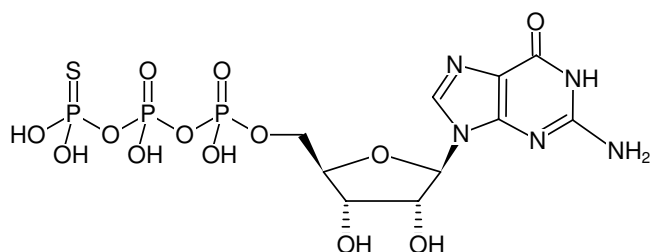


**GTPyS**

Guanosine-5'-(γ-thio)-triphosphate, Tetralithium salt  
 Guanosine-5'-(3-thio)-triphosphate, Guanosine-5'-(3-thiotriphosphate)

Cat. No.	Amount
NU-412-2	2 mg
NU-412-10	10 mg
NU-412-20	20 mg



Structural formula of GTPyS

**For research use only!****Shipping:** shipped on dry ice**Storage Conditions:** store at -20 °C**Shelf Life:** 6 months after date of delivery**Molecular Formula:** C<sub>10</sub>H<sub>16</sub>N<sub>5</sub>O<sub>13</sub>P<sub>3</sub>S (free acid)**Molecular Weight:** 539.25 g/mol (free acid)**Exact Mass:** 538.97 g/mol (free acid)**CAS#:** 94825-44-2**Purity:** ≥ 90 % (HPLC), contains < 10 % GDP**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ<sub>max</sub> 252 nm, ε 13.7 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)**Applications:**Functional assay for GPCRs<sup>[1, 2]</sup>Binding assay<sup>[3]</sup>**Specific Ligands:**Heterotrimeric and monomeric G-proteins<sup>[4]</sup>Phosphodiesterase PDE-6<sup>[5]</sup>for P2Y-like receptor<sup>[6]</sup> and for P2Y<sub>13</sub> receptor<sup>[7]</sup>

**Please note:** For reasons of stability, please make sure that the pH value of a solution of this product never drops below 7.0. This can be achieved by dissolving the nucleotide in a buffer of your choice (50 - 100 mM, pH 7 - 10). Dissolve and adjust concentration photometrically.

**Selected References:**

- [1] Cascio *et al.* (2010) In vitro and in vivo pharmacological characterization of two novel selective cannabinoid CB (2) receptor inverse agonists. *Pharmacological research* **61**:349.
- [2] Obara *et al.* (2010) Agonist-dependent attenuation of micro-opioid receptor-mediated G-protein activation in the dorsal root ganglia of neuropathic rats. *J. Neural Transmission* **117**:421.
- [3] Cascio *et al.* (2010) Evidence that the plant cannabinoid cannabigerol is a highly potent alpha2-adrenoreceptor agonist and moderately potent SHT1A receptor antagonist. *Brit. J. Pharmacol.* **159**:129.
- [4] Koval *et al.* (2010) Europium-labeled GTP as a general nonradioactive substitute for [35S]GTPyS in high-throughput G-protein studies. *Analytical Biochem.* **397**:202.
- [5] Yamazaki *et al.* (2010) Mechanism for the regulation of mammalian cGMP phosphodiesterase 6. 2: Isolation and characterization of the transducin-activated form. *Mol. Cell. Biochem.* **339**:235.
- [6] Daniele *et al.* (2011) Agonist-induced desensitization/resensitization of human G protein-coupled receptor 17: a functional cross-talk between purinergic and cysteinyl-leukotriene ligands. *J. Pharmacol. Exp. Ther.* **338** (2):559.
- [7] Marteau *et al.* (2003) Pharmacological characterization of human P2Y<sub>13</sub> receptor. *Mol. Pharmacol.* **64** (1):104.
- Schmoehl *et al.* (2012) Functional analysis of Rho GTPase activation and inhibition in a bead-based miniaturized format. *Methods. Mol. Biol.* **827**: 271.
- Martin-Galiano AJ *et al.* (2011) Bacterial tubulin distinct loop sequences and primitive assembly properties support its origin from a eukaryotic tubulin ancestor. *J. Bio. Chem.* **286** (22):19789.

**GTPγS**

Guanosine-5'-(γ-thio)-triphosphate, Tetralithium salt  
Guanosine-5'-(3-thio)-triphosphate, Guanosine-5'-(3-thiotriphosphate)

Abenza *et al.* (2010) *Aspergillus* RabB Rab5 integrates acquisition of degradative identity with the long distance movement of early endosomes. *Mol. Biol. Cell.* **21** (15):2756.

Beuve *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.

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.

Kirilenko *et al.* (2002) GTP-induced membrane binding and ion channel activity of annexin VI: Is annexin VI a GTP biosensor? *Biophys. J.* **82** (5):2737.

Shiba *et al.* (2002) Hypocretin stimulates [ (35)]GTP gamma S binding in Hcrt2-transfected cell lines and in brain homogenate. *Biochem. Biophys. Res. Commun.* **294** (3):615.

Aiyar *et al.* (1994) Human At (1) receptor is a single-copy gene - characterization in a stable cell-line. *Mol. Cell. Biochem.* **131** (1):75.

Ashktorab *et al.* (1994) Presence of GTP-binding proteins in the plasma-membrane of the *Phycomyces* sporangiophore. *Exp. Mycol.* **18** (2):139.

Schwemmler *et al.* (1994) The interferon-induced 67-kDa Guanylate-binding protein (HGbp1) is a GTPase that converts GTP to GMP. *J. Biol. Chem.* **269** (15):11299.

Noel *et al.* (1993) The 2.2 Å crystal structure of transducin- alpha complexed with GTP gamma S. *Nature* **366**:654.