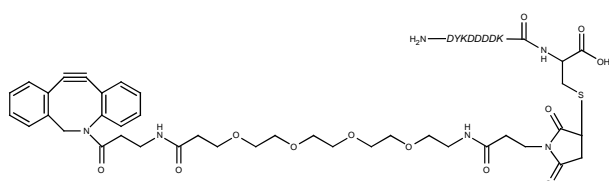


**DBCO-PEG₄-FLAG**Dibenzylcyclooctyne-PEG₄-FLAGtag (DYKDDDDK)

Cat. No.	Amount
CLK-033-S	0,5 µmol
CLK-033-L	5 x 0,5 µmol

Structural formula of DBCO-PEG₄-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:** C₈₀H₁₀₇N₁₅O₃₀S**Molecular Weight:** 1790.85 g/mol**Exact Mass:** 1789.70 g/mol**Purity:** ≥ 95 % (HPLC)**Form:** solid**Color:** white to off-white**Solubility:** water, PBS (up to 30 mM tested)**Spectroscopic Properties:** λ_{max} 308 nm, ε 13.8.0 L mmol⁻¹ cm⁻¹**Description:**

DBCO-PEG₄-FLAG enables the FLAG-tag (DYKDDDDK) attachment to any Azide-functionalized molecule via Cu(I)-free strain-promoted Alkyne-Azide Click Chemistry (SPAAC). 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.

DBCO-FLAG (without PEG-linker) has been successfully used for the detection of cell-surface Azides (final concentration: 100 µM)^[1] or labeling of Azide-functionalized micelles (amount: 100 nmol)^[2]. These concentrations may serve as a starting point for individual assay set-up.

Related Products:Azide-PEG₃-FLAG, #CLK-032**Selected References:**

[1] Patterson *et al.* (2014) Improved cyclopropene reporters for probing protein glycosylation. *Molecular BioSystems* **10**:1693.

[2] Chan *et al.* (2013) Double Click: Dual Functionalized Polymeric Micelles with Antibodies and Peptides. *Bioconjugate Chemistry* **24**:105.