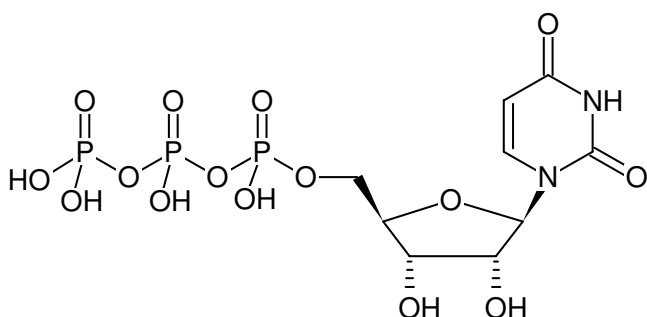


**UTP - Solution**

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

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



Structural formula of UTP - Solution

For in vitro 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. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

Shelf Life: 12 months**Molecular Formula:** C₉H₁₅N₂O₁₅P₃ (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_{max} 262 nm, ε 9.8 L mmol⁻¹ cm⁻¹ (pH 7.0)**Applications:**Activation of purinergic receptors^[1,2,3,4]Cardioprotection against hypoxic damage^[2]Enzyme kinetic parameters^[5]Phosphorylation of EGF-receptor via purinergic receptors^[3]Stimulation of neurogenesis and dopaminergic neurons^[6]**Description:**

Ultrapure UTP supplied as clear aqueous solution.

Specific Ligands:Enterovirus 71 3D RNA polymerase^[7]**Ligand for purinergic receptors:**P2X₁^[4]P2Y₂^[7,8,9,10]P2Y₄^[8,10,11]P2Y₆^[8]**Selected References:**

[1] Raqeeb *et al.* (2011) Purinergic P2Y2 receptors mediate rapid Ca²⁺ 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. *Zoepidemicus*. *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.