



AF488-Picolyl-Azide

Abs/Em = 494/517 nm

also known as Alexa Fluor® 488-Picolyl-Azide

Cat. No.	Amount
CLK-1276-1	1 mg
CLK-1276-5	5 x 1 mg

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: C₃₀H₂₄N₈O₁₁S₂

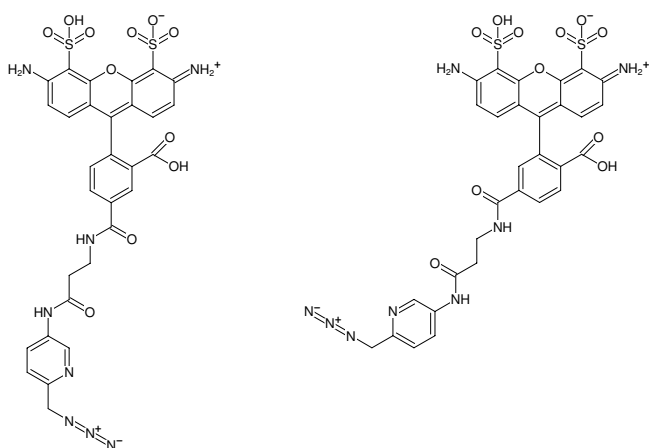
Molecular Weight: 736.69 g/mol

Purity: > 95 % (HPLC)

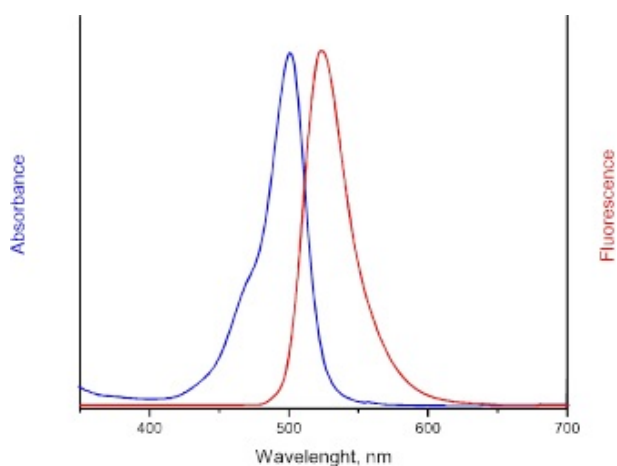
Form: orange to red solid

Solubility: DMSO, DMF, MeOH

Spectroscopic Properties: λ_{exc} 494 nm, λ_{em} 517 nm, ε 73.0 L mmol⁻¹ cm⁻¹



Structural formula of AF488-Picolyl-Azide



excitation and emission spectrum of AF488