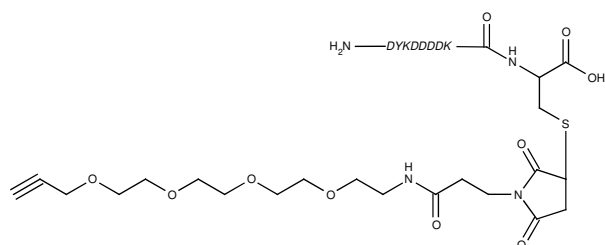




### Alkyne-PEG<sub>4</sub>-FLAG

Alkyne-PEG<sub>4</sub>-FLAGtag (DYKDDDDK)

| Cat. No. | Amount       |
|----------|--------------|
| CLK-088S | 0,5 µmol     |
| CLK-088L | 5 x 0,5 µmol |



Structural formula of Alkyne-PEG<sub>4</sub>-FLAG

#### For general laboratory use.

**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:** C<sub>62</sub>H<sub>91</sub>N<sub>13</sub>O<sub>28</sub>S

**Molecular Weight:** 1498.53 g/mol

**Exact Mass:** 1497.58 g/mol

**Purity:** ≥ 95 % (HPLC)

**Form:** solid

**Color:** white to off-white

**Solubility:** water, PBS (up to 30 mM tested)

#### Description:

Alkyne-PEG<sub>4</sub>-FLAG enables the FLAG-tag (DYKDDDDK) attachment to any Azide-functionalized molecule via Cu(I)-catalyzed terminal Alkyne-Azide Click Chemistry (CuAAC).

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.

#### Related Products:

DBCO-PEG<sub>4</sub>-FLAG, #CLK-033

Azide-PEG<sub>3</sub>-FLAG, #CLK-032