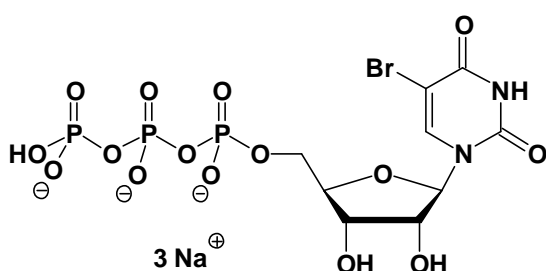


## 5Br-UTP

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

Cat. No.	Amount
NU-121S	50 Units
NU-121L	250 Units



**Cat. No.:** NU-121

**Molecular Formula:**  $C_9H_{11}N_2O_{15}P_3Br$  (Anion)

**Molecular Weight:** 560.01 (Anion)

**Purity:** > 95%, clear aqueous solution, pH 7.5

### Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

### For research use only!

1 unit = 1  $\mu$ l of a 10 mM solution

### Selected References:

Gilles *et al.* (2007) Regulatory Mechanism 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.

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

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

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

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.

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.

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.

Communi *et al.* (1996) Pharmacological characterization of the human P-2Y4 receptor. *Eur. J. Pharmacol.* **317** (2-3):383.

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.