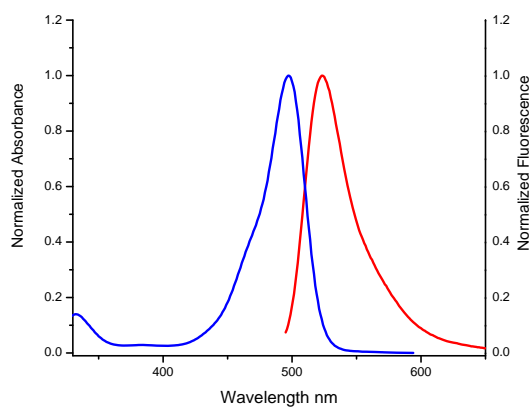
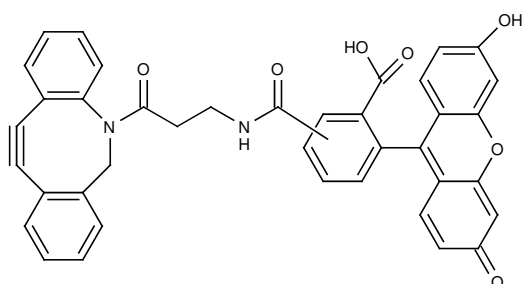


Dibenzylcyclooctyne-Fluor 488

Cat. No.	Amount
CLK-A106N-2	2 mg
CLK-A106N-5	5 mg



Absorption and emission spectrum of Fluor 488

Molecular formula: C₃₉H₂₇N₂O₇

Molecular weight: 635.64 g/mol

Spectroscopic properties:

λ_{abs} 492 nm; λ_{em} 526 nm;

ϵ 79,000 cm⁻¹ M⁻¹ (in phosphate buffer pH 7.4)

Storage conditions: store at -20°C

Purity: >90% (HPLC)

Appearance: yellow solid

Shelf life: 12 months

Solubility: DMSO, DMF, DCM, methanol

Product Features and Benefits: With excitation maximum at 492 nm, the Dibenzylcyclooctyne - Fluorescein conjugate is an excellent match to the intense 488 nm spectral line of the argon ion laser commonly used in many confocal laser-scanning microscopes.

For research use only!

Selected References:

Hao *et al.* (2011) A readily synthesized cyclic pyrrolysine analogue for site-specific protein "click" labeling. *Chem Commun (Camb)* **47(15)**:4502.