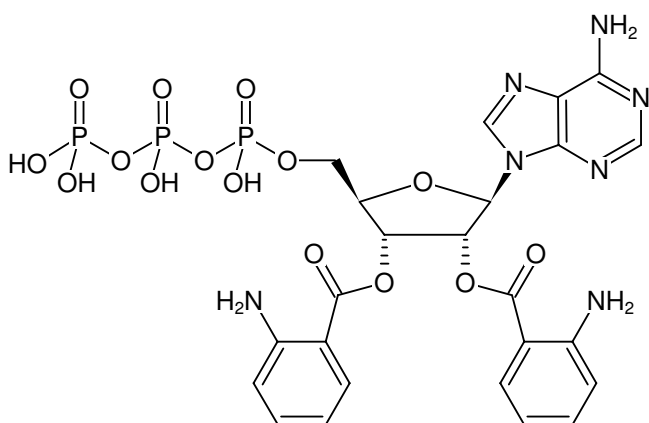


**Bis-Ant-ATP**

2'/3'-O-(Bis-anthraniloyl)-adenosine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-892S	150 µl (10 mM)
NU-892L	5 x 150 µl (10 mM)



Structural formula of Bis-Ant-ATP

For general laboratory use.**Shipping:** shipped on gel packs**Storage Conditions:** store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery**Molecular Formula:** C₂₄H₂₆N₇O₁₅P₃ (free acid)**Molecular Weight:** 745.42 g/mol (free acid)**Exact Mass:** 745.07 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ_{max} 250 nm, ε 20.2 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5), λ_{exc} 337 nm, λ_{em} 423 nm