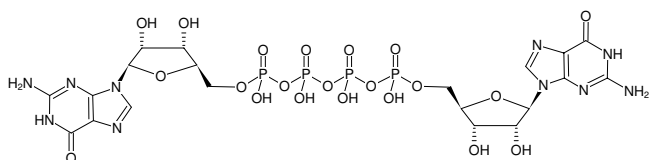


**GP<sub>4</sub>G**P<sup>1</sup>-(5'-Guanosyl)-P<sup>4</sup>-(5'-guanosyl)-tetraphosphate, Sodium salt

Cat. No.	Amount
NU-533S	20 µl (10 mM)
NU-533L	5 x 20 µl (10 mM)

Structural formula of GP<sub>4</sub>G**For general laboratory use.****Shipping:** shipped on gel packs**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>28</sub>N<sub>10</sub>O<sub>21</sub>P<sub>4</sub> (free acid)**Molecular Weight:** 868.39 g/mol (free acid)**Exact Mass:** 868.04 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ<sub>max</sub> 252 nm, ε 25.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)