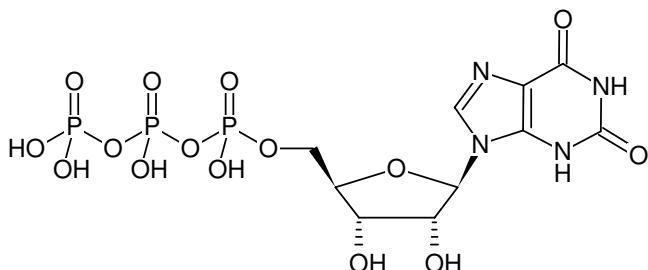




XTP - Sodium salt

Xanthosine-5'-triphosphate, Sodium salt

| Cat. No. | Amount |
|-----------|--------|
| NU-935-2 | 2 mg |
| NU-935-10 | 10 mg |



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: C₁₀H₁₅N₄O₁₅P₃ (free acid)

Molecular Weight: 524.16 g/mol (free acid)

Exact Mass: 523.97 g/mol (free acid)

CAS#: 105931-36-0

Purity: ≥ 95 % (HPLC)

Form: solid

Color: white to off-white

Spectroscopic Properties: λ_{max} 276 nm, ε 9.6 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Applications:

Inhibitor of adenylyl and guanylyl cyclase^[1]

Specific Ligands:

Soluble guanylyl cyclase^[2]

Related Products:

XTP, Triethylammonium salt #NU-602

Selected References:

[1] Spangler et al. (2011) Interaction of the diguanylate cyclase YdeH of Escherichia coli with 2', (3')-substituted purine and pyrimidine nucleotides. *J. Pharmacol. Exp. Ther.* **336** (1):234.

[2] Chang et al. (2005) Nitric Oxide-dependent Allosteric Inhibitory Role of a Second Nucleotide Binding Site in Soluble Guanylyl Cyclase. *J. Biol. Chem.* **280** (12):11513.

Beste et al. (2012) Nucleotidyl cyclase activity of soluble guanylyl cyclase α1β1. *Biochemistry* **51** (1):194.

Spangler et al. (2011) Interaction of the diguanylate cyclase YdeH of Escherichia coli with 2', (3')-substituted purine and pyrimidine nucleotides. *J. Pharmacol. Exp. Ther.* **336** (1):234.

Fulga et al. (2001) SR beta coordinates signal sequence release from SRP with ribosome binding to the translocon. *EMBO J.* **20** (9):238.

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.