
Sirt6 Protein

(Catalog No. E24013)

Description

Human Sirtuin 6 (GenBank Accession No. NM_016539), full length with N-terminal GST tag.

Molecular Weight (kDa)

65 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 an 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.
Sirt6 Protein	100 µg	E24013-1

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