

---

## Sirtuin2 Protein

(Catalog No. E24016)

---

### Description

Human Sirtuin 2 (GenBank Accession No. NM\_030593), amino acids 50-356 with C-terminal His tag

### Molecular Weight (kDa)

35.5 kDa

### Formulation

40mM Tris pH 8.0, 110mM NaCl, 2.1mM KCl, and 20% glycerol.

### Source

Expressed in E.coli expression system.

### Size

100 µg

### Storage Temp & Conditions

>6 months at -80°C

### Purification Notes

Purity: >93%

### Application Notes

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

### Ordering Information:

Products	Size	Cat. No.
Sirtuin2 Protein	100 µg	E24016-1

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