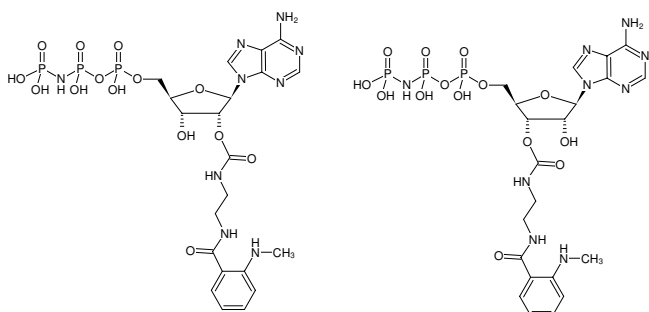




## EDA-AppNHp (EDA-AMPPNP)-MANT

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

Cat. No.	Amount
NU-810-MNT	50 µl (5 mM)



Structural formula of EDA-AppNHp (EDA-AMPPNP)-MANT

### 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>21</sub>H<sub>30</sub>N<sub>9</sub>O<sub>14</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 725.44 g/mol (free acid)

**Exact Mass:** 725.11 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

**Color:** colorless to slightly yellow

**Concentration:** 5.0 mM - 5.5 mM

**pH:** 7.5 ±0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 335 nm, λ<sub>em</sub> 440 nm, ε 2.7 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)