

**5'-Biotin-ApG**

Sodium salt

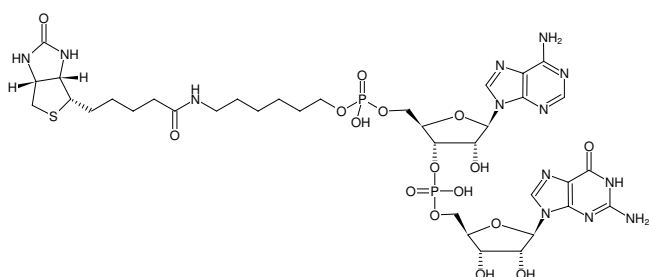
Cat. No.	Amount
NU-888-BIO-S	10 µl (100 mM)
NU-888-BIO-L	5 x 10 µl (100 mM)

Applications:

Initiator dinucleotide for the 5'-labeling of RNA by *in vitro* transcription^[1]

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 5'-Biotin-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:** C₃₆H₅₃N₁₃O₁₆P₂S (free acid)**Molecular Weight:** 1017.90 g/mol (free acid)**Exact Mass:** 1017.29 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_{max} 260 nm, ε 25.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)