

SIRT2 (CT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat SIRT2 (CT)

Instruction Manual

Catalog Number	PK-AB718-4485
Synonyms	SIRT2 Antibody: NAD-dependent deacetylase sirtuin-2, SIR2L, SIR2L2, SIR2-like
Description	The founding member of the sirtuin protein family was the silent information regulator 2 protein (Sir2p) of <i>Saccharomyces cerevisiae</i> , an NAD ⁺ -dependent histone deacetylase (HDAC) that regulates chromatin silencing. The SIR2 family of genes are highly conserved from prokaryotes to eukaryotes. Mammals have seven homologs of Sir2p, SIRT1-7, which are involved in diverse processes ranging from transcriptional regulation, cell cycle progression and DNA-damage repair to aging. SIRT2 is a predominantly cytoplasmic protein that colocalizes with microtubules and can deacetylate α -tubulin and regulate progression through the cell cycle. Most Sirtuins are highly expressed in brain and testis, while Sirt2 expression is higher in fetal relative to adult brain. Recent studies on SIRT2 support the therapeutic utility of inhibitors for the treatment of neurodegenerative diseases such as Parkinson's disease.
Quantity	100 μ g
Source / Host	Rabbit
Immunogen	SIRT2 antibody was raised in rabbits against a 19 amino acid peptide near the carboxy terminus of the human SIRT2.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	
Formulation	Antibody is supplied in PBS containing 0.02% sodium azide.
Reconstitution	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 μ l or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
Storage & Stability	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB, IHC, IF INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. SIRT2 antibody can be used for detection of SIRT2 by Western blot at 1 - 2 μ g/mL. Antibody can also be used for immunohistochemistry starting at 2.5 μ g/mL. For immunofluorescence start at 5 μ g/mL.
Images	Available upon request.
References	Yamamoto H, Schoonjans K, and Auwerx J. Sirtuin functions in health and disease. <i>Mol. Endocrinol.</i> 2007; 21:1745-55. Frye RA. Characterization of five human cDNAs with homology to the yeast SIR2 gene: SIR2-like proteins (sirtuins) metabolize NAD and may have protein ADP-ribosyltransferase activity. <i>Biochem. Biophys. Res. Commun.</i> 1999; 260:273-279. North BJ, Marshall BL, Borra MT, et. al. The human Sir2 ortholog, SIRT2, is an NAD ⁺ -dependent tubulin deacetylase. <i>Mol. Cell</i> 2003; 11:437-44. Inoue T, Hiratsuka M, Osaki M, et al. The molecular biology of mammalian SIRT proteins: SIRT2 in cell cycle regulation. <i>Cell Cycle</i> 2007; 6:1011-8.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4485P; SIRT2 Peptide Cat.No. PK-AB718-1403; Mouse Brain Tissue Lysate

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.