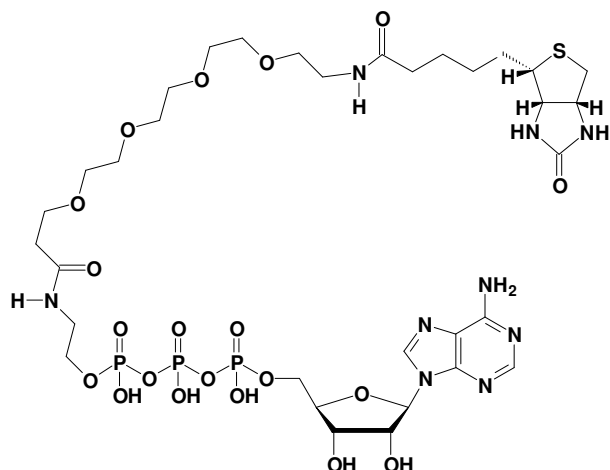


γ -[Ethyl-CONH-(PEG4)]-ATP - Biotin

γ -[Ethyl-CONH-(PEG4)]-Adenosine-5'-triphosphate - Biotin, Triethylammonium salt

| Cat. No. | Amount |
|------------|-------------------|
| NU-927-BIO | 50 μ l / 5 mM |



Cat. No.: NU-927-BIO

Molecular Formula: $C_{33}H_{56}N_9O_{20}P_3S$ (free acid)

Molecular Weight: 1023.83 (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 6 months after date of delivery.

For research use only!