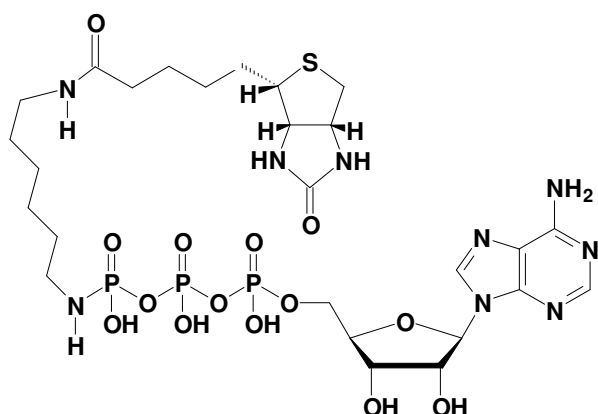


γ -[(6-Aminohexyl)-imido]-ATP - Biotin

γ -[(6-Aminohexyl)-imido]-adenosine-5'-triphosphate - Biotin, Triethylammonium salt

Cat. No.	Amount
NU-814-BIO	50 μ l / 5 mM



Cat. No.: NU-814-BIO

Molecular Formula: C₂₆H₄₄N₉O₁₄P₃S (free acid)

Molecular Weight: 831.66 (free acid)

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

Spectroscopic Properties: λ_{\max} 259 nm;
 ϵ 15,300 M⁻¹cm⁻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!