

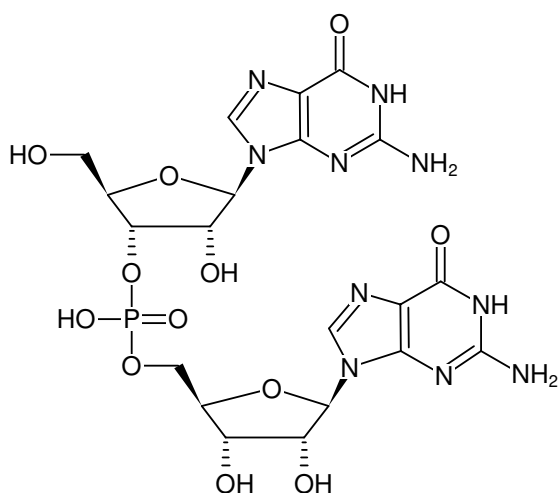
**GpG**

RNA Dinucleotide (5'→3'), Sodium salt

Cat. No.	Amount
NU-950S	20 µl (50 mM)
NU-950L	5 x 20 µl (50 mM)

Selected References:

Salvio *et al.* (2015) Ribonuclease Activity of an Artificial Catalyst That Combines a Ligated Cu(II) Ion and a Guanidinium Group at the Upper Rim of a cone-Calix[4]arene Platform. *J. Org. Chem.* **80** (11):5887.



Structural formula of GpG

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:** C₂₀H₂₅N₁₀O₁₂P (free acid)**Molecular Weight:** 628.45 g/mol (free acid)**Exact Mass:** 628.14 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 50 mM - 55 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ_{max} 260 nm, ε 21.6 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)