



### EDA-ATPyS - ATTO-550

2'/3'-O-(2-Aminoethyl-carbamoyl)-adenosine-5'-(γ-thio)-triphosphate, labeled with ATTO 550, Triethylammonium salt

Cat. No.	Amount
NU-1609-550	50 µl (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:** 6 months after date of delivery

**Molecular Formula:** C<sub>13</sub>H<sub>22</sub>N<sub>7</sub>O<sub>13</sub>P<sub>3</sub>S - ATTO 550 (free acid)

**Molecular Weight:** 1184.33 g/mol (free acid)

**Purity:** ≥ 90 % (HPLC), contains approx. 5 % EDA-ADP-dye

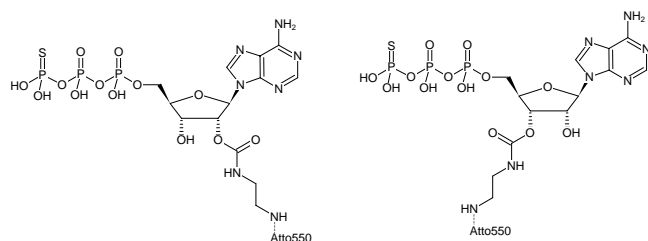
**Form:** red-violet solution in water

**Concentration:** 1.0 mM - 11 mM

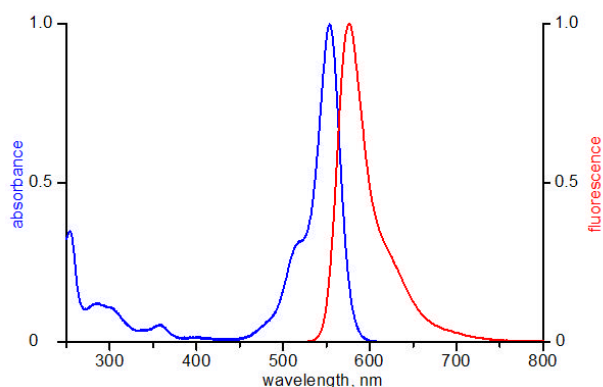
**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 554 nm; λ<sub>em</sub> 576 nm; ε 120.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

The limited stability of labeled ATPyS and GTPyS nucleotides makes it impossible to keep stocks of them and to offer them as catalog products that are available immediately. Therefore, any labeled ATPyS or GTPyS is freshly prepared per order. For this reason the price is higher than for stable nucleotide-dye conjugates which can be synthesized and kept in stock.



Structural formula of EDA-ATPyS - ATTO-550



excitation and emission spectrum of ATTO 550