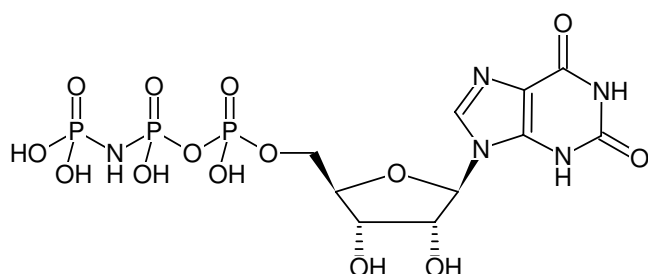


**XppNHp**

(XMPPNP)

Xanthosine-5'-[[β,γ]-imido]triphosphate, Sodium salt

Cat. No.	Amount
NU-996	5 mg



Structural formula of XppNHp

**For general laboratory use.****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>10</sub>H<sub>16</sub>N<sub>5</sub>O<sub>14</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 523.18 g/mol (free acid)**Exact Mass:** 522.99 g/mol (free acid)**CAS#:** 187533-43-3**Purity:** ≥ 90 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ<sub>max</sub> 276 nm, ε 9.6 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)**Please note:**

Due to the limited stability, the price is higher than for stable nucleotides which can be kept in stock. We strongly suggest always to prepare fresh solutions for immediate use. For this, neutral to slightly alkaline buffers and low temperatures are required.

**Selected References:**

Tomar *et al.* (2009) Distinct GDP/GTP bound states of the tandem G-domains of EngA regulate ribosome binding. *Nucleic Acids Research* **37** (7):2359.

Yanachkov *et al.* (1997) Synthesis and effect of nonhydrolyzable xanthosine triphosphate derivatives on prenylation of Rab5 (D136N). *Mol. Pharmacol.* **51** (1):47.