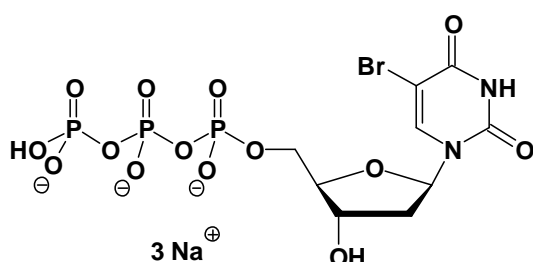


5Br-dUTP

5-Bromo-2'-deoxy-uridine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-122S	50 Units
NU-122L	250 Units



Cat. No.: NU-122

Molecular Formula: $C_9H_{11}N_2O_{14}P_3Br$ (Anion)

Molecular Weight: 544.01 (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 μ l of a 10 mM solution

Selected References:

Ji *et al.* (1991) T4-phage ribonucleotide reductase - allosteric regulation in vivo by thymidine triphosphate. *J. Biol. Chem.* **266** (25):16289.

Srinivasan, L. *et al.* (2002) *Eur. J. Biochem.* **269**:1780

Li, X. *et al.* (1995) *Cell Prolif.* **28-11**:571

Ekstrand, D. H. (1996) *Biotechnol Appl. Biochem.* **23-2**:95

Gorczyca, W. (1996) *Cell Prolif.* **29-10**:539

Mote, J. *et al.* (1998) *J. Biological Chemistry* **273-27**:16843

Pal, M. *et al.* (2002) *Molecular and Cellular Biology* **22-1**:30