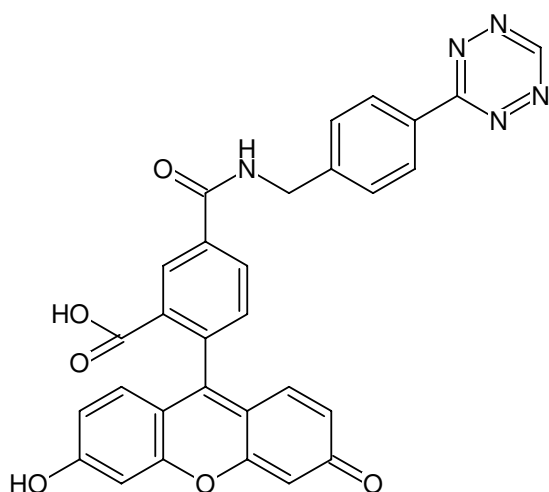


**Tetrazine-5-FAM**

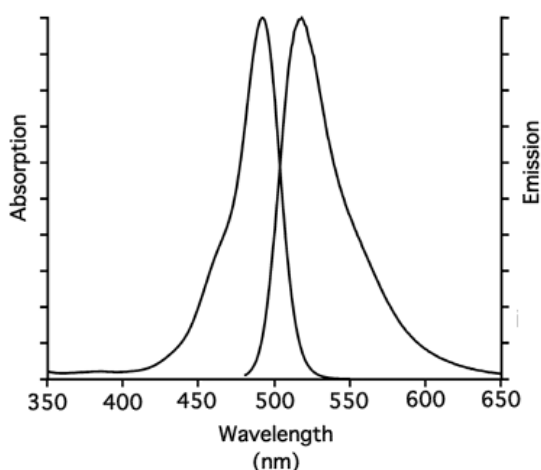
Abs/Em = 492/517 nm

3-(p-Benzylamino)-1,2,4,5-tetrazine – 5-Fluorescein

Cat. No.	Amount
CLK-013-05	0,5 mg



Structural formula of Tetrazine-5-FAM



excitation and emission spectrum of 5-FAM

For research use only!**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:** C₃₀H₁₉N₅O₆ (free acid)**Molecular Weight:** 545.50 g/mol (free acid)**Exact Mass:** 545.13 g/mol (free acid)**Purity:** ≥ 90 % (HPLC)**Form:** solid**Color:** yellow to orange**Solubility:** DCM, DMF, DMSO, MeOH**Spectroscopic Properties:** λ_{exc} 492 nm, λ_{em} 517 nm, ε 83.0 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 speed-vac. Dried aliquots can be stored at -20°C and freshly dissolved immediately before use e.g. with DMSO.