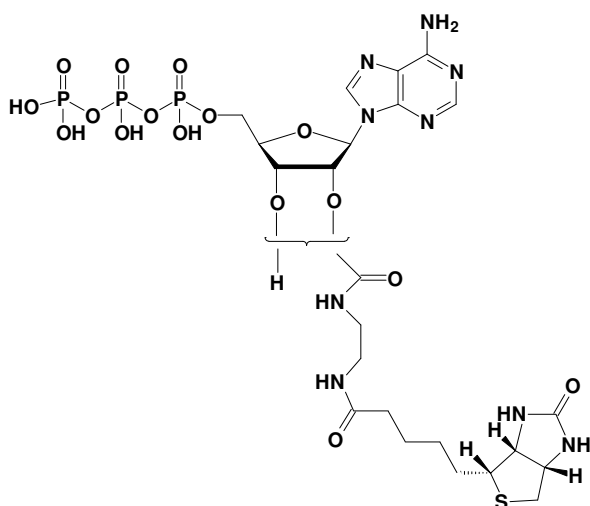


EDA-ATP - Biotin

2'/3'-O-(2-Aminoethyl-carbamoyl)-adenosine-5'-triphosphate - Biotin,
 Triethylammonium salt

Cat. No.	Amount
NU-808-BIO	50 μ l / 5 mM



Cat. No.: NU-808-BIO

Molecular Formula: $C_{23}H_{36}N_9O_{16}P_3S$ (free acid)

Molecular Weight: 819.57 (free acid)

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

Spectroscopic Properties: λ_{max} 259nm;
 ϵ 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°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

For research use only!