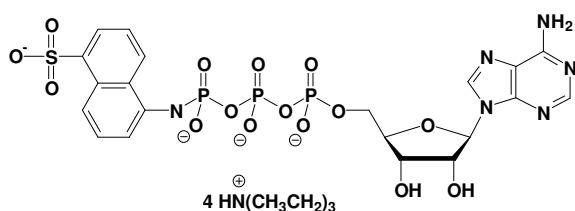


ATP- γ -AmNS

Adenosine-5'-triphosphate- γ -(sulfo-1-naphthyl)amide, Triethylammonium salt

Cat. No.	Amount
NU-1616S	150 μ l / 5 mM
NU-1616L	750 μ l / 5 mM



Cat. No.: NU-1616

Molecular Formula: C₂₀H₁₈N₆O₁₅P₃S (Anion)

Molecular Weight: 707.37 (Anion)

Purity: > 95%

Spectroscopic properties:

λ_{max} 323 nm; ϵ 4200;

λ_{exc} 323nm; λ_{em} 461 nm

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!