



**EDA-AppNHp (EDA-AMPPNP) - ATTO-Thio12**

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-[(β,γ)-imido] triphosphate, labeled with ATTO Thio12, Triethylammonium salt

Cat. No.	Amount
NU-810-THIO12	40 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>42</sub>H<sub>52</sub>N<sub>11</sub>O<sub>15</sub>P<sub>3</sub>S (free acid)

**Molecular Weight:** 1075.91 g/mol (free acid)

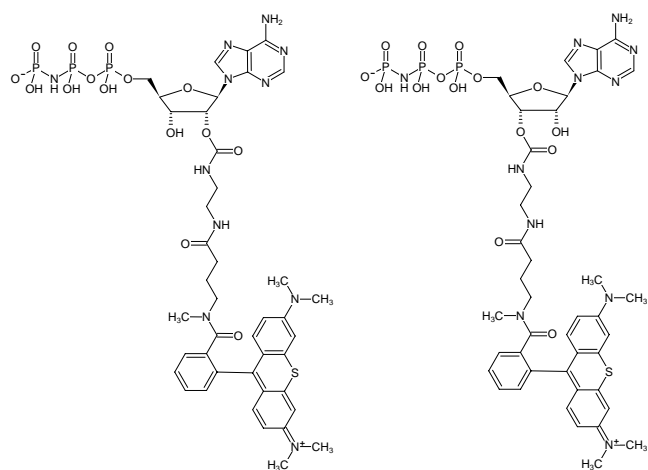
**Purity:** ≥ 95 % (HPLC)

**Form:** clear aqueous solution

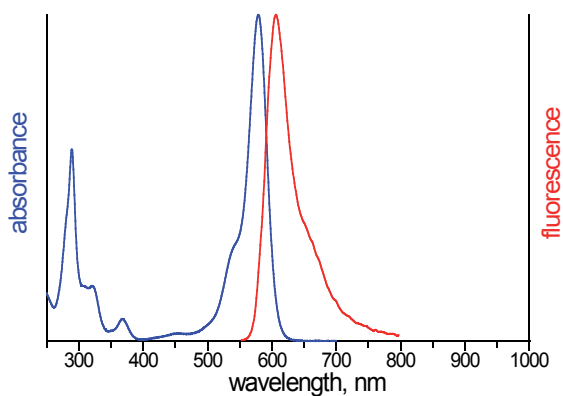
**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 582 nm, λ<sub>em</sub> 607 nm, ε 110.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of EDA-AppNHp (EDA-AMPPNP) - ATTO-Thio12



excitation and emission spectrum of ATTO Thio12