine tissue with SIK1 antibody at 5 µg/ml.

Anti-SIK1

Rabbit Polyclonal Antibody

Catalog Number

S14-63R

Specific Lot Number

O2121-43

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.