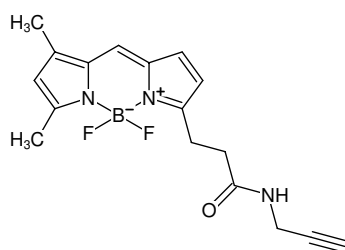


**Alkyne-BDP-FL**

Abs/Em = 503/512 nm

also known as Alkyne-BODIPY® FL

Cat. No.	Amount
CLK-045-1	1 mg
CLK-045-5	5 x 1 mg



Structural formula of Alkyne-BDP-FL

For general laboratory use.**Shipping:** shipped at ambient temperature**Storage Conditions:** store at -20 °C**Additional Storage Conditions:** store dark and dry**Shelf Life:** 12 months after date of delivery**Molecular Formula:** C₁₇H₁₈BF₂N₃O**Molecular Weight:** 329.16 g/mol**Exact Mass:** 329.15 g/mol**Purity:** ≥ 90 % (HPLC)**Form:** solid**Color:** orange**Solubility:** High: DMF, DMSO, DichlormethanLow: H₂O**Spectroscopic Properties:** λ_{exc} 503 nm, λ_{em} 512 nm,
ε 80.0 L mmol⁻¹ cm⁻¹