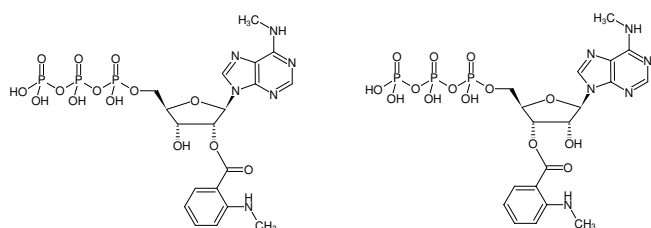




## Mant-N<sup>6</sup>-Methyl-ATP

2'/3'-O-(N-Methyl-anthraniloyl)-N<sup>6</sup>-methyl-adenosine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-216S	30 µl (10 mM)
NU-216L	5 x 30 µl (10 mM)



Structural formula of Mant-N<sup>6</sup>-Methyl-ATP

### 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:** 12 months after date of delivery

**Molecular Formula:** C<sub>19</sub>H<sub>25</sub>N<sub>6</sub>O<sub>14</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 654.36 g/mol (free acid)

**Exact Mass:** 654.06 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** colorless to slightly yellow solution in water

**Concentration:** 10 mM - 11 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>max</sub> 255/355 nm, ε 23.3/5.8 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5), λ<sub>exc</sub> 355 nm, λ<sub>em</sub> 448 nm