



## 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 general laboratory use.**

**Shipping:** shipped on dry ice

**Storage Conditions:** store at -20 °C

**Shelf Life:** 6 months after date of delivery

**Molecular Formula:** C<sub>50</sub>H<sub>63</sub>N<sub>10</sub>O<sub>16</sub>P<sub>3</sub>S (free acid)

**Molecular Weight:** 1185.09 g/mol (free acid)

**Exact Mass:** 1184.34 g/mol (free acid)

**Purity:** ≥ 90 % (HPLC)

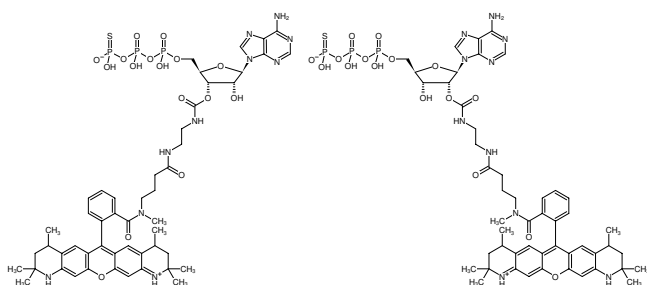
**Form:** solution in water

**Color:** red-violet

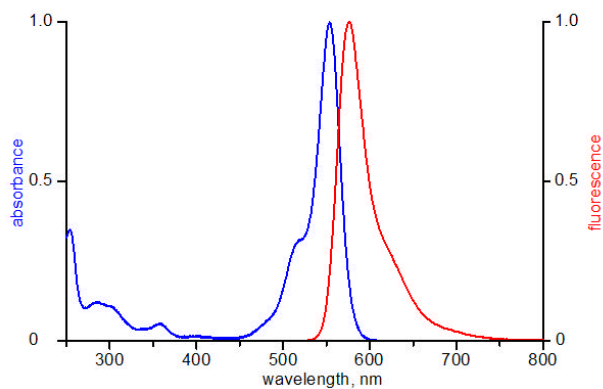
**Concentration:** 1.0 mM - 1.1 mM

**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)



Structural formula of EDA-ATPyS-ATTO-550



excitation and emission spectrum of ATTO 550