

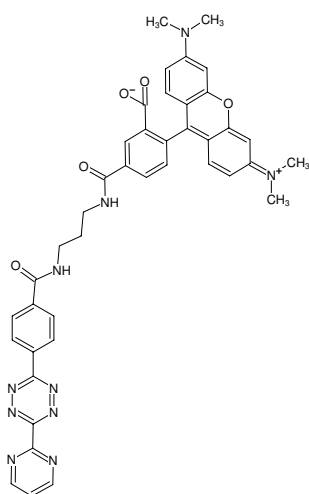


Pyrimidyl-Tetrazine-5-TAMRA

Abs/Em = 545/575 nm

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

Cat. No.	Amount
CLK-097	0,5 mg



Structural formula of Pyrimidyl-Tetrazine-5-TAMRA

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: C₄₁H₃₆N₁₀O₅

Molecular Weight: 748.80 g/mol

Exact Mass: 748.29 g/mol

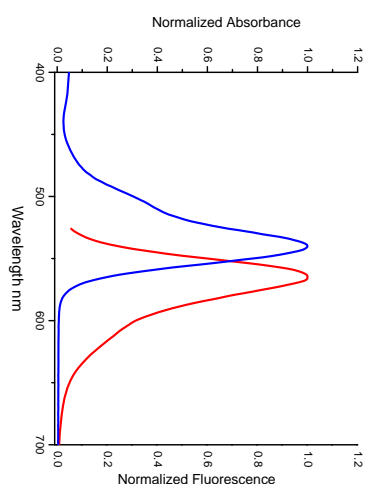
Purity: ≥ 90 % (HPLC)

Form: solid

Color: pink to red

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

Spectroscopic Properties: λ_{exc} 545 nm, λ_{em} 575 nm,
ε 90.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5) (after click reaction)



excitation and emission spectrum of 5-TAMRA