

# PRODUCT INFORMATION



## SIRT6 (human, recombinant)

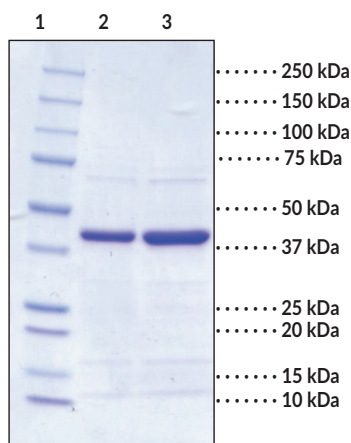
Item No. 10315

### Overview and Properties

<b>Synonyms:</b>	NAD-dependent Deacetylase 6, Silent Information Regulator 6, SIR2L6, SIR2-like Protein 6, Sirtuin 6
<b>Source:</b>	Recombinant N-terminal His-tagged enzyme expressed in <i>E. coli</i>
<b>Amino Acids:</b>	1-355
<b>Uniprot No.:</b>	Q8N6T7
<b>Molecular Weight:</b>	43.7 kDa
<b>Storage:</b>	-80°C (as supplied); avoid freeze/thaw cycles by aliquoting protein
<b>Stability:</b>	≥1.5 years
<b>Purity:</b>	≥80% estimated by SDS-PAGE
<b>Supplied in:</b>	25 mM Tris-HCl, pH 8.0, with 100mM NaCl, 20% glycerol
<b>Protein Concentration:</b>	<i>batch specific</i> mg/ml

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Image



Lane 1: MW Markers  
Lane 2: SIRT6 (2 µg)  
Lane 3: SIRT6 (4 µg)

WARNING  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 08/31/2021

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

---

The sirtuins represent a distinct class of trichostatin A-insensitive lysyl-deacetylases (class III HDACs) and have been shown to catalyze a reaction that couples lysine deacetylation to the formation of nicotinamide and O-acetyl-ADP-ribose from NAD<sup>+</sup> and the abstracted acetyl group.<sup>1-3</sup> There are seven human sirtuins, which have been designated SIRT1-7.<sup>4</sup> SIRT6 associates specifically with telomeres and functions at chromatin to decrease NF-κB signaling.<sup>5,6</sup> Mammalian cells depleted of SIRT6 display abnormal telomere structures similar to defects found in Werner syndrome, a premature ageing disorder, and have a shortened life span.<sup>5,6</sup> Since SIRT6 binds and attenuates NF-κB signaling, it is proposed that activators of SIRT6 may be effective anti-cancer and anti-inflammatory drugs and may increase longevity.<sup>5</sup>

## References

---

1. Imai, S.-I., Armstrong, C.M., Kaeberlein, M., *et al.* Transcriptional silencing and longevity protein Sir2 is an NAD-dependent histone deacetylase. *Nature* **403**, 795-800 (2000).
2. Tanner, K.G., Landry, J., Sternglanz, R., *et al.* Silent information regulator 2 family of NAD-dependent histone/protein deacetylases generates a unique product, 1-O-acetyl-ADP-ribose. *Proc. Natl. Acad. Sci. USA* **97(26)**, 14178-14182 (2000).
3. Tanny, J.C. and Moazed, D. Coupling of histone deacetylation to NAD breakdown by the yeast silencing protein Sir2: Evidence for acetyl transfer from substrate to an NAD breakdown product. *Proc. Natl. Acad. Sci. USA* **98(2)**, 415-420 (2001).
4. Frye, R.A. Phylogenetic classification of prokaryotic and eukaryotic Sir2-like proteins. *Biochem. Biophys. Res. Commun.* **273**, 793-798 (2000).
5. Kawahara, T.L.A., Michishita, E., Adler, A.S., *et al.* SIRT6 links histone H3 lysine 9 deacetylation to NF-κB-dependent gene expression and organismal life span. *Cell* **136**, 62-74 (2009).
6. Michishita, E., McCord, R.A., Berber, E., *et al.* SIRT6 is a histone H3 lysine 9 deacetylase that modulates telomeric chromatin. *Nature* **452**, 492-496 (2008).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM