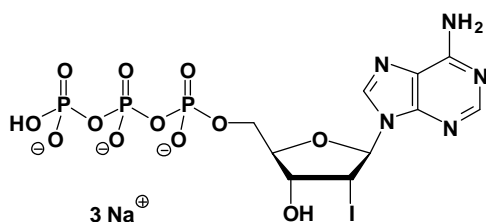


2'-I-ATP

2'-Iodo-adenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-101S	30 Units
NU-101L	150 Units



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.

Naber *et al.* (1995) A novel adenosine-triphosphate analog with a heavy-atom to target the nucleotide-binding site of proteins. *Protein Sci.* **4**:1824.

Cat. No.: NU-101

Molecular Formula: C₁₀H₁₂N₅O₁₂P₃I (Anion)

Molecular Weight: 614.05 (Anion)

Purity: > 95%, clear aqueous solution, pH 7.5

Storage conditions: Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

1 unit = 1 µl of a 10 mM solution

For research use only!