

---

## Sirtuin5 Protein

(Catalog No. E24019)

---

### Description

Human Sirtuin 5 (GenBank Accession No. NM\_012241), full length with N-terminal GST tag.

### Molecular Weight (kDa)

59.8 kDa

### Formulation

45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 18 mM glutathione and 10% 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.
Sirtuin5 Protein	100 µg	E24019 -1

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