

**ApNHpp**

(AMPNPP)

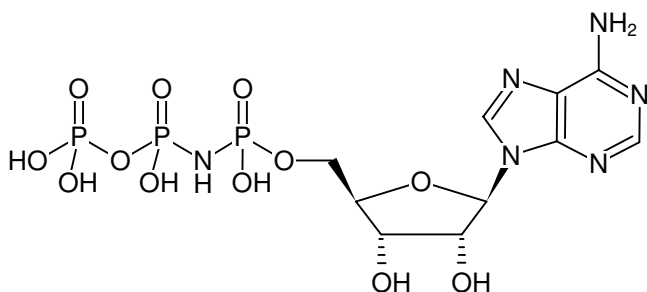
Adenosine-5'-[(α,β)-imido]triphosphate, Sodium salt

Cat. No.	Amount
NU-449-1	1 mg

Selected References:

Niland *et al.* (2021) Mechanism and Inhibition of Human Methionine Adenosyltransferase 2A. *Biochemistry* **60** (10):791-801.

Ma *et al.* (1990) Specificity of S-adenosylmethionine synthetase for ATP analogues mono- and disubstituted in bridging positions of the polyphosphate chain. *Biochemistry* **29** (6):1412.



Structural formula of ApNHpp

For research use only!**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:** C₁₀H₁₇N₆O₁₂P₃ (free acid)**Molecular Weight:** 506.20 g/mol (free acid)**Exact Mass:** 506.01 g/mol (free acid)**CAS#:** 114635-42-6 (free acid), 114635-43-7 (Tetrasodium salt)**Purity:** ≥ 95 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ_{\max} 259 nm, ϵ 15.4 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)