

## 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 Units
NU-533L	100 Units



**Cat. No.:** NU-533

**Molecular Formula:** C<sub>22</sub>H<sub>30</sub>N<sub>10</sub>O<sub>21</sub>P<sub>4</sub> (Anion)

**Molecular Weight:** 894.42 (Anion)

**Purity:** > 95%, clear aqueous solution, pH 7.5

**Storage conditions:**

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

\* 1 unit = 1 µl of a 10 mM solution

**For research use only!**