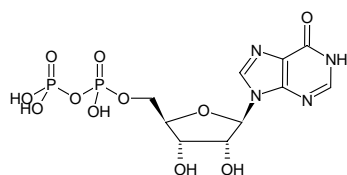


**IDP**

Inosine-5'-diphosphate, Triethylammonium salt

Cat. No.	Amount
NU-1202S	100 µl (10 mM)
NU-1202L	5 x 100 µl (10 mM)



Structural formula of IDP

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. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

Shelf Life: 12 months**Molecular Formula:** C₁₀H₁₄N₄O₁₁P₂ (free acid)**Molecular Weight:** 428.18 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution, pH 7.5 ± 0.5**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:**

AlAli *et al.* (1996) Effects of metal ions on the activity of cytosolic phosphoenolpyruvate carboxykinase from camel kidney. *Arab. Gulf J. Sci. Res.* **14** (3):535.

Vial *et al.* (1995) Purification, partial kinetic characterization and reactive sulfhydryl-groups of the phosphoenolpyruvate carboxykinase from peromyscus-purpuratus adductor muscle. *Comp. Biochem. Phys. B.* **112** (3):451.

Okkuluri *et al.* (1994) Peparation, characterization and crystallization of an antibody fab fragment that recognizes RNA -crystal-structures of native fab and 3 fab-monomucleotide complexes. *J. Mol. Biol.* **243** (2):283.

Jetten *et al.* (1994) Regulation of phospho (enol)-pyruvate and oxaloacetate-converting enzymes in corynebacteriumglutamicum. *Appl. Microbiol. Biot.* **41** (1):47.

Weng *et al.* (1994) Hydrogen-bond interactions of g-proteins with the guanine ring moiety of guanine-nucleotides. *Protein Sci.* **3** (1):22.

Vicente *et al.* (1991) Differentiation between several types of phosphohydrolases in light microsomes of corn roots. *Plant Physiol.* **96** (4):1345.

Tung *et al.* (1991) On the mechanism of nucleotide diphosphate activation of the ATP-sensitive k⁺ channel in ventricular cell of guinea-pig. *J. Physiol.-London* **437**:239.

Schramm *et al.* (1971) Studies on allosteric modification of nucleoside diphosphatase activity by magnesium nucleoside triphosphates and inosine diphosphate. *Biochemistry-US* **10** (12):2272.

Oklik (1955) The formation of nucleoside triphosphate from inosine diphosphate in yeast. *Biochim. Biophys. Acta* **16** (4):610.