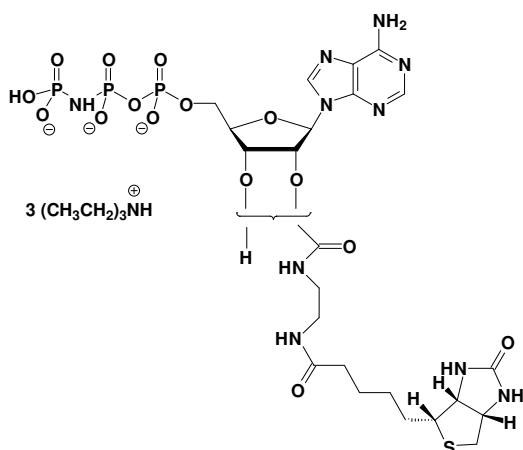


EDA-AppNHp (EDA-AMPPNP) - Biotin

2'/3'-O-(2-Aminoethyl-carbamoyl)-adenosine-5'-[(β,γ)-imido]triphosphate- Biotin,
Triethylammonium salt

Cat. No.	Amount
NU-810-BIO	25 μ l / 5 mM



Cat. No.: NU-810-BIO

Molecular Formula: $C_{23}H_{34}N_{10}O_{15}P_3$ S(Anion)

Molecular Weight: 815.56 (Anion)

Purity: > 95%, clear aqueous solution, 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 6 months after date of delivery.

For research use only!