

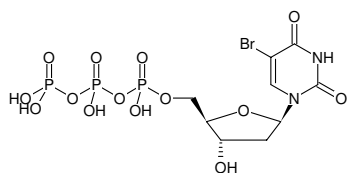
**5-Bromo-dUTP**

(5Br-dUTP)

5-BrdUTP

5-Bromo-2'-deoxyuridine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-122S	50 µl (10 mM)
NU-122L	5 x 50 µl (10 mM)



Structural formula of 5-Bromo-dUTP

For research 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.

Shelf Life: 12 months after date of delivery**Molecular Formula:** C₉H₁₄N₂O₁₄P₃Br (free acid)**Molecular Weight:** 547.04 g/mol (free acid)**Exact Mass:** 545.88 g/mol (free acid)**CAS#:** 102212-99-7**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 278 nm, ε 9.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Selected References:**

Ma *et al.* (2008) The mutagenic properties of BrdUTP in a random mutagenesis process. *Mol. Biol. Rep.* **35**:663.

Pal *et al.* (2002) Strong natural pausing by RNA polymerase II within 10 bases of transcription start may result in repeated slippage and reextension of the nascent RNA. *Mol. Cell Biol.* **22** (1):30.

Srinivasan *et al.* (2002) Characterization of RNA polymerase III transcription factor TFIIIC from the mulberry silkworm, *Bombyx mori*. *Eur. J. Biochem.* **269**(6):1780.

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.

Ekstrand (1996) A sensitive assay for the quantification of reverse transcriptase activity based on the use of carrier-bound template and non-radioactive-product detection, with special reference to human-immunodeficiency-virus isolation. *Biotechnol Appl. Biochem.* **23**-2:95.

Gorczyca (1996) Laser scanning cytometer (LSC) analysis of fraction of labelled mitoses (FLM). *Cell Prolif.* **29** (10):539.

Li *et al.* (1995) Labelling DNA strand breaks with BrdUTP. Detection of apoptosis and cell proliferation. *Cell Prolif.* **28** (11):571.

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