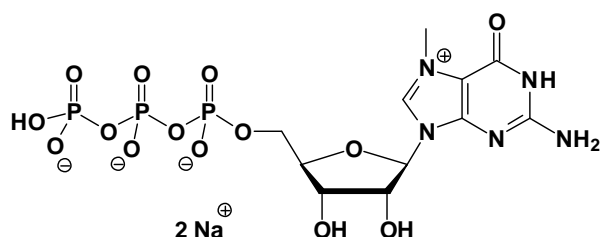


## m<sup>7</sup>GTP

7-methyl-guanosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1122S	100 Units
NU-1122L	500 Units



**Cat. No.:** NU-1122

**Molecular Formula:** C<sub>11</sub>H<sub>16</sub>N<sub>5</sub>O<sub>14</sub>P<sub>3</sub> (Anion)

**Molecular Weight:** 535.19 (Anion)

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

**Spectroscopic properties:** λ<sub>max</sub> 258/280 nm;  
ε 9800/8000

### 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.

### For research use only!

1 unit = 1 μl of a 10 mM solution

### Selected References:

Shen *et al.* (2001) Structural and thermodynamic behavior of eukaryotic initiation factor 4E in supramolecular formation with 4E-binding protein 1 and mRNA cap analogue, studied by spectroscopic methods. *Chem Pharm Bull* **49(10)**:1299.

Carberry *et al.* (1989) A spectroscopic study of the binding of m<sup>7</sup>GTP and m<sup>7</sup>GpppG to human protein synthesis initiation factor 4E. *Biochemistry*. **28(20)**:8078.

Beemon *et al.* (1977) In vitro translation yields a possible Rous sarcoma virus src gene product. *Proc Natl Acad Sci U S A*. **74(8)**:3302.