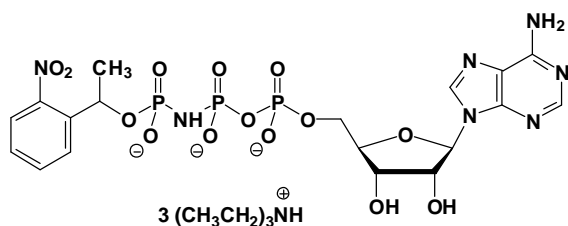


NPE-caged-AppNHp (NPE-caged-AMPPNP)

Adenosine-5'-[(β,γ)-imido]triphosphate, P³-(1-(2-nitrophenyl)-ethyl)-ester,
Triethylammonium salt

Cat. No.	Amount
NU-305S	20 Units
NU-305L	100 Units



Cat. No.: NU-305

Molecular Formula: C₁₈H₂₁N₇O₁₄P₃ (Anion)

Molecular Weight: 652.32 (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. Store in a dark place and protect from light!

For research use only!

1 unit = 1 μ l of a 10 mM solution