

SIRT2 antibody (mAb)

Catalog Nos: 61365, 61366

RRID: AB_2793607

Clone: 7G5

Isotype: IgG2a

Application(s): WB

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Rat

Concentration: 1 µg/µl

Molecular Weight: 40 & 43 kDa

Background: SIRT2 (sirtuin 2) 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 antibody was raised against a synthetic peptide corresponding to human SIRT2 surrounding serine 331. This antibody recognizes both the unphosphorylated and the Ser-331 phosphorylated form of SIRT2.

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

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.