

SIRT1 (human), (recombinant) (His- tag)

Highly active

Citations: 17

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE239-0100	100U
----------------	------

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 1
Formulation	Liquid. In 25 mM TRIS, pH 7.5, 100mM NaCl, 5mM DTT and 10% glycerol.
Gene/Protein Identifier	NM_012238 (RefSeq)
MW	82 kDa
Purity Detail	Partially purified by single-step affinity chromatography and gel filtration.
Source	Produced in <i>E. coli</i> . Recombinant SIRT1 from human cDNA containing an N-terminal His-tag.
Specific Activity	One unit = 1pmol/min at 37°C, 250µM FLUOR DE LYS [®] substrate (Prod. No. BML-K1104), 500µM NAD ⁺ .
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)