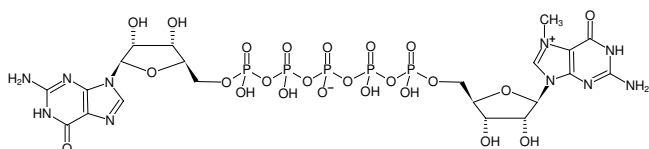


**m⁷GP₅G**P¹-(5'-7-Methyl-Guanosyl)-P⁵-(5'-Guanosyl)-pentaphosphate, Sodium salt

Cat. No.	Amount
NU-532S	20 µl (10 mM)
NU-532L	5 x 20 µl (10 mM)

Structural formula of m⁷GP₅G**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:** C₂₁H₃₁N₁₀O₂₄P₅ (free acid)**Molecular Weight:** 962.39 g/mol (free acid)**Exact Mass:** 962.02 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_{max} 260 nm, ε 21.6 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)