
SUMO2-biotin Protein

(Catalog No. E28003)

Description

SUMO-2 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-2 can be detected using avidin-linked reagents for higher efficiency and sensitivity than with antibodies.

Molecular Weight (kDa)

10.6 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 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.
SUMO2-biotin Protein	50 µg	E28003 -1

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