



**EDA-AppNHp (EDA-AMPPNP)-ATTO-Rho13**

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-[(β,γ)-imido] triphosphate, labeled with ATTO Rho13, Triethylammonium salt

Cat. No.	Amount
NU-810-RHO13	40 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>54</sub>H<sub>68</sub>N<sub>11</sub>O<sub>16</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 1220.12 g/mol (free acid)

**Exact Mass:** 1219.41 g/mol (free acid)

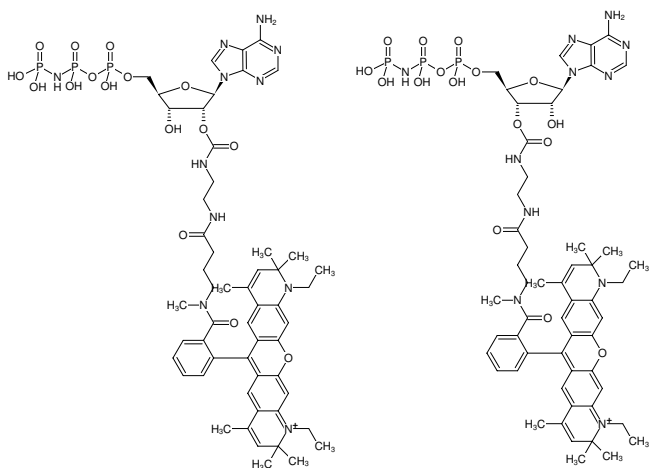
**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

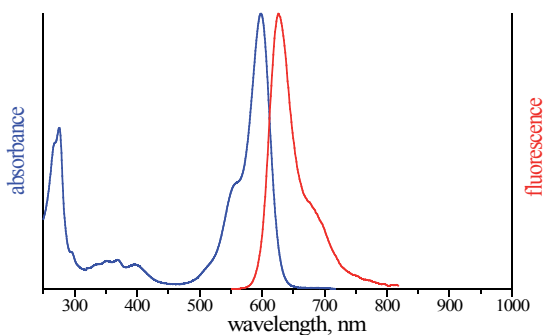
**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 603 nm, λ<sub>em</sub> 627 nm, ε 120.0 L mol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of EDA-AppNHp (EDA-AMPPNP)-ATTO-Rho13



excitation and emission spectrum of ATTO Rho13