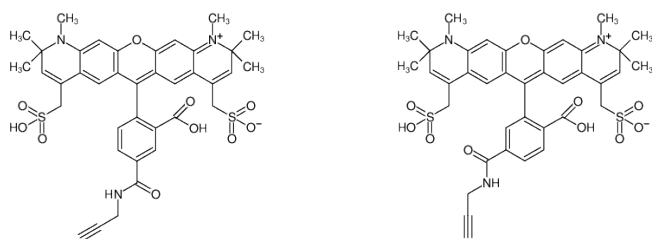




AZDye594-Alkyne

Abs/Em = 590/617 nm
structural analog to Alexa Fluor® 594

Cat. No.	Amount
CLK-1297-AZ-1	1 mg
CLK-1297-AZ-5	5 x 1 mg



Structural formula of AZDye594-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: C₃₈H₃₇N₃O₁₀S₂

Molecular Weight: 759.84 g/mol

Exact Mass: 759.19 g/mol

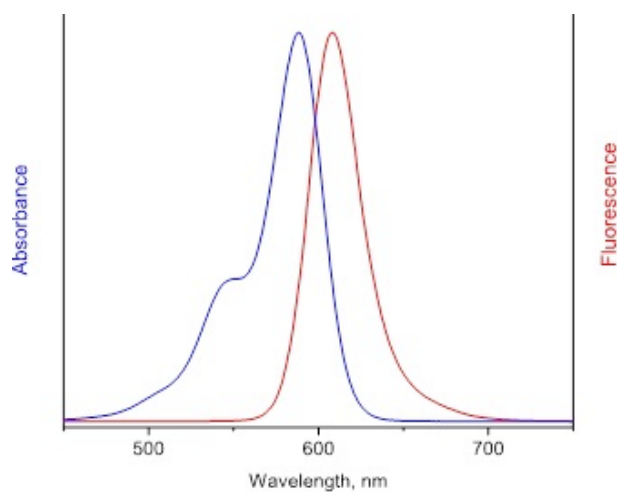
Purity: ≥ 90 % (HPLC)

Form: solid

Color: blue

Solubility: DMSO, DMF, MeOH

Spectroscopic Properties: λ_{exc} 590 nm, λ_{em} 617 nm, ε 92.0 L mmol⁻¹ cm⁻¹



excitation and emission spectrum of AZDye 594