

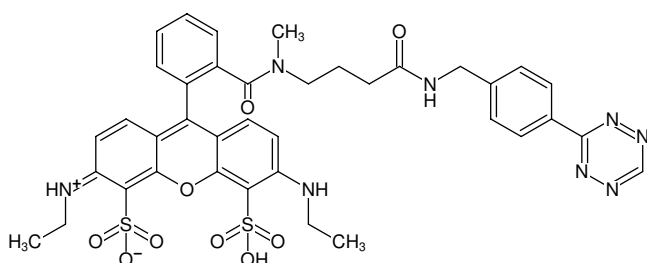


Tetrazine-ATTO-532

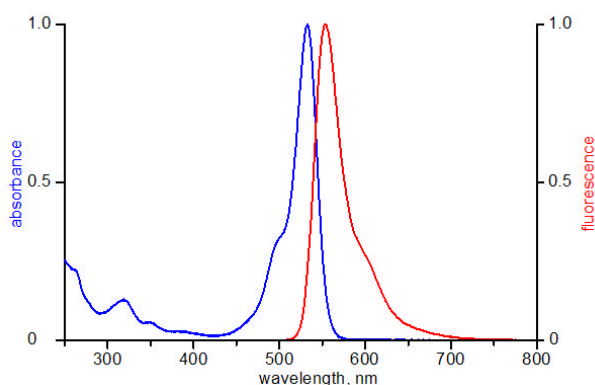
Abs/Em = 532/553 nm

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 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: C₃₈H₃₈N₈O₉S₂

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_{exc} 532 nm, λ_{em} 553 nm, ε 115.0 L mmol⁻¹ cm⁻¹ (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.