

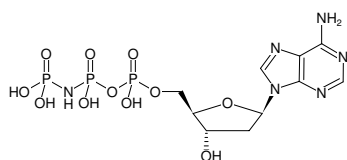
**dAppNHp**

(dAMPPNP)

dAMPPNHP

2'-Deoxyadenosine-5'-[(β,γ)-imido]triphosphate, Lithium salt

Cat. No.	Amount
NU-257	5 mg



Structural formula of dAppNHp

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: 6 months after date of delivery**Molecular Formula:** C₁₀H₁₇N₆O₁₁P₃ (free acid)**Molecular Weight:** 490.20 g/mol (free acid)**Exact Mass:** 490.02 g/mol (free acid)**Purity:** ≥ 90 % (HPLC)**Form:** white to off-white solid**Spectroscopic Properties:** λ_{max} 259 nm, ε 15.4 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Please note: For reasons of stability, please make sure that the pH value of a solution of this product never drops below 7.0. This can be achieved by dissolving the nucleotide in a buffer of your choice (50 - 100 mM, pH 7 - 10). Dissolve and adjust concentration photometrically. When stored at -20 °C, product may hydrolyse, thereby forming dAppNH₂.

The limited stability of dAppNHp makes it impossible to keep it stocked and to offer it as catalog product that is available immediately.

Therefore, dAppNHp is freshly prepared per order. For this reason the price is higher than for stable nucleotides which can be synthesised and kept in stock.