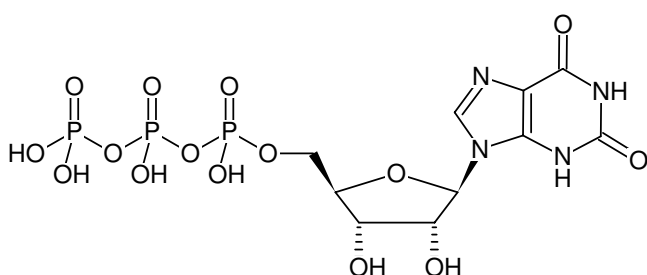




## XTP - Triethylammonium salt

Xanthosine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-602S	15 µl (100 mM)
NU-602L	5 x 15 µl (100 mM)



Structural formula of XTP - Triethylammonium salt

### For research use only!

**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>15</sub>N<sub>4</sub>O<sub>15</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 524.17 g/mol (free acid)

**Exact Mass:** 523.97 g/mol (free acid)

**CAS#:** 6253-56-1

**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

**Color:** colorless to slightly yellow

**Concentration:** 100 mM - 110 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)

### Applications:

Inhibitor of adenylyl and guanylyl cyclase<sup>[1]</sup>

### Specific Ligands:

Soluble guanylyl cyclase<sup>[2]</sup>

### Related Products:

XTP, Sodium salt #NU-935

### 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):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.