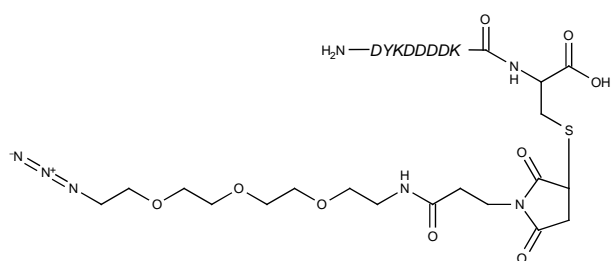


**Azide-PEG₃-FLAG**Azide-PEG₃-FLAGtag (DYKDDDDK)

Cat. No.	Amount
CLK-032-S	0,5 µmol
CLK-032-L	5 x 0,5 µmol

Structural formula of Azide-PEG₃-FLAG**For research use only!****Shipping:** shipped on gel packs**Storage Conditions:** store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery**Molecular Formula:** C₅₉H₈₈N₁₆O₂₇S**Molecular Weight:** 1485.49 g/mol**Exact Mass:** 1484.57 g/mol**Purity:** ≥ 95 % (HPLC)**Form:** solid**Color:** white to off-white**Solubility:** water, PBS (up to 30 mM tested)**Description:**

Azide-PEG₃-FLAG enables the FLAG-tag (DYKDDDDK) attachment to any terminal Alkyne- or Cyclooctyne (e.g. DBCO)-functionalized molecule via Cu(I)-catalyzed terminal Alkyne-Azide Click Chemistry (CuAAC) or Cu(I)-free strain-promoted Alkyne-Azide Click Chemistry (SPAAC), respectively.

The resulting FLAG-tagged molecules can subsequently be detected by an anti-FLAG antibody that is either immobilized onto a matrix (for purification) or coupled with a fluorescent dye or reporter enzyme for direct or indirect detection, respectively. Molecule solubility and efficient FLAG-tag detection is ensured by the integrated PEG-linker.

Azide-FLAG (without PEG-linker) has been successfully used for the detection of Azide-tagged glycoproteins in cell lysates (final concentration: 100 µM)^[1]. This concentrations may serve as a starting point for individual assay set-up.

Related Products:DBCO-PEG₄-FLAG, #CLK-033**Selected References:**

[1] Wang *et al.* (2011) Sulfated Ligands for the Copper (I)-catalyzed Azide-Alkyne Cycloaddition. *Chem Asian J.* **6** (10):2796.