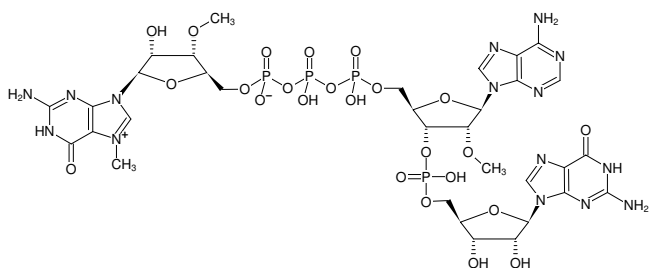



 $m_2^{7,3'-O}GP_3(2'OMe)ApG$ (Cap Analog)
 $m_7(3'OMeG)(5')ppp(5')(2'OMeA)pG$, Sodium salt

Cat. No.	Amount
NU-248S	10 μ l (100 mM)
NU-248L	5 x 10 μ l (100 mM)


 Structural formula of $m_2^{7,3'-O}GP_3(2'OMe)ApG$ (Cap Analog)

For research use only!
Shipping: shipped on blue ice

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: $C_{33}H_{45}N_{15}O_{24}P_4$ (free acid)

Molecular Weight: 1159.70 g/mol (free acid)

Exact Mass: 1159.17 g/mol (free acid)

Purity: $\geq 95\%$ (HPLC)

Form: clear aqueous solution

Concentration: 100 mM - 110 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{max} 255 nm, ϵ 30.5 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)