
SIRT3 Polyclonal Antibody

(Catalog #A73560)

Background

This gene encodes a member of the sirtuin family of proteins, homologs to the yeast Sir2 protein. Members of the sirtuin family are characterized by a sirtuin core domain and grouped into four classes. The functions of human sirtuins have not yet been determined; however, yeast sirtuin proteins are known to regulate epigenetic gene silencing and suppress recombination of rDNA. Studies suggest that the human sirtuins may function as intracellular regulatory proteins with mono-ADP-ribosyltransferase activity. The protein encoded by this gene is included in class I of the sirtuin family. Two alternatively spliced transcript variants that encode different proteins have been described for this gene.

Description

SIRT3 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.05% proclin300, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q9NTG7

Purification

Affinity Purified

Immunogen

A synthetic peptide corresponding to a sequence within amino acids 299-399 of human SIRT3 (NP_036371.1).

Storage

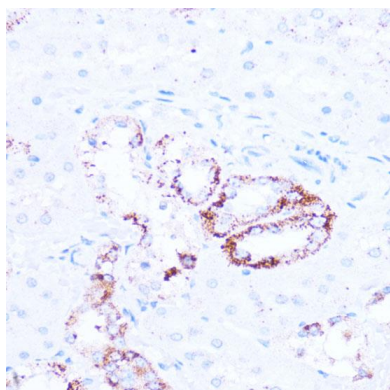
Shipped at 4°C. Store at -20°C. Avoid freeze / thaw cycles.

Alternative Names

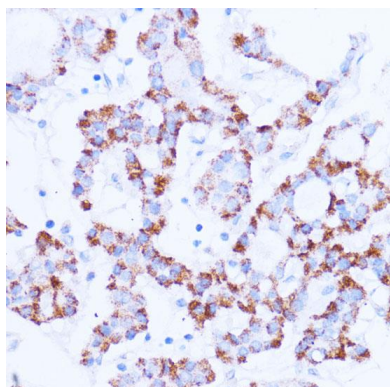
SIRT3; SIR2L3; sirtuin 3, T3

Application

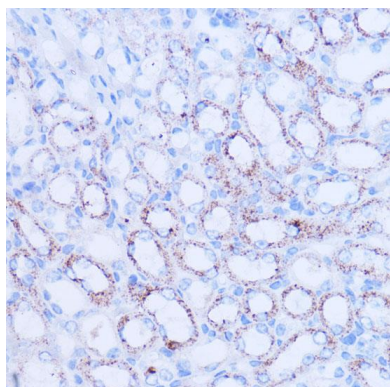
WB, IHC, IF; Recommended Dilutions: WB 1:500 - 1:1000, IHC-P 1:20 - 1:50, IF/ICC 1:50 - 1:200



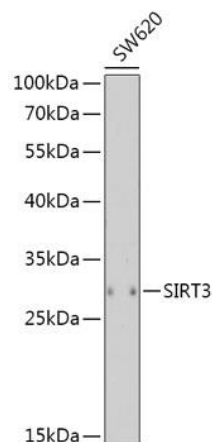
Immunohistochemistry of paraffin-embedded rat kidney using SIRT3 Polyclonal Antibody at dilution of 1:100 (40x lens).



Immunohistochemistry of paraffin-embedded human thyroid cancer using SIRT3 Polyclonal Antibody at dilution of 1:100 (40x lens).



Immunohistochemistry of paraffin-embedded mouse kidney using SIRT3 Polyclonal Antibody at dilution of 1:100 (40x lens).

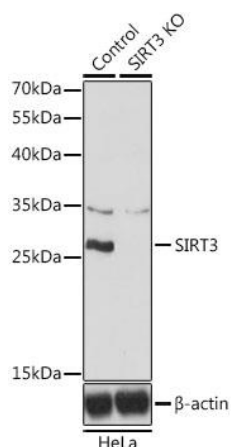


Western blot analysis of extracts of SW620 cells, using SIRT3 Polyclonal Antibody at 1:1000 dilution.

Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.

Lysates/proteins: 25ug per lane.

Blocking buffer: 3% nonfat dry milk in TBST. Exposure time: 180s.



Western blot analysis of extracts from normal (control) and SIRT3 knockout (KO) HeLa cells, using SIRT3 Polyclonal Antibody at 1:500 dilution.

Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.

Lysates/proteins: 25ug per lane.

Blocking buffer: 3% nonfat dry milk in TBST. Exposure time: 3min.