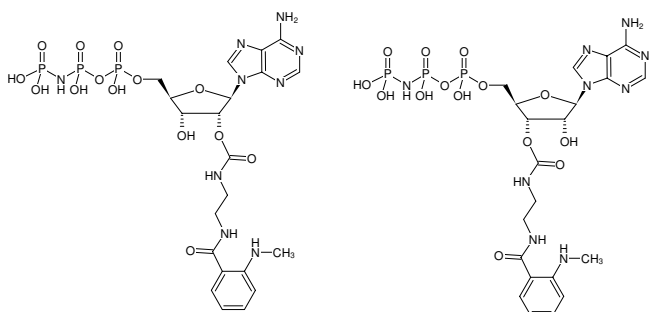



**EDA-AppNHp (EDA-AMPPNP)-MANT**

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 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>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)