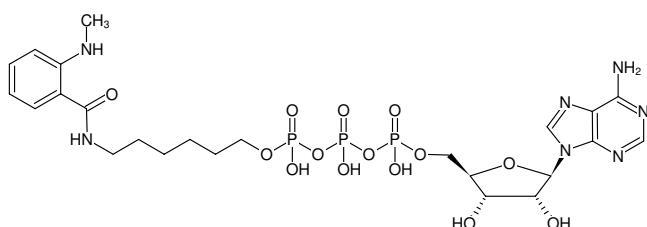




### $\gamma$ -(6-Aminoethyl)-ATP-MANT

$\gamma$ -(6-Aminoethyl)-adenosine-5'-triphosphate, labeled with MANT, Triethylammonium salt

Cat. No.	Amount
NU-833-MNT	200 $\mu$ l (5 mM)



Structural formula of  $\gamma$ -(6-Aminoethyl)-ATP-MANT

**For research use only!**

**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>24</sub>H<sub>36</sub>N<sub>7</sub>O<sub>14</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 739.51 g/mol (free acid)

**Exact Mass:** 739.15 g/mol (free acid)

**Purity:**  $\geq$  95 % (HPLC)

**Form:** solution in water

**Color:** colorless to slightly yellow

**Concentration:** 5.0 mM - 5.5 mM

**pH:** 7.5  $\pm$  0.5

**Spectroscopic Properties:**  $\lambda_{exc}$  335 nm,  $\lambda_{em}$  440 nm,  $\epsilon$  2.7 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)