



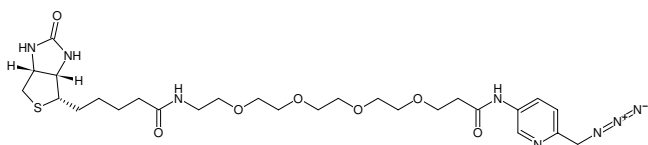
## Picolyl-Azide-PEG<sub>4</sub>-Biotin

Cat. No.	Amount
CLK-1167-5	5 mg
CLK-1167-25	5 x 5 mg
CLK-1167-100	100 mg

### Selected References:

Jiang *et al.* (2014) Monitoring Dynamic Glycosylation in Vivo Using Supersensitive Click Chemistry. *Bioconjug. Chem.* **25** (4):698.

Uttamapinant *et al.* (2012) Fast, cell-compatible click chemistry with copper-chelating azides for biomolecular labeling. *Angew. Chem. Int. Ed. Engl.* **51** (24):5852.



Structural formula of Picolyl-Azide-PEG<sub>4</sub>-Biotin

### For general laboratory use.

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at -20 °C

**Shelf Life:** 12 months after date of delivery

**Molecular Formula:** C<sub>27</sub>H<sub>42</sub>N<sub>8</sub>O<sub>7</sub>S

**Molecular Weight:** 622.74 g /mol

**Exact Mass:** 622.29 g /mol

**Purity:** ≥ 90 % (HPLC)

**Form:** amorphous solid

**Color:** off-white to slightly orange

**Solubility:** DMSO, DMF, MeOH