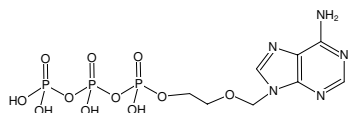


**Acyclo-ATP**

Acycloadenosine-triphosphate, Sodium salt
acyATP, Acyclo-ATP

Cat. No.	Amount
NU-538S	50 µl (100 mM)
NU-538L	5 x 50 µl (100 mM)



Structural formula of Acyclo-ATP

For general laboratory use.

Please centrifuge briefly before opening (volume ≤2 ml).

Shipping: shipped on gel packs

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: C₈H₁₄N₅O₁₁P₃ (free acid)

Molecular Weight: 449.15 g/mol (free acid)

Exact Mass: 448.99 g/mol (free acid)

Purity: ≥ 98 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 100 mM - 110 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{max} 259 nm, ε 15.1 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)