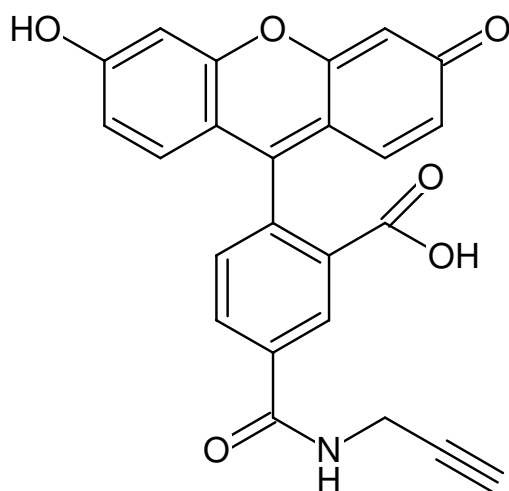


**5-FAM-Alkyne**

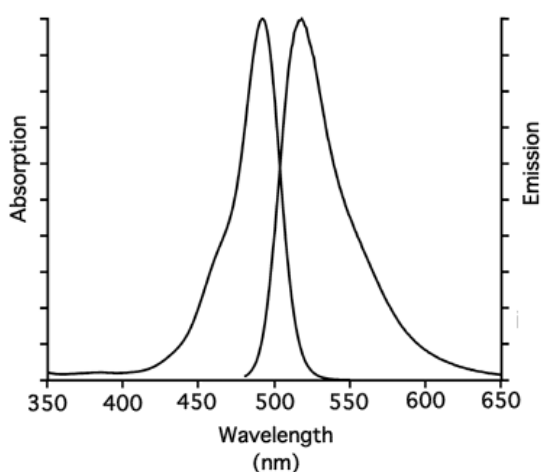
Abs/Em = 490/513 nm

5-Fluorescein-Alkyne

Cat. No.	Amount
CLK-057-5	5 mg
CLK-057-25	5 x 5 mg



Structural formula of 5-FAM-Alkyne



excitation and emission spectrum of 5-FAM

For general laboratory use.**Shipping:** shipped at ambient temperature**Storage Conditions:** store at -20 °C**Additional Storage Conditions:** store dark**Shelf Life:** 12 months after date of delivery**Molecular Formula:** C₂₄H₁₅NO₆**Molecular Weight:** 413.38 g/mol**Exact Mass:** 413.09 g/mol**CAS#:** 510758-19-7**Purity:** ≥ 90 % (HPLC)**Form:** solid**Color:** yellow**Solubility:** aqueous buffer (pH >8), alcohols, DMSO, DMF**Spectroscopic Properties:** λ_{abs} 490 nm, λ_{em} 513 nm, ε 80.0 L mmol⁻¹ cm⁻¹