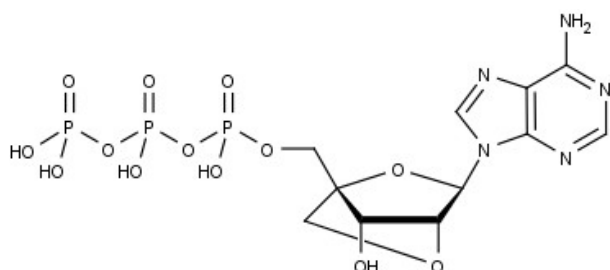


**LNA-ATP**

LNA-adenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-982	10 µl (100 mM)



Structural formula of LNA-ATP

**For *in vitro* 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>11</sub>H<sub>16</sub>N<sub>5</sub>O<sub>13</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 519.19 g/mol (free acid)**Exact Mass:** 519.00 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution**Concentration:** 100 mM - 110 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ<sub>max</sub> 259 nm, ε 15.1 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)**Selected References:**Crouzier *et al.* (2012) Efficient reverse transcription using locked nucleic acid nucleotides towards the evolution of nuclease resistant RNA aptamers. *PLoS One*. **7** (4):e35990.Doessin *et al.* (2012) Amplification and re-generation of LNA-modified libraries. *Molecules*. **17** (11):13087.Veedu *et al.* (2008) Polymerase chain reaction and transcription using locked nucleic acid nucleotide triphosphates. *J. Am. Chem. Soc.* **130** (26):8126.Veedu *et al.* (2007) Enzymatic incorporation of LNA nucleotides into DNA strands. *Chembiochem*. **8** (5):490.