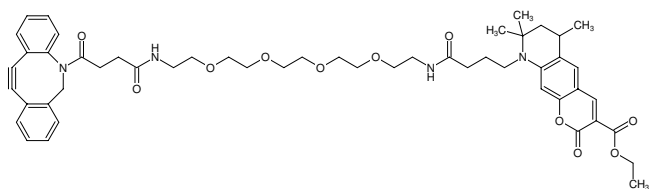
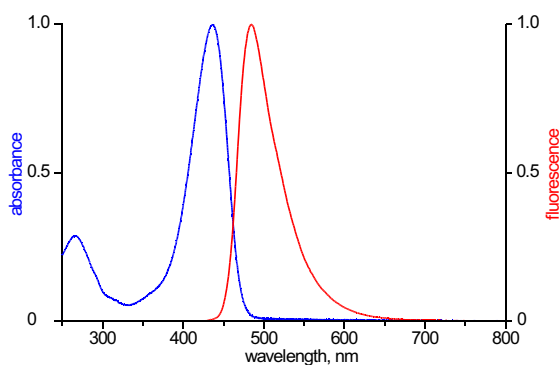


**DBCO-PEG₄-ATTO-425**

Abs/Em = 436/484 nm

Dibenzylcyclooctyne-PEG₄-ATTO-425

Cat. No.	Amount
CLK-085	0,5 mg

Structural formula of DBCO-PEG₄-ATTO-425

excitation and emission spectrum of ATTO 425

For research use only!**Shipping:** shipped at ambient temperature**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:** C₅₁H₆₂N₄O₁₁**Molecular Weight:** 907.06 g/mol**Exact Mass:** 906.44 g/mol**Purity:** ≥ 90 % (HPLC)**Form:** solid**Color:** yellow**Solubility:** DMF, DMSO**Spectroscopic Properties:** λ_{exc} 436 nm, λ_{em} 484 nm, ε 45.0 L mmol⁻¹ cm⁻¹ (after click reaction)**Selected References:**[1] Liu *et al.* (2016) Impact of LytR-CpsA-Psr Proteins on Cell Wall Biosynthesis in *Corynebacterium glutamicum*. *J Bacteriol* **198**:3045.