

SIRT1 (human) polyclonal antibody

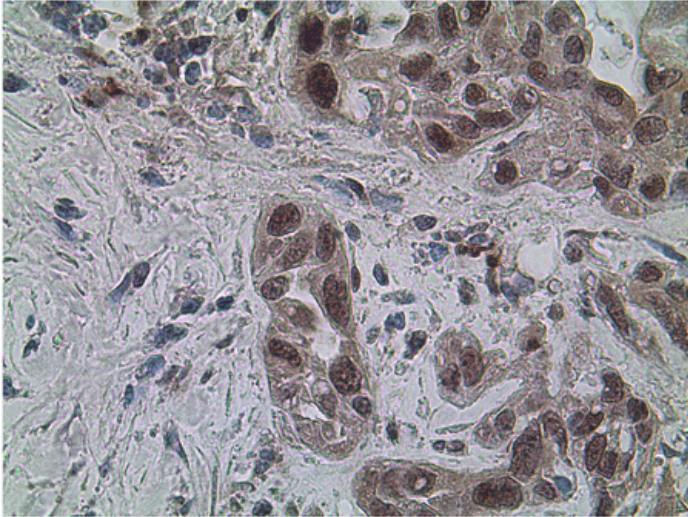
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 2

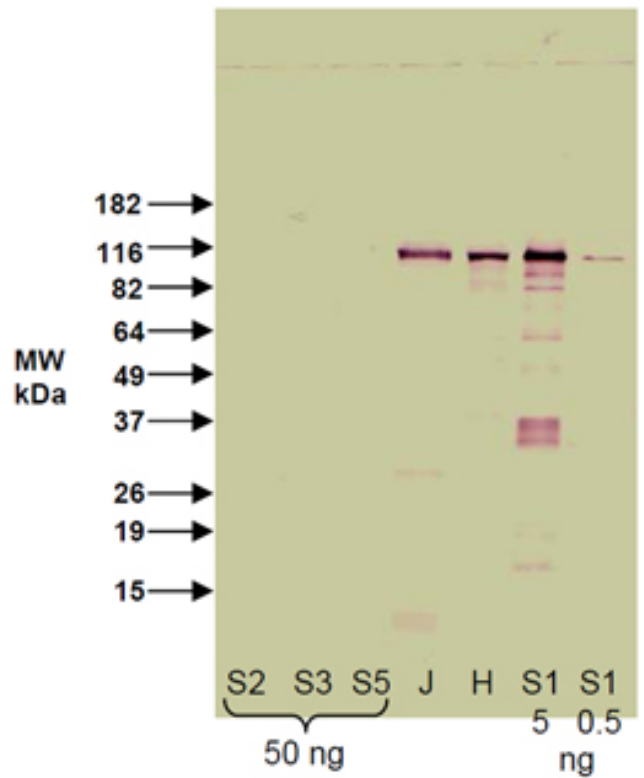
[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry staining of formalin fixed, paraffin-embedded human liver tumor tissue using SIRT1 antibody (Prod. no. BML-SA427) showing nuclear staining of hepatocytes.



Western blot analysis of recombinant human sirtuins, Jurkat cells and HeLa nuclear extract. The nitrocellulose was probed with anti- SIRT1 at 1:1000. The secondary antibody was coupled to alkaline phosphatase and the blot was developed with BCIP/NBT. Arrows indicate positions of molecular weight markers. Lanes: S2, S3, S5 (SIRTs 2, 3 & 5, 50 ng each); J (150,000 Jurkat cells); H (18 μ g HeLa nuclear extract); S1 (SIRT1- 5 ng, 0.5 ng).

Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Sirtuin-1
Application	IHC (PS), WB
Application Notes	Detects a band of ~115 kDa by Western blot. Calculated molecular weight of the full-length protein is 82 kDa.
Crossreactivity	No cross-reactivity observed with recombinant human sirtuins SIRT2, SIRT3 or SIRT5.
Formulation	Liquid. In PBS, pH 7.4.
Host	Rabbit
Immunogen	Recombinant human SIRT1 (sirtuin 1).
Purity Detail	Protein A-affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:1,000) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.
Source	Purified from rabbit serum.
Species Reactivity	Human
Specificity	Recognizes recombinant and native SIRT1.
UniProt ID	Q96EB6

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)