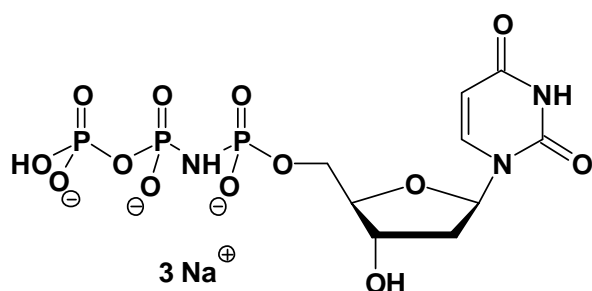


## dUpNHpp (dUMPNPP)

2'-Deoxy-uridine-5'-[( $\alpha,\beta$ )-imido]triphosphate, Sodium salt

Cat. No.	Amount
NU-903S	50 Units
NU-903L	250 Units



**Cat. No.:** NU-903

**Molecular Formula:** C<sub>9</sub>H<sub>13</sub>N<sub>3</sub>O<sub>13</sub>P<sub>3</sub> (Anion)

**Molecular Weight:** 464.13 (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 6 months after date of delivery.

### For research use only!

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

### Selected References:

Vertessy *et al.* (2007) Flexible segments modulate co-folding of dUTPase and nucleocapsid proteins. *Nucleic Acids Research* **35** (2):495.

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.