

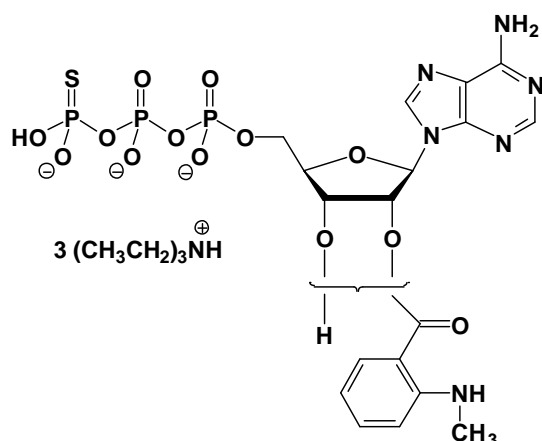
## Mant-ATP $\gamma$ S

2'/3'-O-(N-Methyl-anthraniloyl)-adenosine-5'-( $\gamma$ -thio)-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-232S	20 Units
NU-232L	100 Units

### Selected References:

Gottle *et al.* (2007) Molecular analysis of the interaction of Bordetella pertussis adenylyl cyclase with fluorescent nucleotides. *Molecular Pharmacology Fast Forward*, June 6, 2007; DOI: 10.1124/mol.107.034413 (published online).



**Cat. No.:** NU-232

**Molecular Formula:** C<sub>18</sub>H<sub>20</sub>N<sub>6</sub>O<sub>13</sub>P<sub>3</sub>S (Anion)

**Molecular Weight:** 653.03 (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 6 months after date of delivery.

### For research use only!

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