



## 6-Methyl-Tetrazine-ATTO-488

Abs/Em = 501/523 nm

(4-(6-Methyl-1,2,4,5-tetrazin-3-yl)phenyl)methanamine - ATTO 488

Cat. No.	Amount
CLK-021-02	0,2 mg

**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>35</sub>H<sub>32</sub>N<sub>8</sub>O<sub>9</sub>S<sub>2</sub>

**Molecular Weight:** 772.81 g/mol

**Exact Mass:** 772.17 g/mol

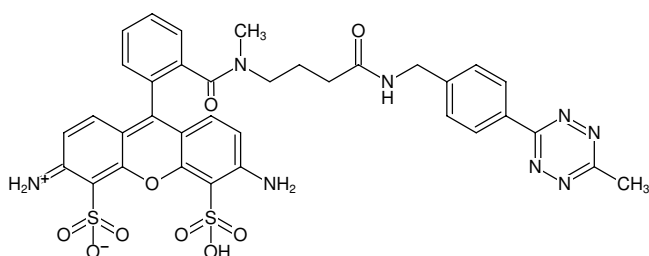
**Purity:** ≥ 90 % (HPLC)

**Form:** solid

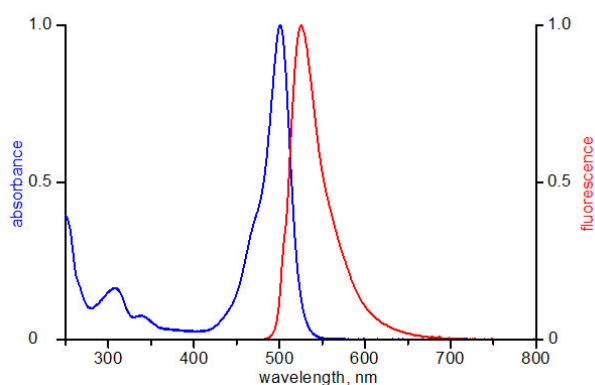
**Color:** orange

**Solubility:** DCM, DMF, MeOH

**Spectroscopic Properties:** λ<sub>exc</sub> 501 nm, λ<sub>em</sub> 523 nm, ε 90.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (after click reaction)



Structural formula of 6-Methyl-Tetrazine-ATTO-488



excitation and emission spectrum of ATTO 488