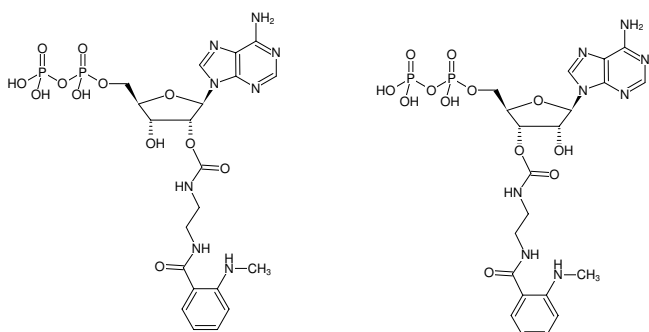


**EDA-ADP-MANT**

2'/3'-O-(2-Aminoethyl-carbamoyl)-adenosine-5'-diphosphate, labeled with MANT, Triethylammonium salt

Cat. No.	Amount
NU-802-MNT	200 µl (5 mM)



Structural formula of EDA-ADP-MANT

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: C₂₁H₂₈N₈O₁₂P₂ (free acid)

Molecular Weight: 646.45 g/mol (free acid)

Exact Mass: 646.13 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 5.0 mM - 5.5 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{exc} 335 nm, λ_{em} 440 nm, ε 2.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)