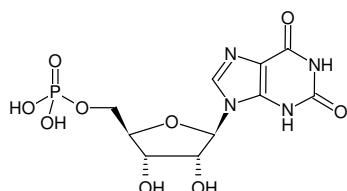


**XMP**

Xanthosine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-603S	150 µl (10 mM)
NU-603L	5 x 150 µl (10 mM)



Structural formula of XMP

**For general laboratory use.****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:** C<sub>10</sub>H<sub>13</sub>N<sub>4</sub>O<sub>9</sub>P (free acid)**Molecular Weight:** 364.21 g/mol (free acid)**Exact Mass:** 364.04 g/mol (free acid)**CAS#:** 25899-70-1**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<sub>max</sub> 276 nm, ε 9.6 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)**Selected References:**

Fulga *et al.* (2001) SR beta coordinates signal sequence release from SRP with ribosome binding to the translocon. *EMBO J.* **20** (9):2338.

Legate *et al.* (2000) Nucleotide-dependent binding of the GTPase domain of the signal recognition particle receptor betasubunit to the alpha-subunit. *J. Biol. Chem.* **275** (35):27439.

Sakash *et al.* (2000) The use of nucleotide analogs to evaluate the mechanism of the heterotropic response of Escherichia coli aspartate transcarbamoylase. *Protein Sci.* **9** (1):53.

Hwang *et al.* (1999) Structure-based identification of a novel NTPase from Methanococcus jannaschii. *Nat. Struct. Biol.* **6** (7):691.

Muraoka *et al.* (1999) Effects of purinenucleotide analogues on microtubule assembly. *Cell Struct. Funct.* **24** (5):305.

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.

Yu *et al.* (1997) Characterization of a Go alpha mutant that binds xanthine nucleotides. *J. Biol. Chem.* **272** (29):18015.

Rybin *et al.* (1996) GTPase activity of Rab5 acts as a timer for endocytic membrane fusion. *Nature* **383** (6597):266.

Powers *et al.* (1995) Reciprocal stimulation of GTP hydrolysis by 2 directly interacting GTPases. *Science* **269** (5229):1422.