

---

## SUMO3-biotin Protein

(Catalog No. E28005)

---

### Description

SUMO-3 modified with biotin via primary amine coupling resulting in modification of lysine residues as well as the N-terminus. Although having a fully functional C-terminus, lysine modification may limit the ability of this reagent to propagate poly-SUMO chains. Biotinylated SUMO-3 can be detected using avidin-linked reagents for higher efficiency and sensitivity than with antibodies.

### Molecular Weight (kDa)

10.5 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.

### Purification Notes

Purity: > 95% by SDS-PAGE

### Application Notes

Enzyme-linked Immunoabsorbent Assay; Western Blot (Recombinant protein; Protein Array; ;Best use within three months from the date of receipt of this protein.

### Ordering Information:

Products	Size	Cat. No.
SUMO3-biotin Protein	50 µg	E28005 -1

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