

Anti-CASK

Rabbit Polyclonal Antibody

Catalog # C19-63R

Lot # O2121-47

Cited Applications

WB, ELISA, IF

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

Specificity

Recognizes the CASK protein

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

CASK antibody was raised against a 16 amino acid synthetic peptide near the amino terminus of human CASK

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

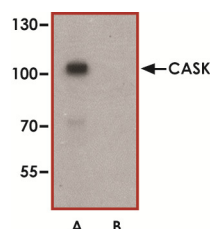
Scientific Background

CASK is a calcium/calmodulin-dependent serine protein kinase which is a MAGUK (membrane-associated guanylate kinase) protein family member (1). CASK is a scaffold protein and the encoded protein is located at synapses in the brain. CASK associates with FG syndrome 4, mental retardation, microcephaly with pontine and cerebellar hypoplasia, and a form of X-linked mental retardation. CASK functions as a cytoskeletal membrane scaffold that coordinates signal transduction pathways within the cortical cytoskeleton (2).

References

1. Atasoy, D.et.al: Deletion of CASK in mice is lethal and impairs synaptic function. Proc. Nat. Acad. Sci. 104: 2525-2530, 2007.
2. Cohen, A. R. et.al: Human CASK/LIN-2 binds syndecan-2 and protein 4.1 and localizes to the basolateral membrane of epithelial cells. J. Cell Biol. 142: 129-138, 1998.

Sample Data



Western blot analysis of CASK in mouse brain tissue lysate with CASK antibody at 1 ug/ml in (A) the absence and (B) the presence of blocking peptide.

Anti-CASK

Rabbit Polyclonal Antibody

Catalog Number

C19-63R

Specific Lot Number

O2121-47

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.