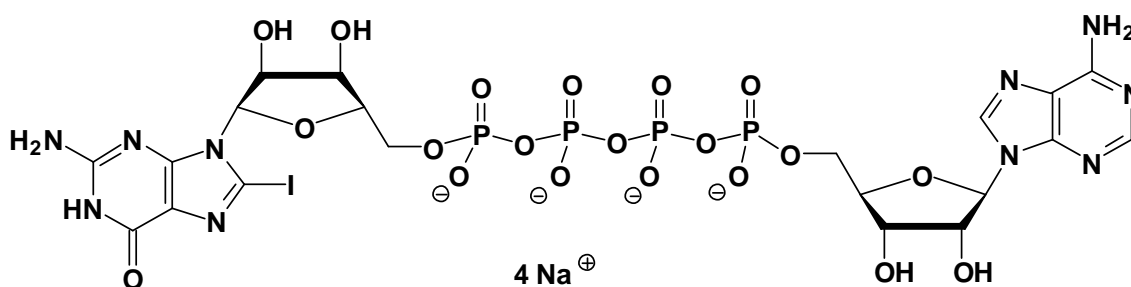


AP₄(8I-G)

P¹-(5'-Adenosyl) P⁴-(5'-(8-iodo)-guanosyl) tetraphosphate, Sodium salt

Cat. No.	Amount
NU-510S	30 Units
NU-510L	150 Units



Cat. No.: NU-510

Molecular Formula: C₂₀H₂₃IN₁₀O₂₀P₄ (Anion)

Molecular Weight: 973.91 (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