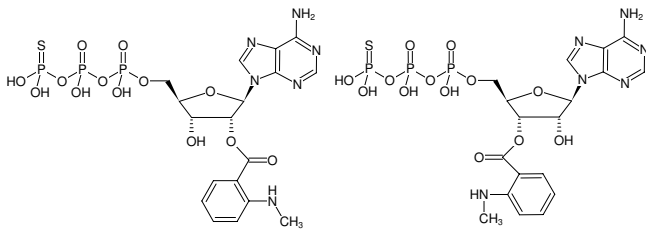




Mant-ATPyS

2'/3'-O-(N-Methyl-anthraniloyl)-adenosine-5'-(γ -thio)-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-232S	20 μ l (10 mM)
NU-232L	5 x 20 μ l (10 mM)



Structural formula of Mant-ATPyS

For research use only!

Shipping: shipped on blue ice

Storage Conditions: store at -20 °C

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

Shelf Life: 6 months after date of delivery

Molecular Formula: C₁₈H₂₃N₆O₁₃P₃S (free acid)

Molecular Weight: 656.39 g/mol (free acid)

Exact Mass: 656.03 g/mol (free acid)

Purity: \geq 90 % (HPLC)

Form: colorless to slightly yellow solution in water

Concentration: 10 mM - 11 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{max} 255/355 nm, ϵ 23.3/5.8 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5), λ_{exc} 355 nm, λ_{em} 448 nm

Applications:

Inhibition of heart adenylyl cyclase AC5^[1]

Inhibition of adenylyl cyclase from *Bordetella pertussis*^[2]

FRET^[2]

Selected References:

[1] Goettle *et al.* (2009) Characterization of Mouse Heart Adenylyl Cyclase. *J. Pharmacol. Exp. Ther.* **329** (3):1156.

[2] Goettle *et al.* (2007) Molecular analysis of the interaction of *Bordetella pertussis* adenylyl cyclase with fluorescent nucleotides. *Molecular Pharmacology* **72** (3):526.

Erdorf *et al.* (2011) Pharmacological characterization of adenylyl cyclase isoforms in rabbit kidney membranes. *Naunyn Schmiedebergs Arch. Pharmacol.* **383** (4):357.