
Sirtuin 7 Protein

(Catalog No. E24014)

Description

Human Sirtuin 7 (GenBank Accession No. NM_016538), full length.

Molecular Weight (kDa)

46 kDa

Formulation

25 mM Tris-HCl, pH 8.0, 100 mM NaCl, 0.05% Tween-20 and 20% glycerol

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.
Sirtuin 7 Protein	100 µg	E24014-1

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