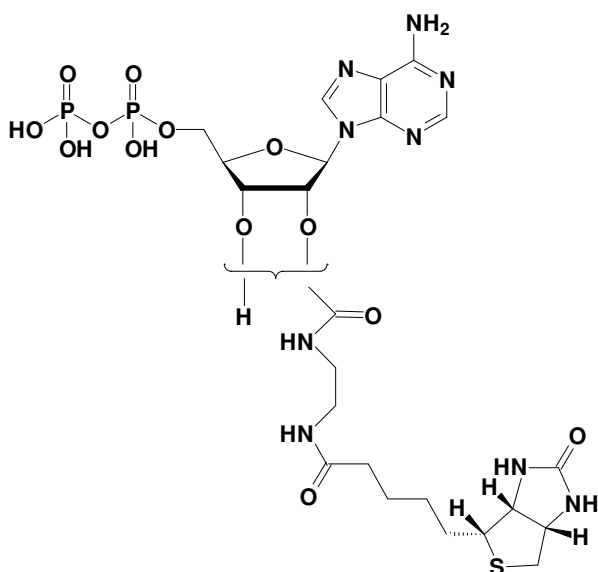


## EDA-ADP-Biotin

### 2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-diphosphate - Biotin

Cat. No.	Amount
NU-802-BIO	50 $\mu$ l / 5 mM



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

**Molecular Formula:** C<sub>24</sub>H<sub>39</sub>N<sub>9</sub>O<sub>13</sub>P<sub>2</sub>S (free acid)

**Molecular Weight:** 755.63 (free acid)

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

**Spectroscopic Properties:**  $\lambda_{\max}$  259nm;  
 $\epsilon$  15.300 M<sup>-1</sup>cm<sup>-1</sup> Tris HCl pH 7.5

**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!**