

Anti-BRSK1

Rabbit Polyclonal Antibody

Catalog # B13-63BR

Lot # O2121-3

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, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

BRSK1 antibody was raised against a 28 amino acid synthetic peptide from near the center of human BRSK1

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

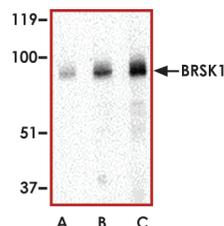
Scientific Background

BRSK1 is serine/threonine kinases 1 which required for presynaptic differentiation in *Caenorhabditis elegans* that are needed for neuronal polarization (1). BRSK1 is highly expressed in all specific adult brain regions followed by fetal brain and adult spinal cord. It is also expressed in adult heart, pancreas, testis, ovary, lung, and kidney, and in fetal liver (2).

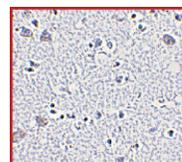
References

1. Kishi, M. et.al: Mammalian SAD kinases are required for neuronal polarization. *Science* 307: 929-932, 2005.
2. Nagase, T.et.al: Prediction of the coding sequences of unidentified human genes. XX. The complete sequences of 100 new cDNA clones from brain which code for large proteins in vitro. *DNA Res.* 8: 85-95, 2001.

Sample Data



Western blot analysis of BRSK1 in human brain tissue lysate with BRSK1 antibody at (A) 0.5, (B) 1 and (C) 2 ug/ml.



Immunohistochemistry of BRSK1 in human brain tissue with BRSK1 antibody at 5 ug/ml.

Anti-BRSK1

Rabbit Polyclonal Antibody

Catalog Number

B13-63BR

Specific Lot Number

O2121-3

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.