

Inosine Kit

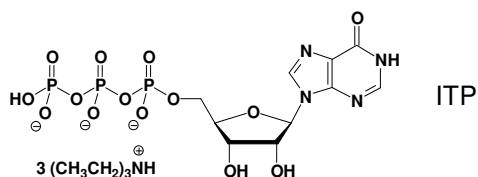
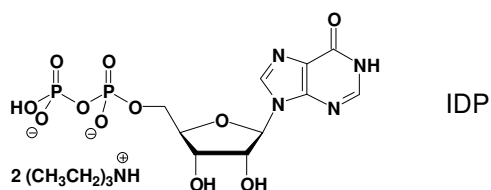
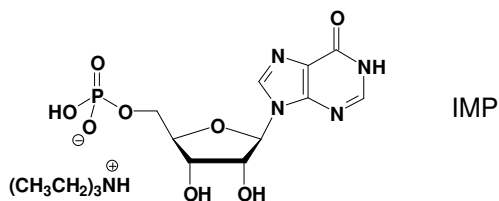
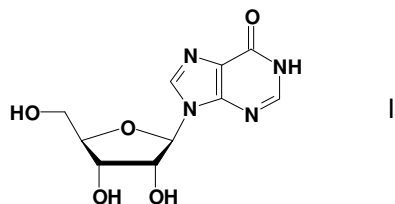
Cat.No.: NK-115

Kit Contents

Inosine Analogs	Cat. No.	Amount
I	N-1049	5 mg
IMP	NU-1201	200 Units
IDP	NU-1202	200 Units
ITP	NU-1203	200 Units

1 Unit = 1 μ l of a 10 mM solution

Structures



Introduction

Inosines are widely used for mechanistic studies on ATP (cATP) or GTP (cGTP) binding proteins.

Kit Description

The Inosine Kit contains a set of 4 typical Inosine analogs (Nucleoside, 5'-Mono,-Di and Triphosphate).

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 (ITP4-) 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.

Storage and Stability

Short term exposure (up to 1 week cumulative) to ambient temperature is possible. Long term storage is recommended at < -20°C. If properly stored, Jena Bioscience guarantees optimal performance of the compounds for 12 months after date of delivery.