



## $\gamma$ -[[6-Aminoethyl)-imido]-ATP $\alpha$ S-ATTO-488

$\gamma$ -[[6-Aminoethyl)-imido]-adenosine-5'-( $\alpha$ -thio)-triphosphate, labeled with ATTO-488, Triethylammonium salt

Cat. No.	Amount
NU-278-488	60 $\mu$ l (1 mM)

**For research use only!**

**Shipping:** shipped on blue ice

**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>41</sub>H<sub>51</sub>N<sub>10</sub>O<sub>20</sub>P<sub>3</sub>S<sub>3</sub> (free acid)

**Molecular Weight:** 1193.01 g/mol (free acid)

**Exact Mass:** 1192.17 g/mol (free acid)

**Purity:**  $\geq$  95 % (HPLC)

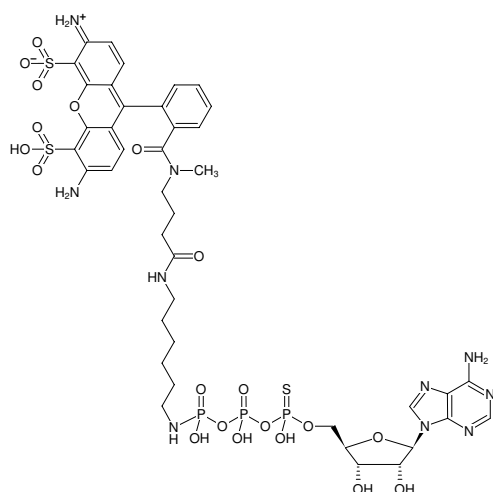
**Form:** solution in 10 mM Tris-HCl

**Color:** yellow

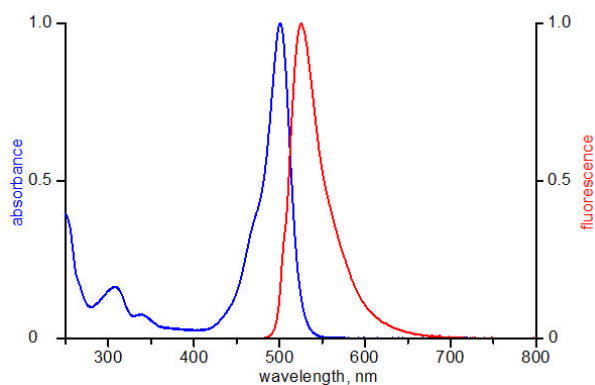
**Concentration:** 1.0 mM - 1.1 mM

**pH:** 8.0  $\pm$  0.5

**Spectroscopic Properties:**  $\lambda_{exc}$  500 nm,  $\lambda_{em}$  520 nm,  $\epsilon$  90.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of  $\gamma$ -[[6-Aminoethyl)-imido]-ATP $\alpha$ S-ATTO-488



excitation and emission spectrum of ATTO 488

**Please note:** This compound contains a phosphoramidate linkage which is hydrolyzed at pH < 7.0.