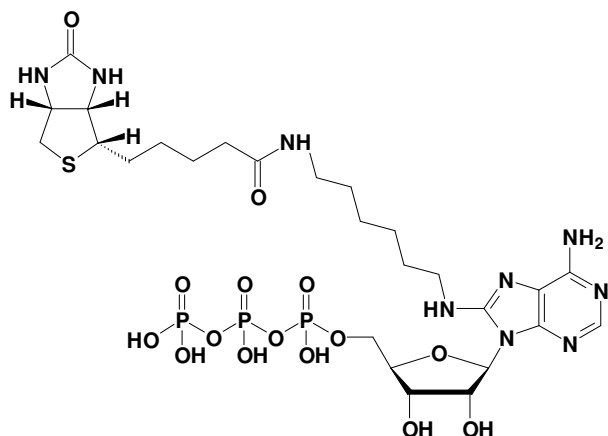


8-[(6-Amino)hexyl]-amino-ATP - Biotin

8-[(6-Amino)hexyl]-amino-adenosine-5'-triphosphate - Biotin, Triethylammonium salt

Cat. No.	Amount
NU-807-BIO	50 μ l / 5 mM



Cat. No.: NU-807-BIO

Molecular Formula: $C_{26}H_{44}N_9O_{15}P_3S$ (free acid)

Molecular Weight: 847.66 (free acid)

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 12 months after date of delivery.

For research use only!