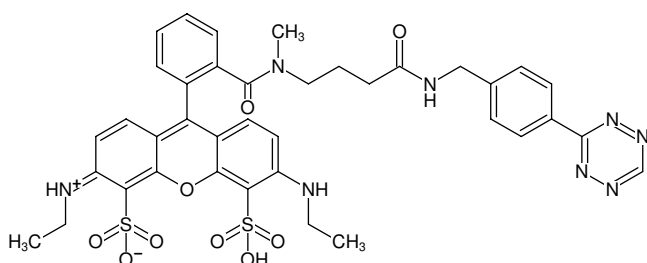


**Tetrazine-ATTO-532**

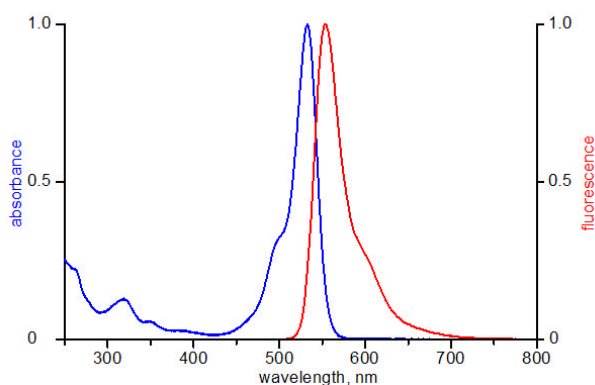
Abs/Em = 532/553 nm

3-(*p*-Benzylamino)-1,2,4,5-tetrazine - ATTO 532

Cat. No.	Amount
CLK-011-02	0,2 mg



Structural formula of Tetrazine-ATTO-532



excitation and emission spectrum of ATTO 532

**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:** C<sub>38</sub>H<sub>38</sub>N<sub>8</sub>O<sub>9</sub>S<sub>2</sub>**Molecular Weight:** 814.89 g/mol**Exact Mass:** 814.22 g/mol**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 speedvac. Dried aliquots can be stored at -20°C and freshly dissolved immediately before use e.g. with DMSO.