

---

## Sumo1-biotin Protein

(Catalog No. E28001)

---

### Description

SUMO-1 modified with biotin via primary amine coupling. This results in multiple biotinylated SUMO-1 species modified at the N-terminus, as well as lysine residues. Although having a fully functional C-terminus, lysine modification may limit the ability of this reagent to propagate poly-SUMO chains. Biotinylated SUMO-1 can be detected using avidin-linked reagents.

### Molecular Weight (kDa)

11.1 kDa

### Formulation

50 mM HEPES pH 8.0, 150 mM NaCl, 1mM DTT, 10% glycerol

### Source

Expressed in E.coli expression system

### Size

50 µg

### Storage Temp & Conditions

Store stock solution at -80°C. Avoid multiple freeze/thaw cycles. Best use within three months from the date of receipt of this protein.

### Purification Notes

Purity: > 95% by SDS-PAGE

### Application Notes

Enzyme-linked Immunoabsorbent Assay; Western Blot (Recombinant protein); Protein Array.

### Ordering Information:

Products	Size	Cat. No.
SUMO1-biotin Protein	50 µg	E28001 -1

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