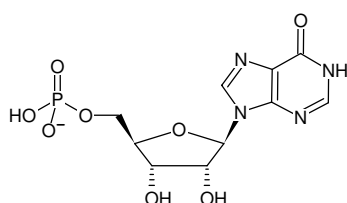


**IMP**

Inosine-5'-monophosphate, Triethylammonium salt

Cat. No.	Amount
NU-1201S	150 µl (10 mM)
NU-1201L	5 x 150 µl (10 mM)



Structural formula of IMP

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:** C₁₀H₁₃N₄O₈P (free acid)**Molecular Weight:** 348.21 g/mol (free acid)**Exact Mass:** 348.05 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ_{max} 249 nm, ε 12.2 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Selected References:**

Bhattacharya *et al.* (2000) Regiospecificity of nucleotide-amino acid mating vs. water dynamics: a key to protein-nucleic acid assemblies: structure of unidecahydrated inosine-5'-monophosphate and L-glutamic acid. *J. Chem. Crystallogr.* **30** (10):655.

Bera *et al.* (1999) Influences of cooperative hydrogen bonding economy on protein-nucleic acid complexation: Structure of unidecahydrated Inosine C5'-monophosphate and L-glutamine cocrystal at atomic resolution. *J. Chem. Crystallogr.* **29** (5):531.

Heroux *et al.* (1999) Crystal structures of the Toxoplasma gondii hypoxanthine-guanine phosphoribosyltransferase-GMP and -IMP complexes: Comparison of purine binding interactions with the XMP complex. *Biochemistry-US* **38** (44):14485.

Tullson *et al.* (1999) IMP degradative capacity in rat skeletal muscle fiber types. *Mol. Cell Biochem.* **199** (1-2): 111.

Bera *et al.* (1998) Hydrogen bonding cooperativity of water to nucleotide recognition: structure of octadecahydrated inosine 5'- monophosphate at atomic resolution. *J. Chem. Crystallogr.* **28** (7):509.

Szabados *et al.* (1998) Inosine-5'-monophosphate analogues as inhibitors of human IMP cyclohydrolase and cellular growth. *Biochem. Mol. Biol. Int.* **44** (3):617.

Mukhopadhyay *et al.* (1995) Role of water-molecules in protein-nucleic acid interactions - visualization of a model highly hydrated complex structure of inosine 5'-monophosphate and L-serine at atomic-resolution. *J. Chem. Crystallogr.* **25** (8):477.

Yu *et al.* (1993) Sulfur-containing volatile compounds generated from the interactions of inosine-5'-monophosphate and the nonvolatile flavor precursors of garlic. *Abstr. Pap. Am. Chem. S.* **206**:138, Part 1.

Greenhaff *et al.* (1993) Energy-metabolism in single human muscle-fibers during intermittent contraction with occluded circulation. *J. Physiol.-London* **460**:443.

Schiffman *et al.* (1987) Inosine-5'-monophosphate and inosine enhance some sweet tastes. *Chem. Senses* **12** (4):694.

Capparelli *et al.* (1983) The site of metal-binding in an iron (II) derivative of inosine 5'-monophosphate - an X-ray and spectral study. *FEBS Lett.* **163** (2):241.

Young *et al.* (1983) An enzymatic method for inosine 5'-monophosphate in the femtomole range. *Anal. Biochem.* **134** (1):11.