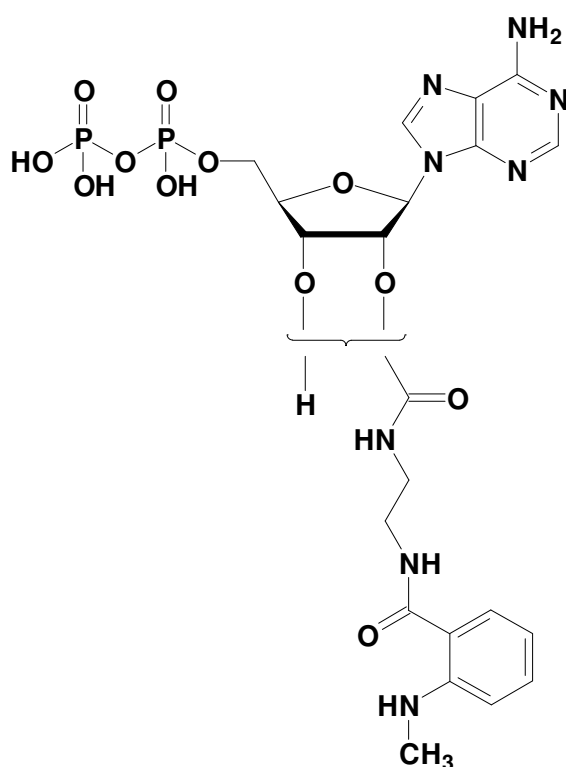


## EDA-ADP - MANT

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-diphosphate, labeled with MANT,  
 Triethylammonium salt

Cat. No.	Amount
NU-802-MNT	200 $\mu$ l (5 mM)



structural formula of EDA-ADP - MANT

**Cat. No.:** NU-802-MNT

**Molecular Formula:** C<sub>21</sub>H<sub>28</sub>N<sub>8</sub>O<sub>12</sub>P<sub>2</sub> (free acid)

**Molecular Weight:** 646.44 (free acid)

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

**Spectroscopic properties:**

$\lambda_{\text{exc}}$  335 nm;  $\lambda_{\text{em}}$  440 nm;  $\epsilon$  2,700 cm<sup>-1</sup> M<sup>-1</sup>

**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!**