

**UpNHpp**

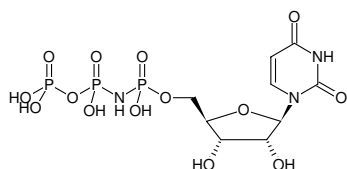
(UMPNPP)

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

Cat. No.	Amount
NU-930S	50 $\mu$ l (10 mM)
NU-930L	5 x 50 $\mu$ l (10 mM)

**Selected References:**

Kroupova *et al.* (2019) Structural basis for acceptor RNA substrate selectivity of the 3' terminal uridylyl transferase Tailor. *Nucleic Acids Res.* **47** (2):1030.



Structural formula of UpNHpp

**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>16</sub>N<sub>3</sub>O<sub>14</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 483.15 g/mol (free acid)**Exact Mass:** 482.98 g/mol (free acid)**Purity:**  $\geq$  95 % (HPLC)**Form:** clear aqueous solution**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)