

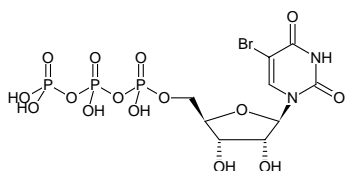
**5-Bromo-UTP**

(5Br-UTP)

5-BrUTP

5-Bromo-uridine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-121S	50 µl (10 mM)
NU-121L	5 x 50 µl (10 mM)



Structural formula of 5-Bromo-UTP

**For research use only!****Shipping:** shipped on blue ice**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:** C<sub>9</sub>H<sub>14</sub>N<sub>2</sub>O<sub>15</sub>P<sub>3</sub>Br (free acid)**Molecular Weight:** 563.03 g/mol (free acid)**CAS#:** 161848-60-8**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution, pH 7.5 ±0.5**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ<sub>max</sub> 278 nm; ε 9.7 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)**Applications:**Agonist for P2Y<sub>2</sub>, P2Y<sub>4</sub> and P2Y<sub>6</sub> receptors<sup>[1,2,3]</sup>**Selected References:**

[1] Wilson *et al.* (1998) Nucleotide-evoked calcium signals and anion secretion in equine cultured epithelia that express apical P2Y (2) receptors and pyrimidine nucleotide receptors. *Brit. J. Pharmacol.* **124** (4):832.

[2] Communi *et al.* (1996) Pharmacological characterization of the human P-2Y<sub>4</sub> receptor. *Eur. J. Pharmacol.* **317** (2-3):383.

[3] El-Tayeb *et al.* (2006) Synthesis and structure-activity relationships of uracil nucleotide derivatives and analogues as agonists at human P2Y<sub>2</sub>, P2Y<sub>4</sub>, and P2Y<sub>6</sub> receptors. *J. Med. Chem.* **49** (24):7076.

Gilles *et al.* (2007) Regulatory Mechanisms Differ in UMP Kinase from Gram-negative and Gram-positive Bacteria. *J. Biol. Chem.* **282** (10):7242.

Dunoyer *et al.* (2001) Peanut clump virus RNA-1-Encoded p15 regulates viral RNA accumulation but is not abundant at viral RNA replication sites. *J. Virology* **75** (4):1941.

Cmarko *et al.* (2000) Ultrastructural analysis of nucleolar transcription in cells microinjected with 5-bromo-UTP. *Histochem. Cell Biol.* **113** (3):181.

Cmarko *et al.* (1999) Ultrastructural analysis of transcription and splicing in the cell nucleus after bromo-UTP microinjection. *Mol. Biol. Cell* **10** (1):211.

Keene *et al.* (1999) Initially transcribed sequences strongly affect the extent of abortive initiation by RNA polymerase II. *J. Biol. Chem.* **274** (17):11526.

Wilson *et al.* (1999) Pyrimidine nucleotide-evoked inhibition of cyclic amp accumulation in equine epithelial cells. *Exp. Physiol.* **84** (4):639.

Mote *et al.* (1998) Recognition of a human arrest site is conserved between RNA polymerase II and prokaryotic RNA polymerases. *J. Biol. Chem.* **273** (27):16843.

Schul *et al.* (1998) Coiled bodies and U2 snRNA genes adjacent to coiled bodies are enriched in factors required for snRNA transcription. *Mol. Biol. Cell* **9** (5):1025.

Tanner *et al.* (1988) Binding interactions between yeast tRNA ligase and a precursor transfer ribonucleic acid containing two photoreactive uridine analogues. *Biochemistry* **27** (24):8852.