

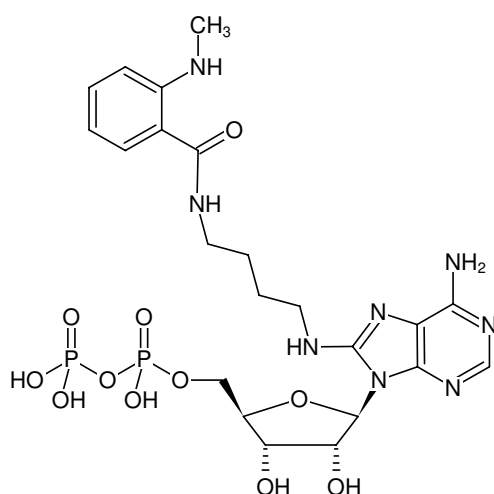


8-[(4-Amino)butyl]-amino-ADP-MANT

(MABA-ADP)

8-[(4-Amino)butyl]-amino-adenosine-5'-diphosphate, labeled with MANT, Triethylammonium salt

Cat. No.	Amount
NU-893-MNT	200 µl (5 mM)



Structural formula of 8-[(4-Amino)butyl]-amino-ADP-MANT

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: 12 months after date of delivery

Molecular Formula: C₂₂H₃₂N₈O₁₁P₂ (free acid)

Molecular Weight: 646.48 g/mol (free acid)

Exact Mass: 646.17 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, ε₂₈₀ 18.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Selected References:

Yan *et al.* (2019) MANF antagonizes nucleotide exchange by the endoplasmic reticulum chaperone BiP. *Nat. Commun.* **10** (1):541.

Leskovar *et al.* (2008) Photophysical properties of popular fluorescent adenosine nucleotide analogs used in enzyme mechanism probing. *Arch. Biochem. Biophys.* **473** (1):16.

Jakob *et al.* (1996) Assessment of the ATP binding properties of Hsp90. *J. Mol. Chem.* **271** (17):10035.

Theysen *et al.* (1996) The second step of ATP binding to DnaK induces peptide release. *J. Mol. Biol.* **263** (5):657.