

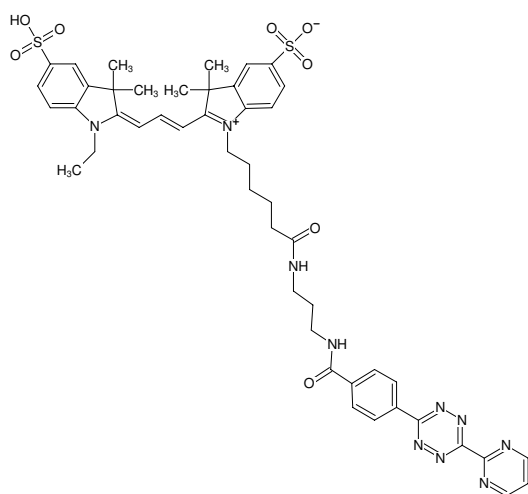


Pyrimidyl-Tetrazine-Cy3

Abs/Em = 550/570 nm

N-(3-Aminopropyl)-4-(6-(pyrimidin-2-yl)-1,2,4,5-tetrazin-3-yl)benzamid - Cy3

Cat. No.	Amount
CLK-100	0,2 mg



Structural formula of Pyrimidyl-Tetrazine-Cy3

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: C₄₇H₅₂N₁₀O₈S₂ (free acid)

Molecular Weight: 949.12 g/mol (free acid)

Exact Mass: 948.34 g/mol (free acid)

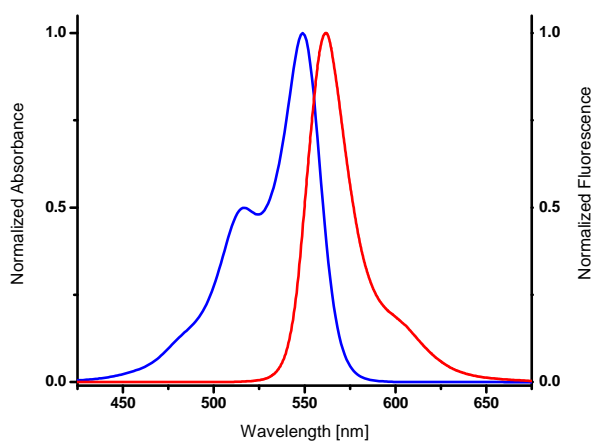
Purity: ≥ 90 % (HPLC)

Form: solid

Color: pink

Solubility: DMF, DMSO, MeOH, water slightly soluble in DCM

Spectroscopic Properties: λ_{exc} 550 nm, λ_{em} 570 nm, ε 150.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5) (after click reaction)



excitation and emission spectrum of Cy3