

RCAN1 Polyclonal Antibody

(Catalog # A72532)

Background

The protein encoded by this gene interacts with calcineurin A and inhibits calcineurin-dependent signaling pathways, possibly affecting central nervous system development. This gene is located in the minimal candidate region for the Down syndrome phenotype, and is overexpressed in the brain of Down syndrome fetuses. Chronic overexpression of this gene may lead to neurofibrillary tangles such as those associated with Alzheimer disease. Alternative splicing results in multiple transcript variants.

Description

RCAN1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P53805

Purification

Affinity Purified

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 60-245 of human RCAN1 (NP_004405.3).

Storage

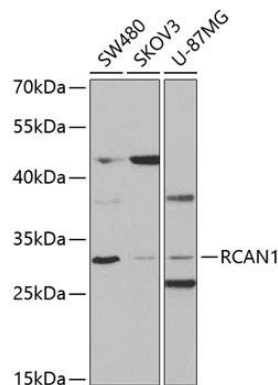
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

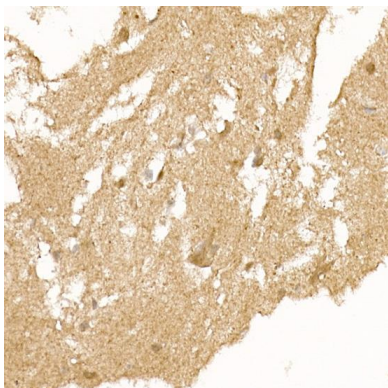
RCAN1; ADAPT78; CSP1; DSC1; DSCR1; MCIP1; RCN1; calcipressin-1

Application

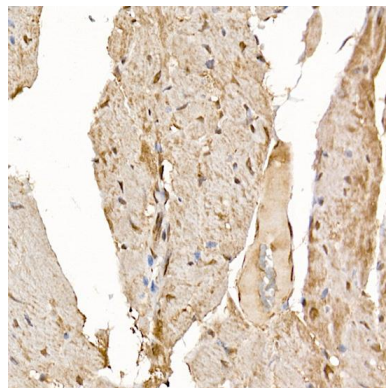
WB, IHC; Recommended dilution: WB 1:500 - 1:2000, IHC 1:50 - 1:200



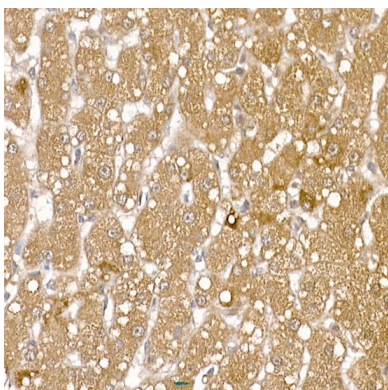
Western blot analysis of extracts of various cell lines, using RCAN1 Polyclonal Antibody at 1:1000 dilution. Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25ug per lane. Blocking buffer: 3% nonfat dry milk in TBST. Exposure time: 90s.



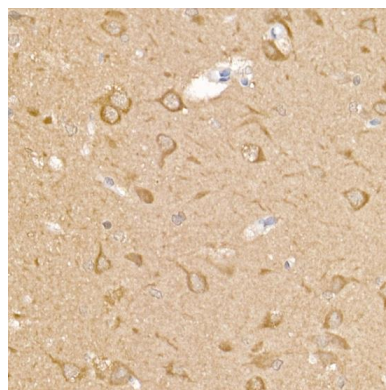
Immunohistochemistry of paraffin-embedded human brain using RCAN1 Polyclonal Antibody at dilution of 1:150 (40x lens). Performed high pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.



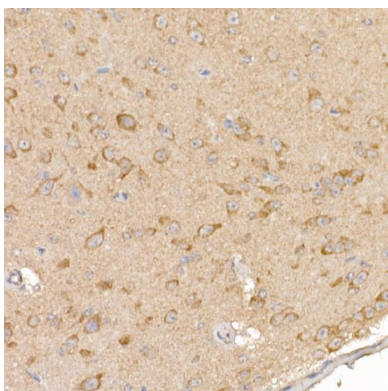
Immunohistochemistry of paraffin-embedded mouse heart using RCAN1 Polyclonal Antibody at dilution of 1:150 (40x lens). Performed high pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.



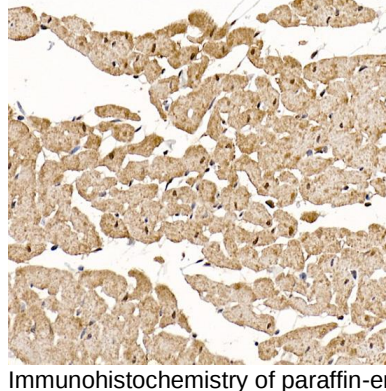
Immunohistochemistry of paraffin-embedded human liver using RCAN1 Polyclonal Antibody at dilution of 1:150 (40x lens). Performed high pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.



Immunohistochemistry of paraffin-embedded rat brain using RCAN1 Polyclonal Antibody at dilution of 1:150 (40x lens). Performed high pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.



Immunohistochemistry of paraffin-embedded mouse brain using RCAN1 Polyclonal Antibody at dilution of 1:150 (40x lens). Performed high pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.



Immunohistochemistry of paraffin-embedded rat heart using RCAN1 Polyclonal Antibody at dilution of 1:150 (40x lens). Performed high pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.