

SIRT6 antibody (pAb)

Catalog Nos: 39911, 39912

RRID: AB_2793392

Isotype: IgG

Application(s): IP, WB

Reactivity: Human

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Concentration: 0.36 μ g/ μ l

Molecular Weight: 46 kDa

Background: SIRT6 (Sir2 homologue 6) is an NAD⁺ dependant histone deacetylase (HDAC) related to SIRT1 and to the yeast Sir2 protein. SIRT6, originally identified as a tubulin deacetylase, it was recently found to be able to deacetylate histone H3 acetylated at lysine 9.

Immunogen: This SIRT6 antibody was raised against a peptide within the C-terminus of human SIRT6.

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:

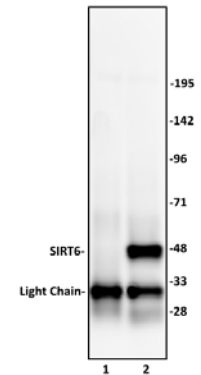
IP: 10 μ l per IP

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

*Note: many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western Blot.

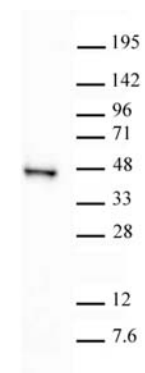
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product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

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



SIRT6 antibody (pAb) tested by Immunoprecipitation.

10 μ l of SIRT6 antibody was used to immunoprecipitate SIRT6 from 500 μ g of HeLa nuclear 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 SIRT6 antibody at a dilution of 1:1,000.



SIRT6 antibody (pAb) tested by Western blot.

HeLa cells nuclear extract (25 μ g) probed with SIRT6 antibody (1:1,000).