

**ApG**

Sodium salt

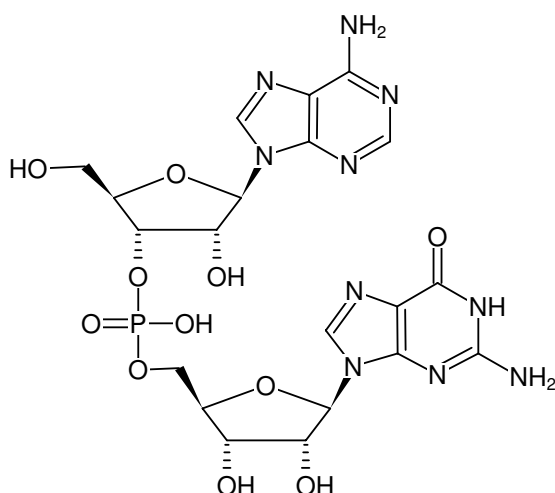
Cat. No.	Amount
NU-888S	10 µl (100 mM)
NU-888L	5 x 10 µl (100 mM)

**Applications:**

Initiator dinucleotide for the 5'-labeling of RNA by *in vitro* transcription<sup>[1]</sup>

**Selected References:**

[1] Pitulle *et al.* (1992) Initiator oligonucleotides for the combination of chemical and enzymatic RNA synthesis. *Gene*. **112 (1)**:101.



Structural formula of ApG

**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>20</sub>H<sub>25</sub>N<sub>10</sub>O<sub>11</sub>P (free acid)**Molecular Weight:** 612.45 g/mol (free acid)**Exact Mass:** 612.14 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** sterile solution in water**Color:** colorless to slightly yellow**Concentration:** 100 mM - 110 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ<sub>max</sub> 260 nm, ε 25.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)