

SIRT5 polyclonal antibody

SIRT5 is the only human class III sirtuin and as such is the closest homolog to bacterial (e.g. *E. coli* CobB) and archaeal sirtuins. SIRT5 is an imported mitochondrial protein with a calculated precursor molecular weight of 34kDa. In various human, rat, mouse and bovine cells and tissues, the major reactive band migrates with an apparent molecular weight of ~29kDa. This presumably reflects removal of an N-terminal transit sequence due to mitochondrial import and processing. No cross-reactivity observed with recombinant human sirtuins SIRT1, SIRT2 or SIRT3.

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

Citations: 1

[View Online »](#)

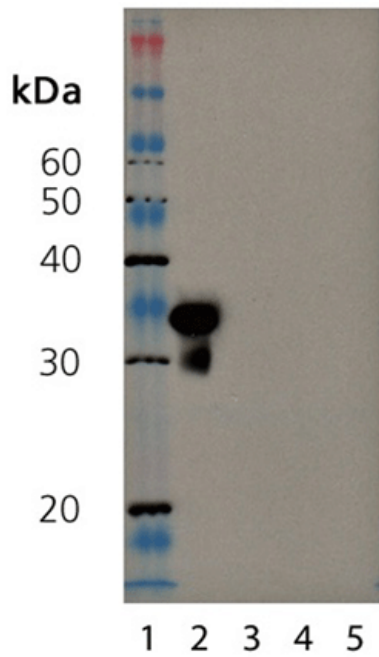
Ordering Information

[Order Online »](#)

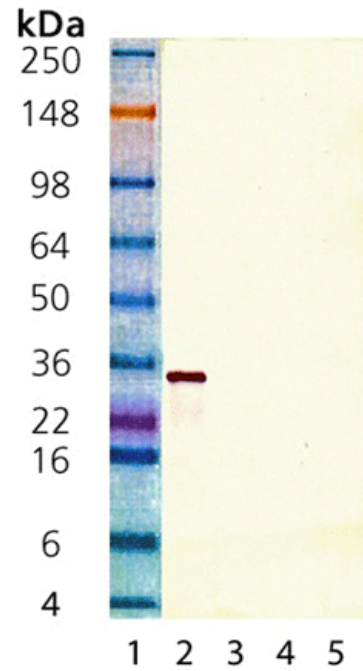
BML-SA464-0100	100µl
----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of SIRT5, pAb (Prod. No. BML-SA464): Lane 1: MW marker, Lane 2: SIRT5 (human), (recombinant) (His-tag) (Prod. No. BML-SE555), Lane 3: SIRT3 (human), (recombinant) (His-tag) (Prod. No. BML-SE270), Lane 4: SIRT2 (human), (recombinant) (His-tag) (Prod. No. BML-SE251), Lane 5: SIRT1 (human), (recombinant) (His-tag) (Prod. No. BML-SE239).



Western blot analysis of SIRT5, pAb (Prod. No. BML-SA464): Lane 1: MW marker, Lane 2: SIRT5 (human), (recombinant) (His-tag) (Prod. No. BML-SE555), Lane 3: SIRT3 (human), (recombinant) (His-tag) (Prod. No. BML-SE270), Lane 4: SIRT2 (human), (recombinant) (His-tag) (Prod. No. BML-SE251), Lane 5: SIRT1 (human), (recombinant) (His-tag) (Prod. No. BML-SE239).

Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Sirtuin-5
Application	WB
Application Notes	Detects a band of ~29 kDa by Western blot.
Formulation	Liquid. In PBS.
Host	Rabbit
Immunogen	Recombinant human SIRT5.
Purity Detail	Partially purified from rabbit serum with immunodepletion.
Recommendation Dilutions/Conditions	Western Blot: (1:2,500, Colorimetric, using secondary antibody coupled to alkaline phosphatase and BCIP/NBT as substrate. / 1:3,000, ECL, using secondary antibody coupled HRP)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	Bovine, Human, Mouse, Rat
UniProt ID	Q9NXA8
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .



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)