

## Anti-BRSK2

Rabbit Polyclonal Antibody

**Catalog # B14-63R**

Lot # O2121-5

### Cited Applications

WB, ELISA, IF, IHC

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

### Specificity

Recognizes the BRSK1 protein

### Cross Reactivity

Human and Mouse

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

BRSK2 antibody was raised against a 23 amino acid synthetic peptide from near the carboxy terminus of human BRSK2.

### Formulation

PBS + 0.02% sodium azide

### Stability

1yr at -20°C from date of shipment

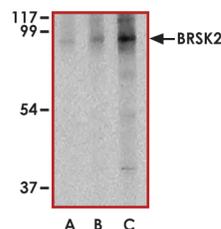
### Scientific Background

BRSK2 is a brain-selective serine/threonine kinase 2 that is mainly expressed in the brain, with weaker expression in testis and pancreas. BRSK2 expressed in insect cells specifically phosphorylates WEE1A, CDC25C and CDC25B in an in vitro assay. DNA damage induced by ultraviolet (UV) irradiation or methyl methane sulfonate, but not by ionizing radiation, enhanced endogenous BRSK2 kinase activity in a caffeine-sensitive manner and caused translocation of BRSK2 from the cytoplasm to the nucleus (1). Overexpression of BRSK2 induces G2/M arrest in HeLa cells while small interfering RNA against BRSK2 partly abrogated UV-induced G2/M arrest. Mammalian BRSK2 kinases are required for neuronal polarization (2).

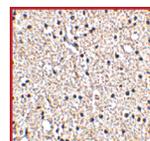
### References

1. Lu, R. et.al: Human SAD1 kinase is involved in UV-induced DNA damage checkpoint function. J. Biol. Chem. 279: 31164-31170, 2004.
2. Kishi, M. et.al: Mammalian SAD kinases are required for neuronal polarization. Science 307: 929-932, 2005.

### Sample Data



Western blot analysis of BRSK2 in human brain tissue lysate with BRSK2 antibody at (A) 0.5, (B) 1 and (C) 2 ug/ml.



Immunohistochemistry of BRSK2 in human brain tissue with BRSK2 antibody at 5 ug/ml.

## Anti-BRSK2

Rabbit Polyclonal Antibody

Catalog Number

B14-63R

Specific Lot Number

O2121-5

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.**