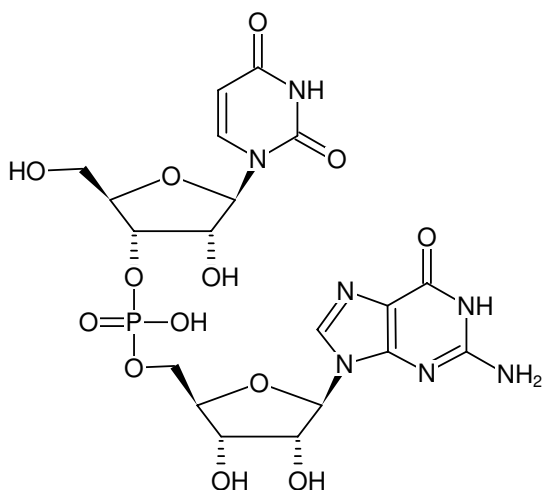


**UpG**

RNA Dinucleotide (5'→3'), Sodium salt

Cat. No.	Amount
NU-898S	10 µl (100 mM)
NU-898L	5 x 10 µl (100 mM)



Structural formula of UpG

**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>19</sub>H<sub>24</sub>N<sub>7</sub>O<sub>13</sub>P (free acid)**Molecular Weight:** 589.41 g/mol (free acid)**Exact Mass:** 589.12 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** 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 20.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)