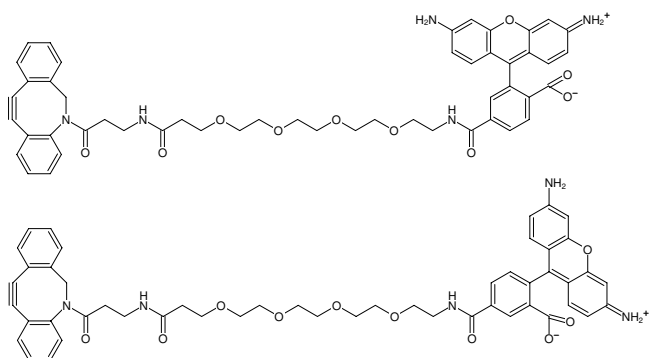
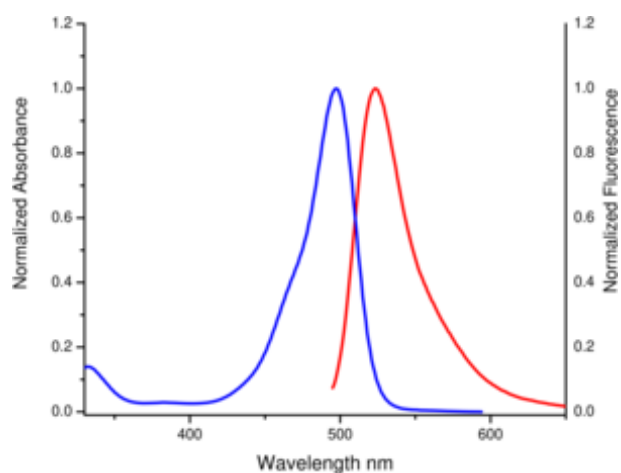



DBCO-PEG₄-5/6-Carboxyrhodamine 110

Abs/Em = 501/526 nm

Dibenzylcyclooctyne-PEG₄-5/6-Carboxyrhodamine 110DBCO-PEG₄-Rhodamine Green

Cat. No.	Amount
CLK-A127-1	1 mg
CLK-A127-5	5 x 1 mg
CLK-A127-25	25 mg

For research use only!**Shipping:** shipped at ambient temperature**Storage Conditions:** store at -20 °C**Shelf Life:** 12 months after date of delivery**Molecular Formula:** C₅₀H₄₉N₅O₁₀**Molecular Weight:** 879.95 g/mol**Exact Mass:** 879.34 g/mol**Purity:** ≥ 90 % (HPLC)**Form:** solid**Color:** red**Solubility:** DMF, DMSO**Spectroscopic Properties:** λ_{abs} 501 nm, λ_{em} 526 nm, ε 74.0 L mmol⁻¹ cm⁻¹Structural formula of DBCO-PEG₄-5/6-Carboxyrhodamine 110

excitation and emission spectrum of 5/6-Carboxyrhodamine 110