

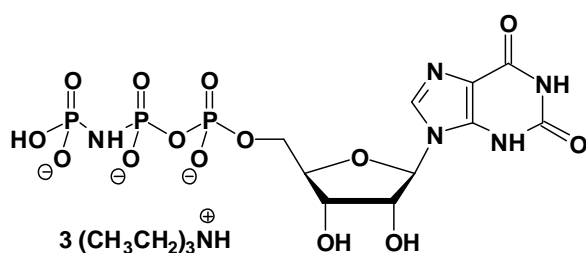
XppNHp (XMPPNP)

Xanthosine-5'-[(β , γ)-imido]triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-403S	50 Units
NU-403L	250 Units

Selected References:

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



Cat. No.: NU-403

Molecular Formula: C₁₀H₁₃N₅O₁₄P₃ (Anion)

Molecular Weight: 520.15 (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 μ l of a 10 mM solution