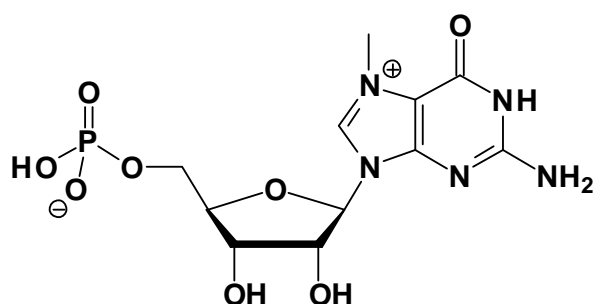


m⁷GMP

7-methyl-guanosine-5'-monophosphate

Cat. No.	Amount
NU-1135S	100 Units
NU-1135L	500 Units



Carberry *et al.* (1989) A spectroscopic study of the binding of m⁷GTP and m⁷GpppG 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.

Cat. No.: NU-1135

Molecular Formula: C₁₁H₁₆N₅O₈P

Molecular Weight: 377.25

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

Spectroscopic properties: λ_{max} 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.