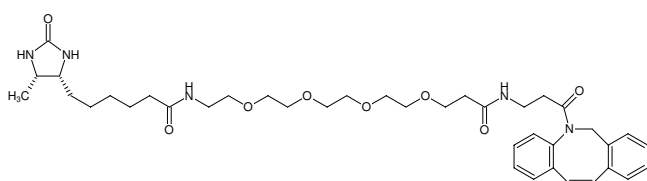




## DBCO-PEG<sub>4</sub>-Desthiobiotin

Dibenzylcyclooctyne-PEG<sub>4</sub>-Desthiobiotin

Cat. No.	Amount
CLK-1108-10	10 mg
CLK-1108-100	100 mg



Structural formula of DBCO-PEG<sub>4</sub>-Desthiobiotin

### 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>39</sub>H<sub>53</sub>N<sub>5</sub>O<sub>8</sub>

**Molecular Weight:** 719.88 g/mol

**Exact Mass:** 719.39 g/mol

**Purity:** ≥ 90 % (HPLC)

**Form:** oil

**Color:** colorless to slightly yellow

**Solubility:** DMSO, DMF, MeOH, DCM, THF

### Description:

DBCO-PEG<sub>4</sub>-Desthiobiotin is suitable for the introduction of a desthiobiotin moiety to Azide-labeled biomolecules via Cu(I)-free strain-promoted Alkyne-Azide Click Chemistry (SPAAC) reaction.

The hydrophilic PEG<sub>4</sub> linker reduces or eliminates aggregation and precipitation during the labeling process by increasing the hydrophilicity of the target molecule. It furthermore enhances the accessibility of the desthiobiotin moiety and thus the detection efficiency of the desthiobiotinylated molecule via fluorescent or HRP-labeled streptavidin or its affinity purification via streptavidin agarose.

Desthiobiotin binds less tightly to streptavidin ( $K_D = 10^{-11}$  M) than Biotin ( $K_D = 10^{-15}$  M). Desthiobiotinylated molecules are therefore easily eluted from the complex in the presence of excess Biotin.