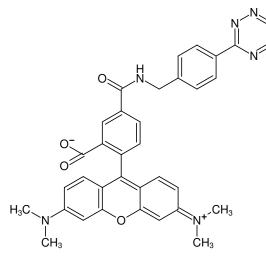




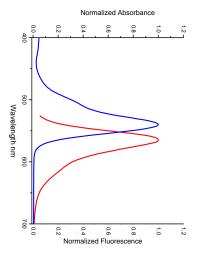
Tetrazine-5-TAMRA

Abs/Em = 545/575 nm 3-(p-Benzylamino)-1,2,4,5-tetrazine - 5-Carboxytetramethylrhodamine

Cat. No.	Amount
CLK-017-05	0,5 mg



Structural formula of Tetrazine-5-TAMRA



excitation and emission spectrum of 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: C₃₄H₂₉N₇O₄

Molecular Weight: 599.64 g/mol

Exact Mass: 599.23 g/mol

Purity: ≥ 90 % (HPLC)

Form: solid

Color: pink to red

Solubility: DMF, DMSO slightly soluble in DCM, MeOH insoluble in Water

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)

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.

Selected References:

Liu *et al.* (2012) Diels-Alder cycloaddition for fluorophore targeting to specific proteins inside living cells. *J. Am. Chem. Soc.* **134 (2)**:792.

