

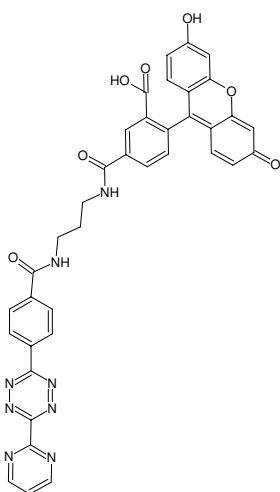


## Pyrimidyl-Tetrazine-5-FAM

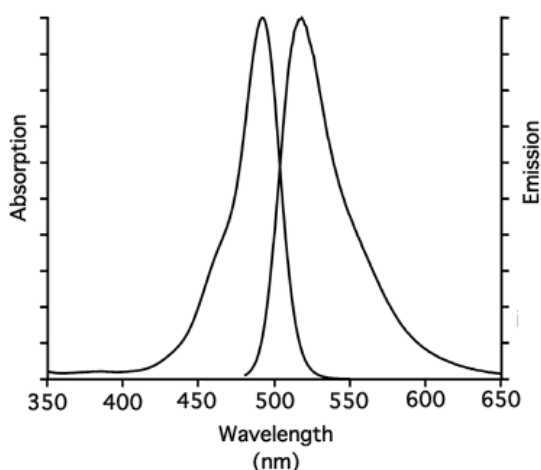
Abs/Em = 492/517 nm

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

Cat. No.	Amount
CLK-096	0,5 mg



Structural formula of Pyrimidyl-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:** C<sub>37</sub>H<sub>26</sub>N<sub>8</sub>O<sub>7</sub> (free acid)

**Molecular Weight:** 694.66 g/mol (free acid)

**Exact Mass:** 694.19 g/mol (free acid)

**Purity:** ≥ 90 % (HPLC)

**Form:** solid

**Color:** yellow to orange

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

**Spectroscopic Properties:**  $\lambda_{exc}$  492 nm,  $\lambda_{em}$  517 nm,  
 $\epsilon$  83.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5) (after click reaction)