

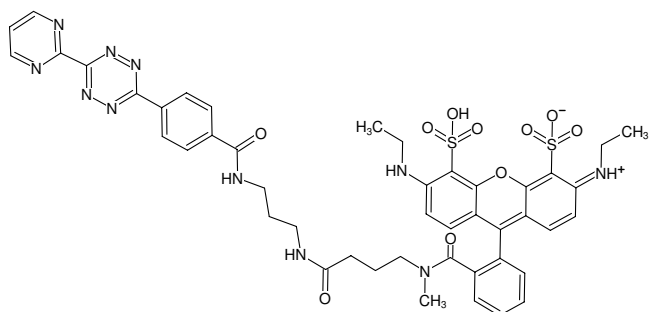


## Pyrimidyl-Tetrazine-ATTO-532

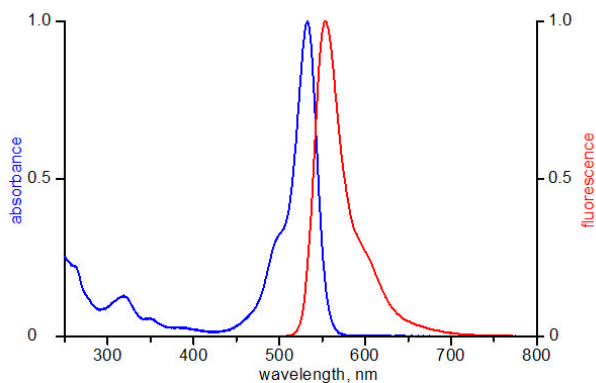
Abs/Em = 532/553 nm

N-(3-Aminopropyl)-4-(6-(pyrimidin-2-yl)-1,2,4,5-tetrazin-3-yl)benzamid - ATTO 532

Cat. No.	Amount
CLK-105	0,2 mg



Structural formula of Pyrimidyl-Tetrazine-ATTO-532



excitation and emission spectrum of ATTO 532

### 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>45</sub>H<sub>45</sub>N<sub>11</sub>O<sub>10</sub>S<sub>2</sub> (free acid)

**Molecular Weight:** 964.04 g/mol (free acid)

**Exact Mass:** 963.28 g/mol (free acid)

**Purity:** ≥ 90 % (HPLC)

**Form:** solid

**Color:** pink

**Solubility:** DCM, DMF, MeOH

**Spectroscopic Properties:** λ<sub>exc</sub> 532 nm, λ<sub>em</sub> 553 nm, ε 115.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (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.