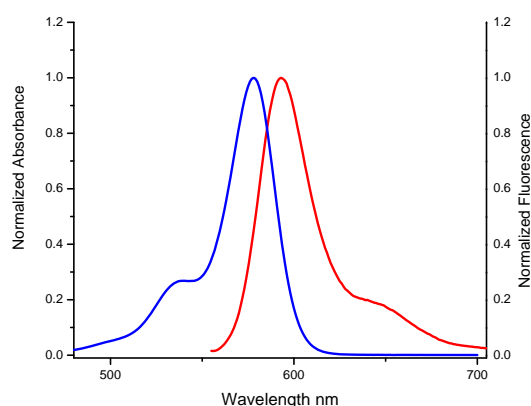
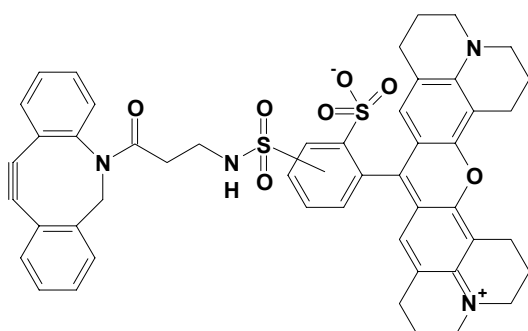


Dibenzylcyclooctyne-Fluor 585

Cat. No.	Amount
CLK-A111-2	2 mg
CLK-A111-5	5 mg



Absorption and emission spectrum Fluor 585

Molecular formula: C₄₉H₄₄N₄O₇S₂

Molecular weight: 865.03 g/mol

Spectroscopic properties:

λ_{abs} 584 nm; λ_{em} 603 nm, ϵ 110000 cm⁻¹ M⁻¹ (in MeOH)

Storage conditions: store at -20°C

Purity: >90% (H NMR)

Appearance: dark red solid

Shelf life: 12 months

Solubility: DMSO, DMF, DCM, Chloroform, MeOH

Product Features and Benefits:

This Sulforhodamine 101 (commonly known as Texas Red) based Dibenzylcyclooctyne-Fluor 585 probe emits at a longer wavelength than do either tetramethylrhodamine or Lissamine™ rhodamine B conjugates. Unlike other rhodamine based probes, the Dibenzylcyclooctyne-Fluor 585 exhibits very little spectral overlap with fluorescein. With peak absorption at 584 nm, Dibenzylcyclooctyne-Fluor 585 probe is particularly well suited for excitation by the 568 nm spectral line of the Ar – Kr mixed gas laser commonly used in many confocal laser-scanning microscopes or the 594 nm spectral line of the orange He-Ne laser. The fluorescence quantum yield of the Dibenzylcyclooctyne-Fluor 585 probe is higher than that of tetramethylrhodamine or Lissamine™ rhodamine B. Usually, conjugates of Dibenzylcyclooctyne-Fluor 585 exhibit higher fluorescence intensity than conjugates of other rhodamine based probes.

For research use only!