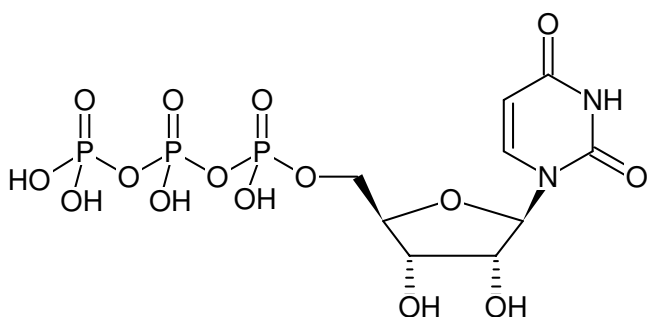


**UTP - Solution**

100 mM Sodium salt solution  
Uridine 5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1013	1 ml (100 mM)
NU-1013-100ML	100 ml (100 mM)



Structural formula of UTP - Solution

**For general laboratory use.**

**Shipping:** shipped on gel packs

**Storage Conditions:** store at -20 °C

**Shelf Life:** 12 months

**Molecular Formula:** C<sub>9</sub>H<sub>15</sub>N<sub>2</sub>O<sub>15</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 484.14 g/mol (free acid)

**CAS#:** 19817-92-6

**Purity:** ≥ 99 % (HPLC)

**Form:** clear aqueous solution

**Concentration:** 100 mM ±2 %

**pH:** 8.0 ±0.2 (22 °C)

**Spectroscopic Properties:** λ<sub>max</sub> 262 nm, ε 9.8 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.0)

**Applications:**

Activation of purinergic receptors<sup>[1,2,3,4]</sup>

Cardioprotection against hypoxic damage<sup>[2]</sup>

Enzyme kinetic parameters<sup>[5]</sup>

Phosphorylation of EGF-receptor via purinergic receptors<sup>[3]</sup>

Stimulation of neurogenesis and dopaminergic neurons<sup>[6]</sup>

**Description:**

Ultrapure UTP supplied as clear aqueous solution.

**Specific Ligands:**

Enterovirus 71 3D RNA polymerase<sup>[7]</sup>

**Ligand for purinergic receptors:**

P2X<sub>1</sub><sup>[4]</sup>  
P2Y<sub>2</sub><sup>[7,8,9,10]</sup>  
P2Y<sub>4</sub><sup>[8,10,11]</sup>  
P2Y<sub>6</sub><sup>[8]</sup>

**Selected References:**

- [1] Raqeeb *et al.* (2011) Purinergic P2Y2 receptors mediate rapid Ca<sup>2+</sup> mobilization, membrane hyperpolarization and nitric oxide production in human vascular endothelial cells. *Cell Calcium* **49**:240.
- [2] Golan *et al.* (2011) Extracellular nucleotide derivatives protect cardiomyocytes against hypoxic stress. *Biochemical Pharmacology* **81**:1219.
- [3] Boucher *et al.* (2011) Distinct activation of epidermal growth factor receptor by UTP contributes to epithelial cell wound repair. *American Journal Pathology* **178**:1092.
- [4] Sugihara *et al.* (2011) Dual signaling pathway of arterial constriction by extracellular uridine-5-triphosphate in the rat. *J. Pharmacological Sciences (Japan)* **115**:293.
- [5] Ma *et al.* (2011) Molecular cloning and analysis of the UDP-glucose pyrophosphorylase in *Streptococcus equi* subsp. *Zooepidemicus*. *Molecular Biology Reports* **38**:2751.
- [6] Delic *et al.* (2011) Nucleotides affect neurogenesis and dopaminergic differentiation of mouse fetal midbrain-derived neural precursor cells. *Purinergic Signalling* **6**:417.
- [7] Jiang *et al.* (2011) Biochemical characterization of enterovirus 71 3D RNA polymerase. *Biochim. Biophys. Acta, Gene Regulatory Mechanisms* **1809**:211.
- [8] Pendergast *et al.* (2001) Synthesis and P2Y receptor activity of a series of uridine dinucleoside 5'-polyphosphates. *Bioorg. Med. Chem. Lett.* **11** (2):157.

**UTP - Solution**

100 mM Sodium salt solution

Uridine 5'-triphosphate, Sodium salt

[9] Shaver *et al.* (1997) 4-substituted uridine 5'-triphosphates as agonists of the P2Y2 purinergic receptor. *Nucleosides and Nucleotides* **16 (7)**:1099.

[10] Kim *et al.* (2002) Methanocarba modification of uracil and adenine nucleotides: High potency of northern ring conformation at P2Y1, P2Y2, P2Y4 and P2Y11 but not P2Y6 receptors. *J. Med. Chem.* **45**:208.

[11] Nguyen *et al.* (1995) Cloning, expression, and chromosomal localization of human uridine nucleotide receptors. *J. Biol. Chem.* **270 (52)**:30845.