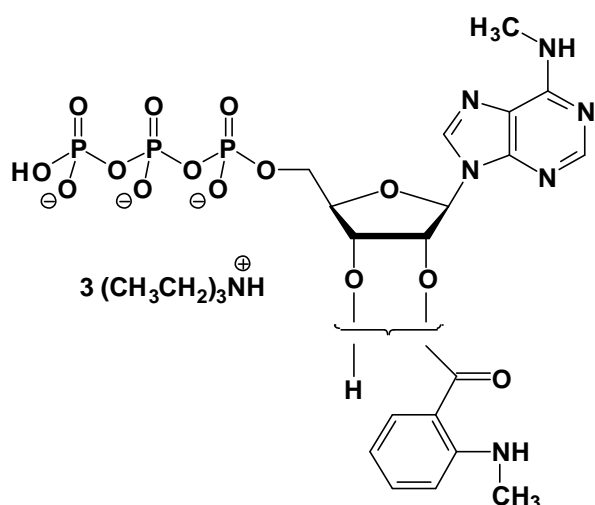


## mant-N<sup>6</sup>-Methyl-ATP

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

Cat. No.	Amount
NU-216S	30 Units
NU-216L	150 Units



**Cat. No.:** NU-216

**Molecular Formula:** C<sub>19</sub>H<sub>21</sub>N<sub>6</sub>O<sub>14</sub>P<sub>3</sub> (Anion)

**Molecular Weight:** 650.32 (Anion)

**Purity:** > 95%, clear aqueous solution, pH 7.5

**Spectroscopic properties:**

$\lambda_{\max}$  255/355 nm;  $\epsilon$  23300/5800;  $\lambda_{\text{exc}}$  355nm;

$\lambda_{\text{em}}$  448 nm

**Storage conditions:**

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

**For research use only!**

1 unit = 1  $\mu$ l of a 10 mM solution