

**2'-Iodo-dAppNHp**

(2'-dAppNHp, 2'-I-dAMPPNP)

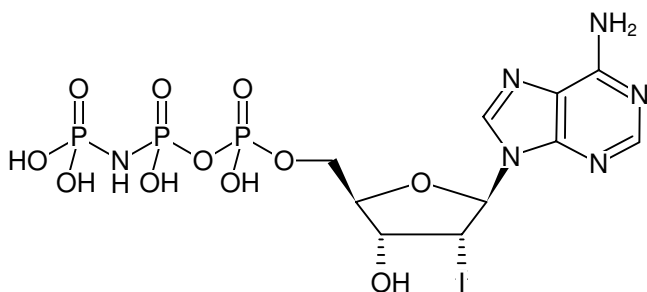
2'-Iodo-2'-deoxyadenosine-5'-[(β,γ)-imido]triphosphate, Sodium salt

Cat. No.	Amount
NU-103	200 μ l (10 mM)

Selected References:

Gruen *et al.* (1999) 2'-Halo-ATP and -GTP analogues: Rational phasing tools for protein crystallography. *Protein Sci.* **8**:2524.

Gruen *et al.* (1999) Synthesis of 2'-iodo- and 2'-bromo-ATP and GTP analogues as potential phasing tools for X-ray crystallography. *Nucleos. Nucleot.* **18** (1):137.



Structural formula of 2'-Iodo-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:** C₁₀H₁₆N₆O₁₁P₃I (free acid)**Molecular Weight:** 616.09 g/mol (free acid)**Exact Mass:** 615.91 g/mol (free acid)**CAS#:** 221092-56-4**Purity:** \geq 95 % (HPLC)**Form:** clear aqueous solution**Concentration:** 10 mM - 11 mM**pH:** 7.5 \pm 0.5**Spectroscopic Properties:** λ_{\max} 259 nm, ϵ 15.1 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.0)