

## SIRT2 Protein

Full-length recombinant protein expressed in E. coli cells

**Catalog # S36-30EG**

Lot # M310-3

### Product Description

Full-length recombinant human SIRT2 was expressed in E. coli cells using an N-terminal GST tag. The gene accession number is [NM\\_030593](#).

### Gene Aliases

SIR2L, SIR2L2, Sirtuin 2

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

### Storage and Stability

Store product at  $-70^{\circ}\text{C}$ . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

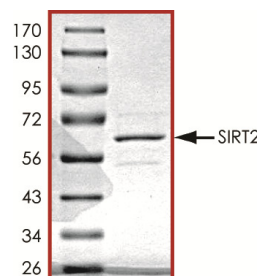
### Scientific Background

SIRT2 is a member of the sirtuin family of proteins which are homologs to the yeast Sir2 protein. Sirtuin family contain a sirtuin core domain and are grouped into four classes with SIRT2 being a member of class I. Inhibition of SIRT2 can lead to neuroprotection in cellular and invertebrate models of Huntington's disease (1). Huntington's disease is characterized by increased sterols synthesis in neuronal cells and this process is reversed by SIRT2 inhibition. SIRT2 can deacetylate lys40 of alpha-tubulin both *in vitro* and *in vivo* (2). Knockdown of SIRT2 via small interfering RNA results in tubulin hyperacetylation.

### References

1. Luthi-Carter, R. Et al: SIRT2 inhibition achieves neuroprotection by decreasing sterol biosynthesis. Proc Natl Acad Sci U S A. 2010 Apr 27;107(17):7927-32.
2. North, B. J. et al: The human Sir2 ortholog, SIRT2, is an NAD(+)-dependent tubulin deacetylase. Molec. Cell 11: 437-444, 2003.

### Purity



The purity was determined to be **>85%** by densitometry. Approx. MW **64kDa**.

## SIRT2 Protein

Full-length recombinant protein expressed in E. coli cells

Catalog Number	S36-30EG
Specific Lot Number	M310-3
Purity	>85%
Concentration	0.2µg/µl
Stability	1yr at $-70^{\circ}\text{C}$ from date of shipment
Storage & Shipping	Store product at $-70^{\circ}\text{C}$ . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**