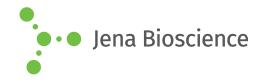
DATA SHEET

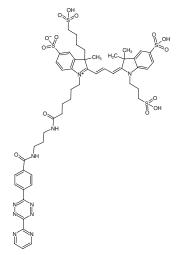




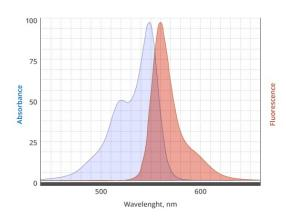
Pyrimidyl-Tetrazine-AF555

Abs/Em = 555/572 nm N-(3-Aminopropyl)-4-(6-(pyrimidin-2-yl)-1,2,4,5-tetrazin-3-yl)benzamid - AF555

Cat. No.	Amount
CLK-098	0,2 mg



Structural formula of Pyrimidyl-Tetrazine-AF555



excitation and emission spectrum of AF555

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

 $\begin{aligned} & \textbf{Molecular Formula:} \ \ C_{51}H_{60}N_{10}O_{14}S_4 \\ & \textbf{Molecular Weight:} \ \ 1165.34 \ g/mol \end{aligned}$

Exact Mass: 1164.32 g/mol **Purity:** ≥ 90 % (HPLC)

Form: solid
Color: red

Solubility: DMF, DMSO, MeOH, water

slightly soluble in DCM

Spectroscopic Properties: λ_{exc} 555 nm, λ_{em} 572 nm, ϵ 155.0 L mmol⁻¹

cm⁻¹

Please note:

Product can not be stored in solution.

Dissolve the product in methanol, split into aliquots of sufficient amounts for one experiment and dry the product e.g. with a speedvac. Dried aliquots can be stored at -20°C and freshly dissolved immediately before use e.g. with DMSO.