

## SIRT5 Protein

Full-length recombinant protein expressed in E. coli cells

**Catalog # S39-30EG**

Lot # M131-5

### Product Description

Full-length recombinant human SIRT5 was expressed in E. coli cells using an N-terminal GST tag. The gene accession number is [NM\\_012241](#).

### Gene Aliases

SIR2L5, Sirtuin 5

### 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

SIRT5 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 SIRT5 being a member of class III. SIRT5 consists of eight exons and is found in two isoforms which encode a 310 aa and a 299 aa protein, respectively. Human SIRT5 is most predominantly expressed in heart muscle cells and in lymphoblasts. Fluorescence in situ hybridization analysis localized the human SIRT5 gene to chromosome 6p23. SIRT5 can deacetylate cytochrome c, a protein of the mitochondrial intermembrane space with a central function in oxidative metabolism as well as apoptosis initiation (1).

### References

- Schlicker C. et al: Substrates and regulation mechanisms for the human mitochondrial sirtuins Sirt3 and Sirt5. *J Mol Biol.* 2008 Oct 10;382(3):790-801.
- Mahlknecht, U. et al: Assignment of the NAD-dependent deacetylase sirtuin 5 gene (SIRT5) to human chromosome band 6p23 by in situ hybridization. *Cytogenet Genome Res.* 2006;112(3-4):208-12.

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)

### Purity



The purity was determined to be **>95%** by densitometry. Approx. MW **60kDa**.

## SIRT5 Protein

Full-length recombinant protein expressed in E. coli cells

Catalog Number	S39-30EG
Specific Lot Number	M131-5
Purity	>95%
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.