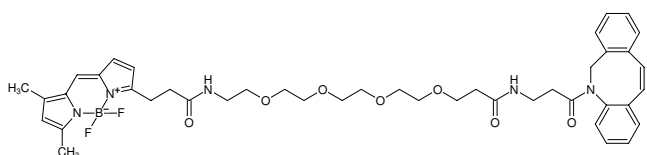


**BDP-FL-PEG₄-DBCO**also known as BODIPY® FL-PEG₄-DBCO

Abs/Em = 503/512 nm

Cat. No.	Amount
CLK-040-05	0,5 mg

Structural formula of BDP-FL-PEG₄-DBCO**For general laboratory use.****Shipping:** shipped on gel packs**Storage Conditions:** store at -20 °C**Additional Storage Conditions:** store dry

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

Shelf Life: 12 months after date of delivery**Molecular Formula:** C₄₃H₅₀BF₂N₅O₇**Molecular Weight:** 797.69 g/mol**Exact Mass:** 797.38 g/mol**Purity:** ≥ 90 % (HPLC)**Form:** solid**Color:** orange**Solubility:** DCM, DMF, DMSO, MeOH**Spectroscopic Properties:** λ_{exc} 503 nm, λ_{em} 512 nm, ε 80.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5 / Methanol 1:1)