

# 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 »](#)

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 <i>E. coli</i> . Contains a N-terminal His-tag.
Specific Activity	One unit will deacetylate 1pmol/min of FLUOR DE LYS <sup>®</sup> -SIRT2 substrate (Prod. No. BML-KI179).
UniProt ID	Q8IXJ6

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)