

## SIRT6 Protein

Full-length recombinant protein expressed in E. coli cells

**Catalog # S40-30EG**

Lot # M358-3

### Product Description

Full-length recombinant human SIRT6 was expressed in E. coli cells using an N-terminal GST tag. The gene accession number is [NM\\_016539](#).

### Gene Aliases

SIR2L6, Sirtuin 6

### 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°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.

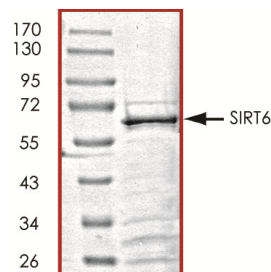
### Scientific Background

SIRT6 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 SIRT6 being a member of class IV. Human SIRT6 protein is a NAD(+)-dependent histone H3 lysine-9 deacetylase that modulates telomeric chromatin (1). SIRT6 associates specifically with telomeres and SIRT6 depletion leads to telomere dysfunction with end-to-end chromosomal fusions and premature cellular senescence. SIRT6 <sup>-/-</sup> mouse cells show that SIRT6 promotes resistance to DNA damage and suppresses genomic instability in association with a role in base excision repair (2).

### References

1. Michishita, E. et al: SIRT6 is a histone H3 lysine 9 deacetylase that modulates telomeric chromatin. Nature 452: 492-496, 2008.
2. Mostoslavsky, R. et al: Genomic instability and aging-like phenotype in the absence of mammalian SIRT6. Cell 124: 315-329, 2006.

### Purity



The purity was determined to be **>85%** by densitometry. Approx. MW **66kDa**.

## SIRT6 Protein

Full-length recombinant protein expressed in E. coli cells

Catalog Number **S40-30EG**

Specific Lot Number **M358-3**

Purity	>85%
Concentration	0.2µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°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.**