

**AF555-Alkyne**

Abs/Em = 555/572 nm
structural analog to Alexa Fluor® 555

Cat. No.	Amount
CLK-1289-1	1 mg
CLK-1289-5	5 x 1 mg

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at -20 °C

Shelf Life: 12 months after date of delivery

Molecular Formula: C₃₇H₄₇N₃O₁₃S₄

Molecular Weight: 870.03 g/mol

Exact Mass: 869.20 g/mol

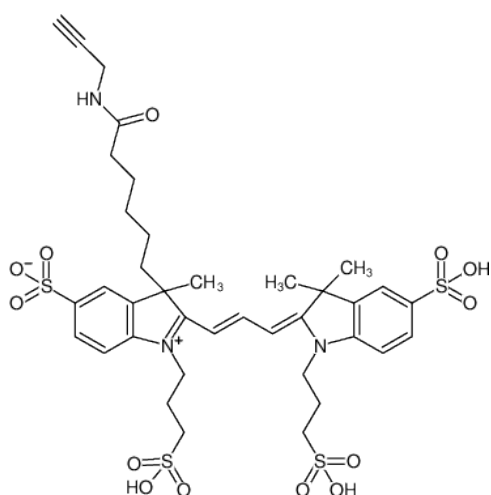
Purity: ≥ 90 % (HPLC)

Form: solid

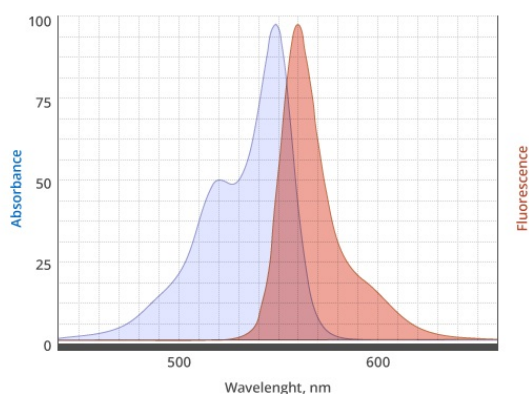
Color: red

Solubility: DMSO, DMF

Spectroscopic Properties: λ_{exc} 555 nm, λ_{em} 572 nm, ε 155.0 L mmol⁻¹ cm⁻¹



Structural formula of AF555-Alkyne



excitation and emission spectrum of AF555