

**cyclic-tri-AMP**

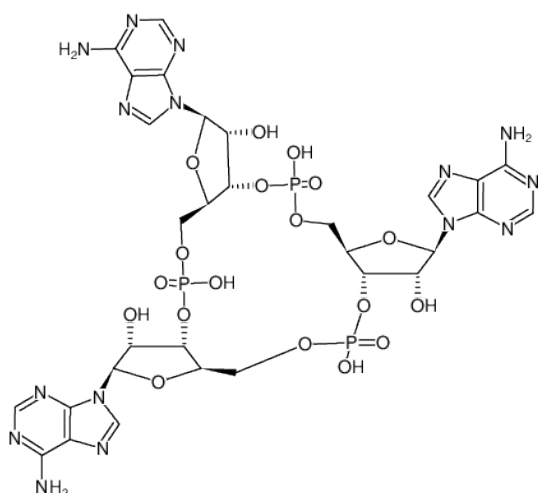
3',3',3'-cyclic AMP-AMP-AMP, c-tri-AMP

3',3',3'-Cyclic adenosine monophosphate-adenosine monophosphate-adenosine monophosphate, Sodium salt

Cat. No.	Amount
NU-458S	1 $\mu$ mol
NU-458L	5 x 1 $\mu$ mol

**Selected References:**

Wenzl *et al.* (2024) How enzyme-centered approaches are advancing research on cyclic oligo-nucleotides. *FEBS Lett.* **598**(8): 839.



Structural formula of cyclic-tri-AMP

**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>30</sub>H<sub>36</sub>N<sub>15</sub>O<sub>18</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 987.63 g/mol (free acid)**Exact Mass:** 987.16 g/mol (free acid)**CAS#:** 54447-85-7**Purity:**  $\geq$  98 % (HPLC) /  $\leq$  2.5 EU/ml Endotoxin (LAL test)**Form:** solid (lyophilized from filtered (10 kDa) solution in water)**Color:** white to off-white**Spectroscopic Properties:**  $\lambda_{\max}$  259 nm,  $\epsilon$  40.5 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)