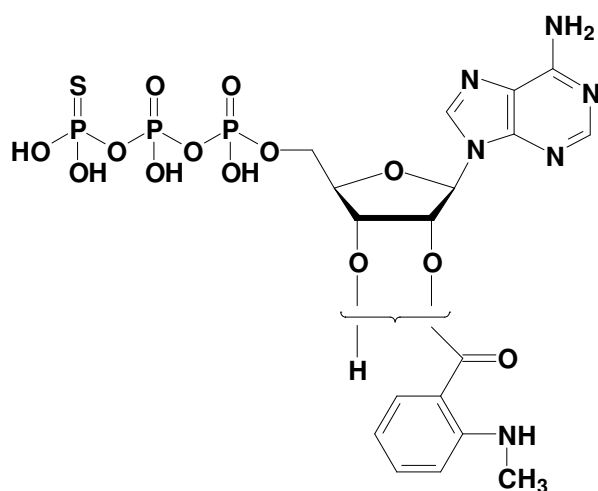


Mant-ATP- γ S

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

Cat. No.	Amount
NU-232S	20 Units
NU-232L	100 Units



Inhibition of heart adenylyl cyclase AC5^[1]

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

FRET^[2]

Selected References:

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

[2] Göttele *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 Schmiedeberg's Arch. Pharmacol.* **383** (4):357.

Cat. No.: NU-232

Molecular Formula: C₁₈H₂₃N₆O₁₃P₃S (free acid)

Molecular Weight: 656.39 (free acid)

Purity: > 90%, clear aqueous solution, pH 7.5

Spectroscopic properties:

λ_{\max} 255/355 nm; ϵ 23300/5800; λ_{exc} 355nm;
 λ_{em} 448 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 6 months after date of delivery.

For research use only!

* 1 unit = 1 μ l of a 10 mM solution

Applications: