

SIRT1 antibody (mAb)

Catalog Nos: 39353, 39354

RRID: AB_2615002

Clone: 2G1/F7

Isotype: IgG1

Application(s): ICC, IF, IP, WB

Reactivity: Human

Quantities: 200 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 120 kDa

Background: SIRT1 (Sirtuin 1, SIR2L1) is a member of the Sirtuin family of proteins related to the yeast Sir2 protein. Like Sir2 and other Sirtuins, SIRT1 is an NAD-dependent histone and protein deacetylase. SIRT1 catalyzes the removal of acetyl groups from a variety of protein substrates (NBS1, p53, Ku70, *etc.*) as well as from Histone H4 Lys16. SIRT1 is involved in the regulation of a number of cellular processes, including transcription, metabolism, DNA repair and aging, and increased SIRT1 is associated with extended lifespan of *C. elegans*.

Immunogen: This SIRT1 antibody was raised against a synthetic peptide corresponding to amino acids 581-630 of human SIRT1.

Buffer: 10 mM sodium phosphate pH 7.5, 150 mM NaCl, 30% glycerol, 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

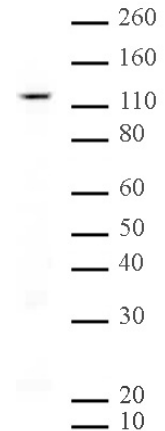
Applications Validated by Active Motif:

WB: 0.5 - 2.0 µg/ml 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.



SIRT1 mAb (Clone 2G1/F7) tested by Western blot.

Nuclear extract (20 µg) of CMSL0 cells stained with SIRT1 mAb (Clone 2G1/F7) at a dilution of 1 µg/ml.