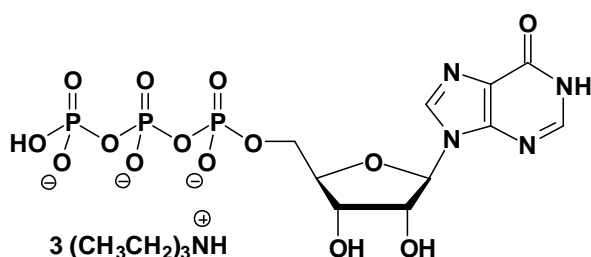


## ITP

### Inosine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-1203S	150 Units
NU-1203L	750 Units



**Cat. No.:** NU-1203

**Molecular Formula:** C<sub>10</sub>H<sub>12</sub>N<sub>4</sub>O<sub>14</sub>P<sub>3</sub> (Anion)

**Molecular Weight:** 505.14 (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.

#### For research use only!

1 unit = 1 µl of a 10 mM solution

#### Selected References:

Noji *et al.* (2001) Purine but not pyrimidine nucleotides support rotation of F(1)-ATPase. *J. Biol. Chem.* **276** (27):25480.

Bianchi *et al.* (2001) Intramolecular equilibria in metal ion complexes of guanosine 5'-triphosphate (GTP(4-)) and inosine 5'-triphosphate (ITP(4-)) in aqueous solution. *J. Inorg. Biochem.* **86** (1):148.

Chakrabarti *et al.* (2000) Nucleoside triphosphate specificity of tubulin. *Biochemistry* **39** (33):10269.

Jacob *et al.* (2000) Involvement of asparagine 118 in the nucleotide specificity of the catalytic subunit of protein kinase CK2. *FEBS Lett.* **466** (2-3):363.

Seifert *et al.* (1999) Effects of guanine, inosine, and xanthine nucleotides on beta(2)-adrenergic receptor/G(s) interactions: evidence for multiple receptor conformations. *Mol. Pharmacol.* **56** (2):348.

Wang *et al.* (1999) Identification of residues of Escherichia coli phosphofructokinase that contribute to nucleotide binding and specificity. *Biochemistry* **38** (14):4313.

Sasaki *et al.* (1998) Identification of stable RNA hairpins causing band compression in transcriptional sequencing and their elimination by use of inosine triphosphate. *Gene* **222** (1):17.

Nakahara *et al.* (1998) Inosine 5'-triphosphate can dramatically increase the yield of NASBA products targeting GC-rich and intramolecular base-paired viroid RNA. *Nucleic Acids Res.* **26** (7):1854.

Klinker *et al.* (1997) Functionally nonequivalent interactions of guanosine 5'-triphosphate, inosine 5'-triphosphate, and xanthosine 5'-triphosphate with the retinal G-protein, transducin, and with G(i)-proteins in HL-60 leukemia cell membranes. *Biochem. Pharmacol.* **54** (5):551.

Pollardknight *et al.* (1987) Kinetics of Hexokinase-D (Glucokinase) with inosine triphosphate as phosphate donor - loss of kinetic cooperativity with respect to glucose. *Biochem. J.* **245** (3):625.

Chanda *et al.* (1983) In vitro synthesis of genome length complementary RNA of vesicular stomatitis-virus in the presence of inosine 5'-triphosphate. *Virology* **129** (1):225.

Morgan *et al.* (1980) Initiation of reovirus transcription by inosine 5'-triphosphate and properties of 7-methylinosine-capped, inosinesubstituted messenger ribonucleic-acids. *Biochemistry-US* **19** (26):5960.

West (1970) Adenosine triphosphate and inosine triphosphate dependent conformational changes of adenosine diphosphate-G-actin. *Biochemistry-US* **9** (20):3847.

Azuma *et al.* (1963) Kinetic studies on hydrolysis of adenosine triphosphate and inosine triphosphate by Myosin A. *Biochim. Biophys. Acta* **73** (3):499.