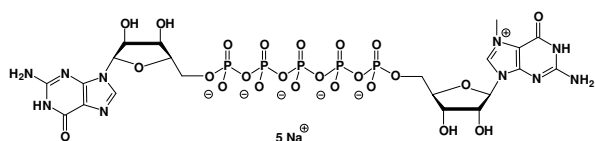


## m<sup>7</sup>GP<sub>5</sub>G

P<sup>1</sup>-(5'-7-Methyl-Guanosyl)-P<sup>5</sup>-(5'-Guanosyl)-pentaphosphate, Sodium salt

Cat. No.	Amount
NU-532S	20 Units
NU-532L	100 Units



**Cat. No.:** NU-532

**Molecular Formula:** C<sub>23</sub>H<sub>33</sub>N<sub>10</sub>O<sub>24</sub>P<sub>5</sub> (Anion)

**Molecular Weight:** 988.43 (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!**