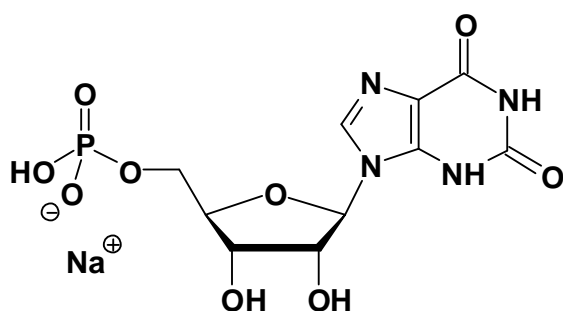


XMP

Xanthosine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-603S	150 Units
NU-603L	750 Units



Cat. No.: NU-603

Molecular Formula: C₁₀H₁₂N₄O₉P (Anion)

Molecular Weight: 363.20 (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:

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.

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.

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

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.