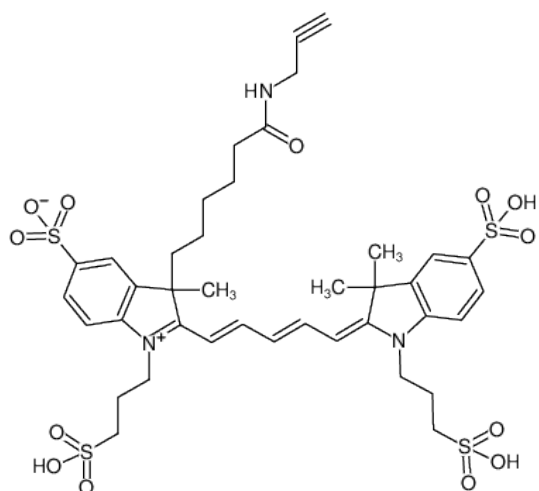




AZDye647A-Alkyne

Abs/Em = 648/671 nm
structural analog to Alexa Fluor® 647

Cat. No.	Amount
CLK-1301A-AZ-1	1 mg
CLK-1301A-AZ-5	5 x 1 mg



Structural formula of AZDye647A-Alkyne

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: C₃₉H₄₉N₃O₁₃S₄

Molecular Weight: 896.07 g/mol

Exact Mass: 895.21 g/mol

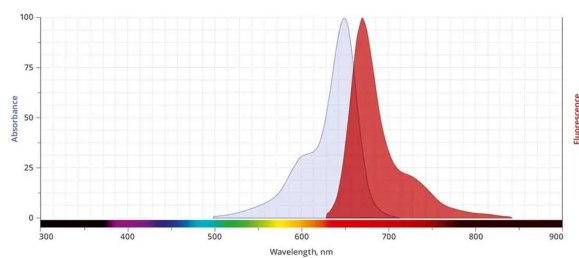
Purity: ≥ 90 % (HPLC)

Form: solid

Color: blue

Solubility: DMSO, DMF

Spectroscopic Properties: λ_{exc} 648 nm, λ_{em} 671 nm, ε 270.0 L mmol⁻¹ cm⁻¹



excitation and emission spectrum of AZDye 647A