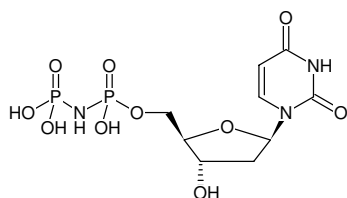


**dUpNHp**

(dUMPNP)

2'-Deoxyuridine-5'-[( $\alpha,\beta$ )-imido]diphosphate, Sodium salt

Cat. No.	Amount
NU-902S	100 $\mu$ l (10 mM)
NU-902L	5 x 100 $\mu$ l (10 mM)



Structural formula of dUpNHp

**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.

**Shelf Life:** 12 months after date of delivery**Molecular Formula:** C<sub>9</sub>H<sub>15</sub>N<sub>3</sub>O<sub>10</sub>P<sub>2</sub> (free acid)**Molecular Weight:** 387.17 g/mol (free acid)**CAS#:** 178809-71-7**Purity:**  $\geq$  95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5  $\pm$  0.5**Spectroscopic Properties:**  $\lambda_{\max}$  262 nm,  $\epsilon$  10.2 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)**Selected References:**

Kovári *et al.* (2008) Methylene substitution at the  $\alpha,\beta$  bridging position within the phosphate chain of dUDP profoundly perturbs ligand accommodation into the dUTPase active site. *J. Biol. Chem.* **283**:308.

Barabás *et al.* (2004) Structural Insights into the Catalytic Mechanism of Phosphate Ester Hydrolysis by dUTPase. *J. Biol. Chem.* **279** (41):42907.

Persson *et al.* (1996) Synthesis of 2'-deoxyuridine 5'-( $\alpha,\beta$ -imido)triphosphate: A substrate analogue and potent inhibitor of dUTPase. *Bioorgan. Med. Chem.* **4** (4):553.