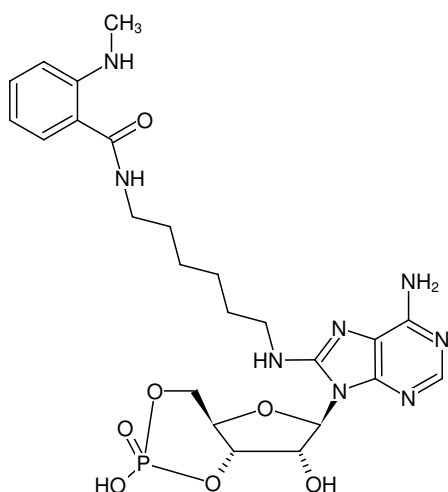




### 8-(6-Aminoethyl)-amino-cAMP-MANT

8-(6-Aminoethyl)-amino-adenosine-3',5'-cyclic monophosphate, labeled with MANT, Triethylammonium salt

| Cat. No.   | Amount        |
|------------|---------------|
| NU-851-MNT | 200 µl (1 mM) |



Structural formula of 8-(6-Aminoethyl)-amino-cAMP-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:** C<sub>24</sub>H<sub>33</sub>N<sub>8</sub>O<sub>7</sub>P (free acid)

**Molecular Weight:** 576.55 g/mol (free acid)

**Exact Mass:** 576.22 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

**Color:** colorless to slightly yellow

**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 335 nm, λ<sub>em</sub> 440 nm, ε 2.7 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)