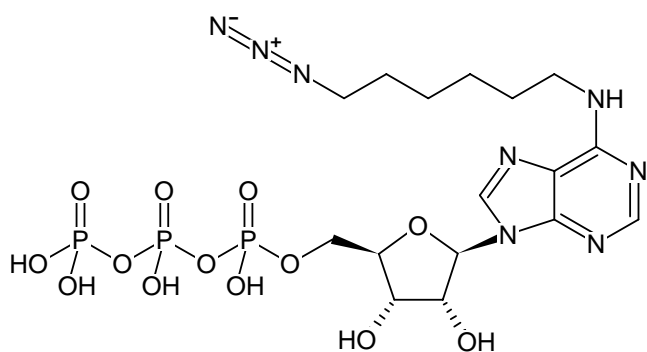




N⁶-(6-Azido)hexyl-ATP

N⁶-(6-Azido)hexyl-adenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
CLK-NU-003-1	1 mg
CLK-NU-003-5	5 mg



Structural formula of N⁶-(6-Azido)hexyl-ATP

For research use only!

Shipping: shipped on blue ice

Storage Conditions: store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: C₁₆H₂₇N₈O₁₃P₃ (free acid)

Molecular Weight: 632.35 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: lyophilised

Spectroscopic Properties: λ_{max} 259 nm; ε 15.3 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)