

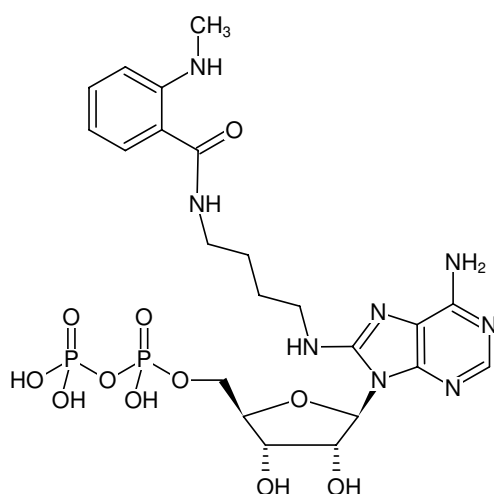


## 8-[(4-Amino)butyl]-amino-ADP-MANT

(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 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:** C<sub>22</sub>H<sub>32</sub>N<sub>8</sub>O<sub>11</sub>P<sub>2</sub> (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<sub>exc</sub> 335 nm, λ<sub>em</sub> 440 nm, ε<sub>280</sub> 18.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (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.