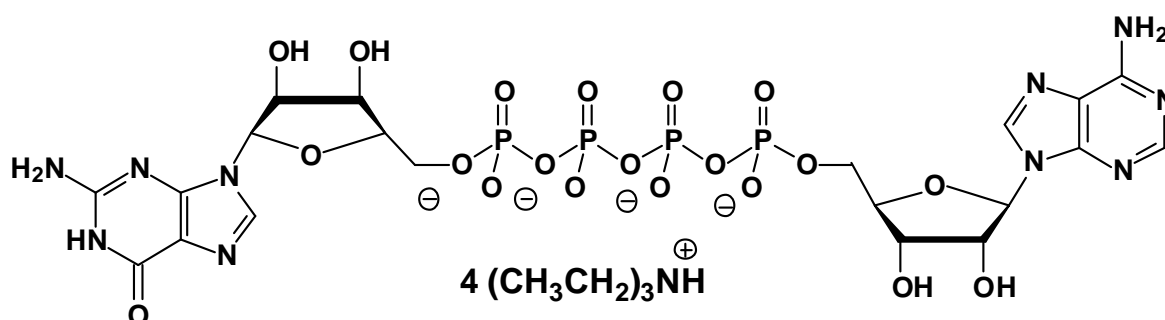


AP₄G

P¹-(5'-Adenosyl) P⁴-(5'-guanosyl) tetrphosphate, Triethylammonium salt

Cat. No.	Amount
NU-503S	50 Units
NU-503L	250 Units



Cat. No.: NU-503

Molecular Formula: C₂₀H₂₄N₁₀O₂₀P₄ (Anion)

Molecular Weight: 848.36 (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 μl of a 10 mM solution

Selected References:

- Safrany *et al.* (2007) Characterisation of a bis(5'-nucleosyl)-tetraphosphatase (asymmetrical) from *Drosophila melanogaster*. *Int. J. Biochem. Cell Biol.* **39** (5):943.
- Ortiz *et al.* (1993) specific synthesis of adenosine (5')tetraphospho(5')nucleoside and adenosine- (5')oligophospho(5')adenosine (n-greaterthan- 4) catalyzed by firefly luciferase. *Eur. J. Biochem.* **212** (1):263.
- Palfi *et al.* (1991) Alterations in the accumulation of adenylylated nucleotides in heavy-metal-ion-stressed and heat-stressed *Synechococcus* sp strain pcc-6301, a cyanobacterium, in light and dark. *Biochem. J.* **276**:487.
- Brevet *et al.* (1991) Isolation and characterization of a dinucleoside triphosphatase from *Saccharomyces-cerevisiae*. *J. Bacteriol.* **173** (17):5275.
- Sillero *et al.* (1991) Synthesis of dinucleoside polyphosphates catalyzed by firefly luciferase. *Eur. J. Biochem.* **202** (2):507.
- Avila *et al.* (1991) A paradoxical increase of a metabolite upon increased expression of its catabolic enzyme - the case of diadenosine tetraphosphate (AP4A) and AP4A phosphorylase-I in *Saccharomyces-cerevisiae*. *J. Bacteriol.* **173** (24):7875.
- Garrison *et al.* (1984) Assay of adenosine 5'-P1-tetraphospho-P4-5'''-adenosine and adenosine 5'-P1-tetraphospho-P4-5'''-guanosine in *Physarum-polycephalum* and other eukaryotes - an isocratic high-pressure liquid-chromatography method. *Biochem. J.* **217** (3):805.
- Barnes *et al.* (1984) Measurement of AP4A and AP4G during the cell-cycle of *Physarumpolycephalum*. *H-S Z. Physiol. Chem.* **365** (6):608.