

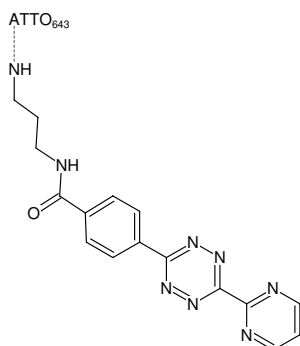


Pyrimidyl-Tetrazine-ATTO-643

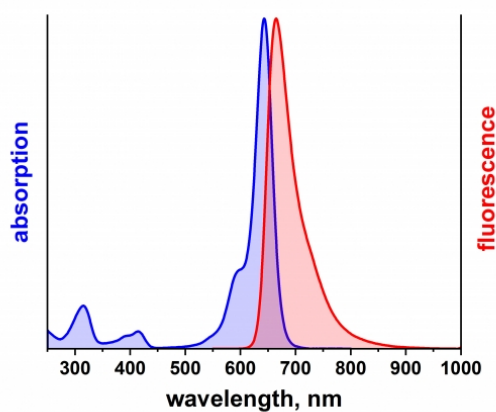
Abs/Em = 643/665 nm

N-(3-Aminopropyl)-4-(6-(pyrimidin-2-yl)-1,2,4,5-tetrazin-3-yl)benzamid - ATTO-643

Cat. No.	Amount
CLK-101	0,2 mg



Structural formula of Pyrimidyl-Tetrazine-ATTO-643



excitation and emission spectrum of ATTO 643

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: C₁₆H₁₅N₈O - ATTO 643

Molecular Weight: 1154 g/mol

Purity: ≥ 90 % (HPLC)

Form: solid

Color: blue

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

Spectroscopic Properties: λ_{exc} 643 nm, λ_{em} 665 nm, ε 150.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5) (after click reaction)