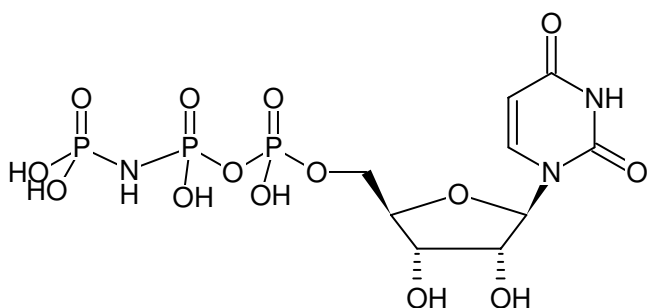


**UppNHp - Solid**

(UMPPNP)

Uridine-5'-[(β,γ)-imido]triphosphate, Sodium salt

Cat. No.	Amount
NU-415-5	5 mg



Structural formula of UppNHp - Solid

**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:** 6 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)**CAS#:** 82145-58-2**Purity:** ≥ 95 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ<sub>max</sub> 262 nm, ε 10.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)