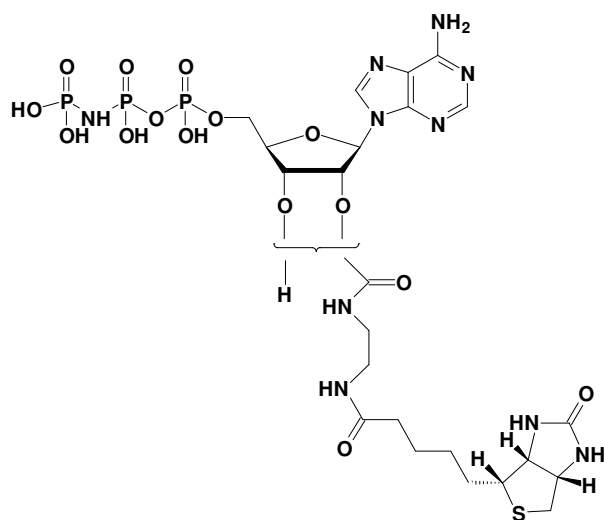


## EDA-AppNHp (EDA-AMPPNP) - Biotin

2'/3'-O-(2-Aminoethyl-carbamoyl)-adenosine-5'-[( $\beta,\gamma$ )-imido]triphosphate- Biotin,  
 Triethylammonium salt

Cat. No.	Amount
NU-810-BIO	25 $\mu$ l / 5 mM



**Cat. No.:** NU-810-BIO

**Molecular Formula:**  $C_{23}H_{37}N_{10}O_{15}P_3$  S (free acid)

**Molecular Weight:** 818.58 (free acid)

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

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

### Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at  $< -20^{\circ}C$ . If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 6 months after date of delivery.

**For research use only!**