

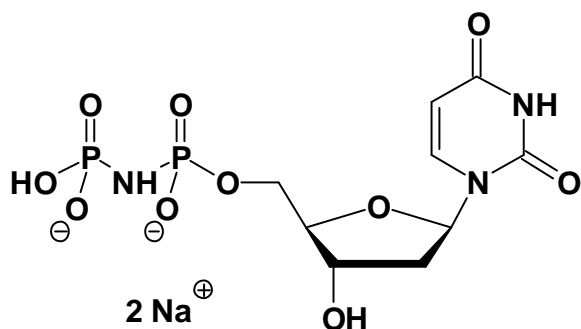
## dUpNHp (dUMPNP)

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

Cat. No.	Amount
NU-902S	100 Units
NU-902L	500 Units

### Selected References:

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.



**Cat. No.:** NU-902

**Molecular Formula:** C<sub>9</sub>H<sub>13</sub>N<sub>3</sub>O<sub>10</sub>P<sub>2</sub> (Anion)

**Molecular Weight:** 385.16 (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