

SIRT2 antibody (pAb)

Catalog Nos: 39903, 39904

RRID: AB_2615001

Isotype: IgG

Application(s): IP, WB

Reactivity: Human

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 40 & 45 kDa

Background: SIRT2 is a member of the SIR2 / Sirtuin family of NAD-dependant deacetylase enzymes. It has multiple targets for deacetylation, including tubulin, P53, HOXA10, FOXO3A, P300, FOXO1. SIRT2 suppresses adipogenesis by deacetylating FOXO1, promoting the binding of FOXO1 to PPAR γ and subsequent repression of PPAR γ mediated transcriptional activation. Similar to SIRT1, SIRT2 also deacetylates H4AcK16 similar to SIRT1, SIRT2 deacetylates histone H4 at lysine 16. SIRT2 is present in the cytoplasm during most of the cell cycle except in the G2/M transition and in mitosis, where it is found in the nucleus and in contact with chromatin and required for exit from mitosis.

Immunogen: This SIRT2 antibody was raised against a synthetic peptide derived from the C-terminal region of human SIRT2.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

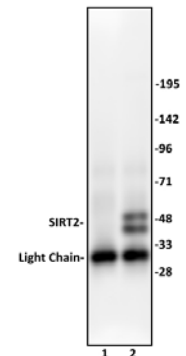
IP: 10 μ l per IP

WB: 1:500 - 1:2,000 dilution

The addition of 0.05% Tween 20 in the blocking buffer and primary antibody incubation buffer is recommended to aid in detection by Western blot. Individual optimization may be required.

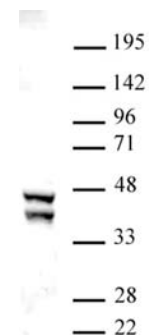
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



SIRT2 antibody (pAb) tested by Immunoprecipitation.

10 μ l of SIRT2 antibody was used to immunoprecipitate SIRT2 from 250 μ g of K-562 cytoplasmic cell extract (lane 2). 10 μ l of rabbit IgG was used as a negative control (lane 1). The immunoprecipitated protein was detected by Western blotting using the SIRT2 antibody at a dilution of 1:500.



SIRT2 antibody (pAb) tested by Western blot.

Cytoplasmic extract of K-562 cells (25 μ g per lane) probed with SIRT2 antibody (1:500).