
Sirtuin4 Protein

(Catalog No. E24018)

Description

Human Sirtuin 4 (GenBank Accession No.NM_012240), amino acids 25-314 with N-terminal GST tag.

Molecular Weight (kDa)

59 kDa

Formulation

40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 16 mM glutathione, and 20% glycerol.

Source

Expressed in E.coli expression system

Size

100 µg

Storage Temp & Conditions

>6 months at -80°C

Purification Notes

Purity: >90%

Application Notes

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

Ordering Information:

Products	Size	Cat. No.
Sirtuin4 Protein	100 µg	E24018-1

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