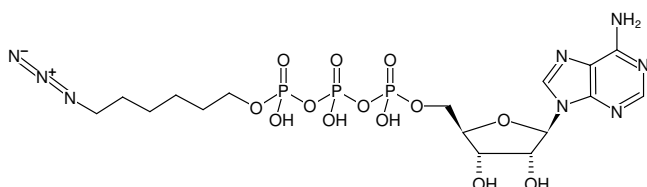




γ -(6-Azidoheptyl)-ATP

γ -(6-Azidoheptyl)-adenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1702S	50 μ l (10 mM)
NU-1702L	5 x 50 μ l (10 mM)



Structural formula of γ -(6-Azidoheptyl)-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)

Exact Mass: 632.09 g/mol (free acid)

Purity: \geq 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 10 mM - 11 mM

pH: 7.5 \pm 0.5

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