

SIRT2 (human), (recombinant) (His- tag)

Due to their roles in gene silencing, aging, and oxidative stress responses, it is believed that sirtuins (NAD-dependent lysine deacetylases) play a crucial role in cell metabolism. Human SIRT2 is a class I sirtuin, although not as closely related to Sir2 and human SIRT1. Its catalytic core has the same basic two-domain architecture and central groove as other sirtuins and an interaction with the homeobox transcription factor HOXA10.

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE251-0500	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 2

Formulation Liquid. In 25mM TRIS, pH 7.5, containing 100mM NaCl, 5mM DTT and 10% glycerol.

MW 43 kDa

Purity Detail Purified by multi-step chromatography.

Source Produced in *E. coli*. Contains a N-terminal His-tag.

Specific Activity One unit will deacetylate 1pmol/min of FLUOR DE LYS[®]-SIRT2 substrate (Prod. No. BML-KI179).

UniProt ID Q8IXJ6



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)