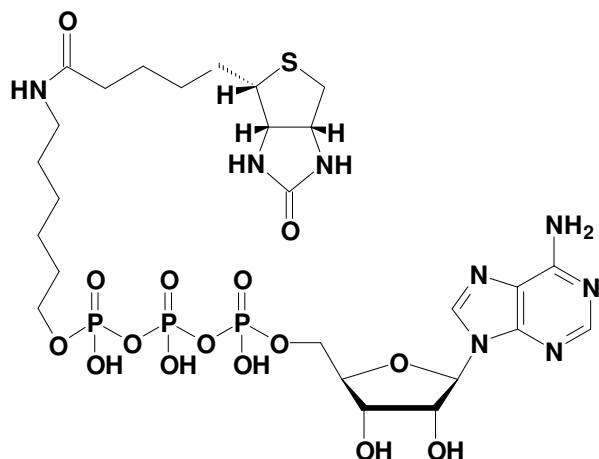


## $\gamma$ -[6-Aminoheptyl]-ATP - Biotin

$\gamma$ -[6-Aminoheptyl]-adenosine-5'-triphosphate - Biotin, Triethylammonium salt

Cat. No.	Amount
NU-833-BIO	50 $\mu$ l / 5 mM



**Cat. No.:** NU-833-BIO

**Molecular Formula:**  $C_{26}H_{43}N_8O_{15}P_3S$  (free acid)

**Molecular Weight:** 832.65 (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!**