

Anti-SIRT2

Rabbit Polyclonal Antibody

Catalog # **S36-463BR**

Lot # O2121-12

Cited Applications

WB, ELISA, IF, IHC

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

Specificity

Recognizes the SIRT2 protein

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

SIRT2 antibody was raised against a 17 amino acid synthetic peptide near the amino terminus of the human SIRT2

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

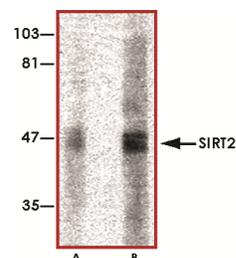
Scientific Background

SIRT2 is a member of the sirtuin family of proteins which are homologs to the yeast Sir2 protein. Sirtuin family contain a sirtuin core domain and are grouped into four classes with SIRT2 being a member of class I. Inhibition of SIRT2 can lead to neuroprotection in cellular and invertebrate models of Huntington's disease (1). Huntington's disease is characterized by increased sterol synthesis in neuronal cells and this process is reversed by SIRT2 inhibition. SIRT2 can deacetylate lys40 of alpha-tubulin both in vitro and in vivo (2). Knockdown of SIRT2 via small interfering RNA results in tubulin hyperacetylation.

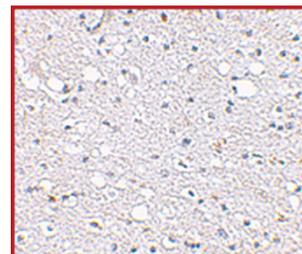
References

1. Luthi-Carter, R. Et al: SIRT2 inhibition achieves neuroprotection by decreasing sterol biosynthesis. Proc Natl Acad Sci U S A. 2010 Apr 27;107(17):7927-32.
2. North, B. J. et al: The human Sir2 ortholog, SIRT2, is an NAD(+)-dependent tubulin deacetylase. Molec. Cell 11: 437-444, 2003.

Sample Data



Western blot analysis of SIRT2 in human brain lysate with SIRT2 antibody at (A) 2.5 and (B) 5 ug/ml.



Immunohistochemical staining of human brain tissue using SIRT2 antibody at 2.5 ug/ml.

Anti-SIRT2

Rabbit Polyclonal Antibody

Catalog Number

S36-463BR

Specific Lot Number

O2121-12

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.