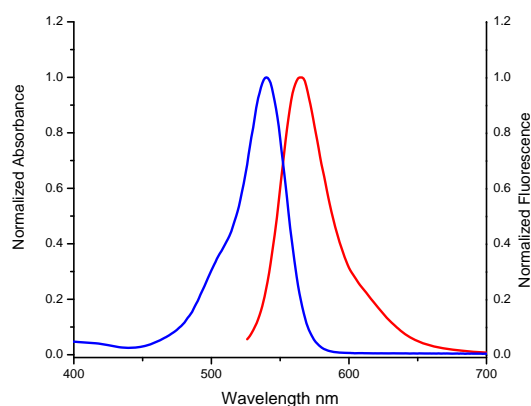
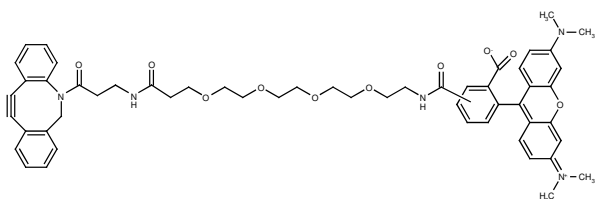


Dibenzylcyclooctyne-PEG4-Fluor 545

Cat. No.	Amount
CLK-A110N-2	2 mg
CLK-A110N-5	5 mg



Absorption and emission spectrum Fluor 545

Molecular formula: C₅₄H₅₉N₅O₁₀

Molecular weight: 938.07 g/mol

Spectroscopic properties:

λ_{abs} 546 nm; λ_{em} 565 nm; ϵ 89,000 cm⁻¹ M⁻¹ (in MeOH)

Storage conditions: store undissolved at -20°C, for use prepare a fresh solution

Purity: >90% (HPLC)

Appearance: dark red solid

Shelf life: 12 months (undissolved)

Solubility: DMSO, DMF, DCM, MeOH

Product Features and Benefits:

With excitation maximum at 546 nm, the dibenzylcyclooctyne - conjugate is an excellent match to the intense 546 nm spectral line of the mercury-arc lamps used in most fluorescence microscopes. This probe is also efficiently excited by the 543 nm spectral line of the Ar - Kr mixed gas laser commonly used in many confocal laser-scanning microscopes.

The fluorescence of the dibenzylcyclooctyne - tetramethylrhodamine conjugate as well as of many other rhodamine based probes is pH insensitive between 4 and 9, and exhibits excellent photostability.

For research use only!