

## Anti-STK39 (STLK3)

Rabbit Polyclonal Antibody

**Catalog # S51-63R**

Lot # O2121-16

### Cited Applications

WB, ELISA, IF, IHC

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

### Specificity

Recognizes the STK39 (STLK3) protein

### Cross Reactivity

Human and Mouse

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

STK39 antibody was raised against a 19 amino acid synthetic peptide from near the carboxy terminus of human STK39

### Formulation

PBS + 0.02% sodium azide

### Stability

1yr at -20°C from date of shipment

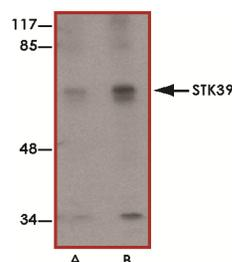
### Scientific Background

STK39 (also known as serine threonine kinase 39) is involved in the cellular stress response pathway. STK39 is activated in response to hypotonic stress leading to phosphorylation of several cation-chloride-coupled co-transporters. STK39 activates the p38 MAP kinase pathway and its interaction with p38 decreases during cellular stress (1). STK39 acts as an intermediate in the response to cellular stress. STK39 is also an independent risk factor for hypertension in men and its intragenic SNPs can interact and function in the control of blood pressure (2).

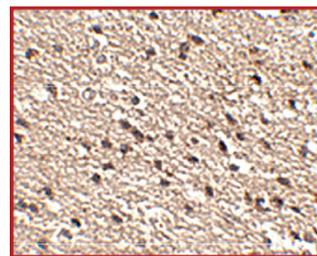
### References

1. Johnston, A. M. et.al: SPAK, a STE20/SPS1-related kinase that activates the p38 pathway. *Oncogene* 19: 4290-4297, 2000.
2. Chen LY. et.al: STK39 is an independent risk factor for male hypertension in Han Chinese. *Int J Cardiol*, 2010 Oct 1.

### Sample Data



Western blot analysis of STK39 in SK-N-SH cell lysate with STK39 antibody at (A) 1 and (B) 2 ug/ml.



Immunohistochemistry of STK39 in human brain tissue with STK39 antibody at 2.5 ug/ml.

## Anti-STK39 (STLK3)

Rabbit Polyclonal Antibody

Catalog Number

S51-63R

Specific Lot Number

O2121-16

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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

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