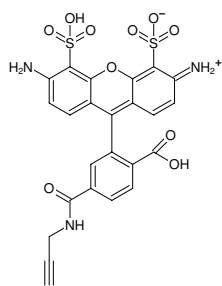
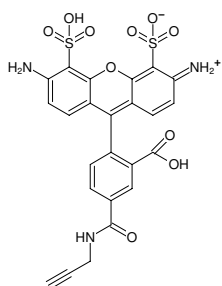


**AF488-Alkyne**

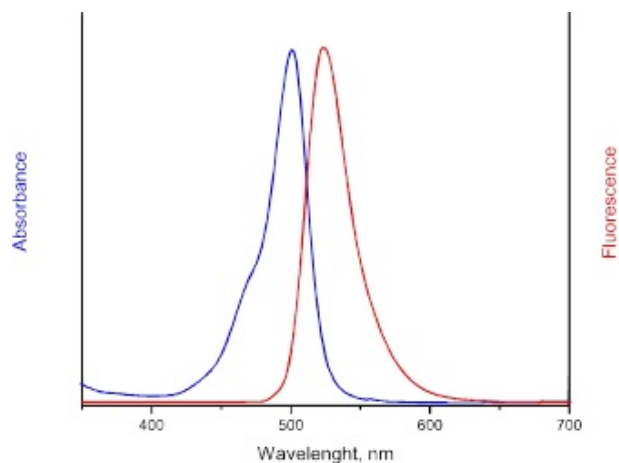
Abs/Em = 494/517 nm

also known as Alexa Fluor® 488-Alkyne

Cat. No.	Amount
CLK-1277-1	1 mg
CLK-1277-5	5 x 1 mg



Structural formula of AF488-Alkyne

For research use only!**Shipping:** shipped at ambient temperature**Storage Conditions:** store at -20 °C**Shelf Life:** 12 months after date of delivery**Molecular Formula:** C₂₄H₁₇N₃O₁₀S₂**Molecular Weight:** 571.54 g/mol**Exact Mass:** 571.04 g/mol**Purity:** > 95 % (HPLC)**Form:** solid**Color:** orange to red**Solubility:** DMSO, DMF, MeOH**Spectroscopic Properties:** λ_{exc} 494 nm, λ_{em} 517 nm, ε 73.0 L mmol⁻¹ cm⁻¹

excitation and emission spectrum of AF488