

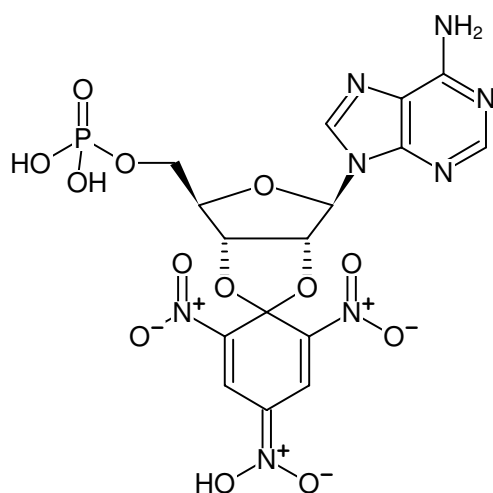
**TNP-AMP**

2',3'-O-Trinitrophenyl-adenosine-5'-monophosphate, Triethylammonium salt

Cat. No.	Amount
NU-224S	2 μ mol
NU-224L	5 x 2 μ mol

Selected References:

Sprang *et al.* (2006) Broad Specificity of Mammalian Adenylyl Cyclase for Interaction with 2',3'-Substituted Purine- and Pyrimidine Nucleotide Inhibitors. *Mol. Pharmacol.* **70**:878.



Structural formula of TNP-AMP

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:** C₁₆H₁₅N₈O₁₃P**Molecular Weight:** 558.31 g/mol**Exact Mass:** 558.05 g/mol**Purity:** \geq 95 % (HPLC)**Form:** solid**Color:** orange**Spectroscopic Properties:** λ_{exc} 408/470 nm, λ_{em} 552 nm, $\epsilon_{408\text{ nm}}$ 24.4 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)