
Sirtuin1 Protein

(Catalog No. E24015)

Description

Human Sirtuin 1 (GenBank) Accession No. NM_012238), amino acids 193- 741 with N-terminal GST tag.

Molecular Weight (kDa)

87.2 kDa

Formulation

50mM Tris pH 7.4, 250mM NaCl, 90mM (NH₄)₂SO₄, 10% glycerol, and 2mM DTT.

Source

Expressed in E. coli expression system

Size

100 µg

Storage Temp & Conditions

>6 months at -80°C

Purification Notes

Purity: >75%

Application Notes

Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Ordering Information:

Products	Size	Cat. No.
Sirtuin1 Protein	100 µg	E24015-1

This product is for research purposes only. Not intended for use in diagnostic procedures.