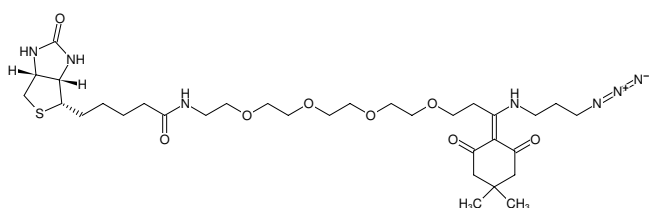




Dde Biotin-Azide

Cat. No.	Amount
CLK-1136-5	5 mg
CLK-1136-25	25 mg



Structural formula of Dde Biotin-Azide

For research use only!**Shipping:** shipped at ambient temperature**Storage Conditions:** store at -20 °C**Shelf Life:** 12 months after date of delivery**Molecular Formula:** C₃₂H₅₃N₇O₈S**Molecular Weight:** 695.87 g/mol**CAS#:** 1802907-93-2**Purity:** ≥ 95 % (HPLC)**Form:** amorphous solid**Solubility:** Chloroform, DCM, DMF, DMSO, THF**Description:**

Dde Biotin-Azide is suitable for the introduction of a biotin moiety to terminal Alkyne- and strained Alkyne (e.g. DBCO)-labeled biomolecules via Cu(I)-catalyzed Alkyne-Azide (CUAAC) or Cu(I)-free strain-promoted Alkyne-Azide Click Chemistry (SPAAC) reaction, respectively.

The biotinylated molecule can be selectively captured through conventional affinity biotin - streptavidin purification and released under mild buffer conditions with hydrazine solution for subsequent identification by mass spectrometry, ELISA, dot blot or Western blot applications.

Important Product Information

If possible, avoid reducing agents in reaction buffers, this can interfere with the azide stability.

General Procedure for Release

Incubate beads in 100 mM phosphate buffer solution (pH 7.4) containing 2%(v/v) hydrazine for 30 - 120 min at room temperature. Reference [1] might be used as starting point for the set up of your individual assay.

Related Products:Copper (II)-Sulphate (CuSO₄), #CLK-MI004

Tris(3-hydroxypropyltriazolylmethyl)amine (THPTA), #CLK-1010

Sodium Ascorbate (Na-Ascorbate), #CLK-MI005

Selected References:

[1] Yang *et al.* (2013) Cleavable Trifunctional Biotin Reagents for Protein Labeling, Capture and Release. *Chem. Commun.* **49**:5366.