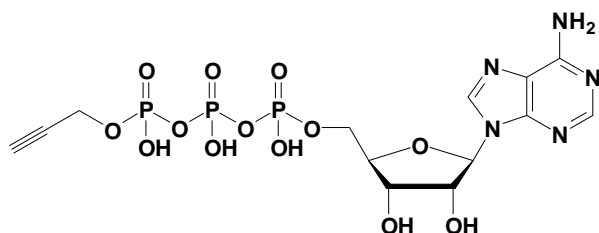


γ -Propargyl-ATP

Adenosine-5'-[γ -(propargyl)]triphosphate, Sodium salt

Cat. No.	Amount
CLK-T10-1	1 mg



Cat. No.: CLK-T10

Molecular formula: $C_{13}H_{18}N_5O_{13}P_3$ (free acid)

Molecular weight: 545.23 (free acid)

Purity: > 95% (HPLC)

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!