
Sirtuin3 Protein

(Catalog No. E24017)

Description

Human Sirtuin 3 (GenBank Accession No. NM_012239), amino acids 102- 399 with N-terminal GST tag.

Molecular Weight (kDa)

59 kDa

Formulation

45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 20 mM glutathione, 10% glycerol.

Source

Expressed in E.coli expression system

Size

100 µg

Storage Temp & Conditions

>6 months at -80°C

Purification Notes

Purity: >50%

Application Notes

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

Ordering Information:

Products	Size	Cat. No.
Sirtuin3 Protein	100 µg	E24017-1

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