

## Biotin-PEG11-Azide

**Catalog #:** 27436

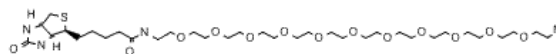
**Lot #:** 141126

**Size:** 100 mg

**CAS Registry #:** NA

**Structure:**

**Purity:** ≥96%



**Chemical Formula:** C<sub>34</sub>H<sub>64</sub>N<sub>6</sub>O<sub>13</sub>S

**Molecular Weight:** 797.0 Da

**Description:** Biotinylation reagent with Azide, and a long 11 PEG spacer. Labels terminal alkynes via a copper-catalyzed “click reaction”. Useful for labeling fluorescent dyes with triple bond alkyne groups.

**Appearance:** White to off-white solid. May appear sticky.

**Solubility:** Soluble in DMSO. Limited solubility in water. Preferably, stock solutions should be made in DMSO, then diluted into aqueous solutions.

**Application:** Useful for biotinylation of fluorescent dyes or other compounds with triple bond alkynes via click chemistry.

**Storage/Stability:** Store at or below ≤4°C. Solid form is stable at least 2 years from date of receipt, when stored as directed. *Protect from moisture, store with dessicant.*

### References:

1. Parrish, B., *et al.*, *J. Am. Chem. Soc.* 2005; **127(20)**: 7404-7410.