

SIRT3 (human), (recombinant) (His- tag)

Highly active

Sirtuins, which are NAD-dependent lysine deacetylases, have been implicated in the control of longevity in yeast and *C. elegans* and, in yeast, have been shown to be necessary for the longevity increase effected by caloric restriction. Human SIRT3, one of seven human sirtuins, has been assigned, along with SIRT1 and SIRT2, to the class I sirtuin homology group.

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE270-0500	500U
----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Sirtuin 3
Formulation	Liquid. In 25mM TRIS, pH 7.5, containing 100mM NaCl, 5mM DTT, and 10% glycerol.
MW	32.7 kDa
Purity Detail	Partially purified by single-step affinity chromatography and gel filtration.
Source	Produced in <i>E. coli</i> . Active SIRT3 (aa 102-399). Contains a N-terminal His-tag.
Specific Activity	One unit will deacetylate 1pmol/min of FLUOR DE LYS [®] -SIRT2 substrate (Prod. No. BML-KI179) at 37°C , using 500µM FLUOR DE LYS [®] -SIRT2 substrate, 500µM NAD ⁺ .
UniProt ID	Q9NTG7

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)