



$m_2^{7,3'-O}GP_3G$ (ARCA Cap Analog) - Solution

$m_2^{7,3'-O}G(5')ppp(5')G$

$P^1-(5'-(3'-O-methyl)-7-methyl-guanosyl) P^3-(5'-(guanosyl))triphosphate$, Sodium salt

Cat. No.	Amount
NU-855S	10 μ l (100 mM)
NU-855L	5 x 10 μ l (100 mM)

Related Products:

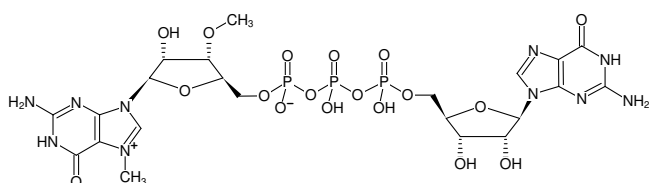
HighYield T7 RNA Synthesis Kit, #RNT-101

Selected References:

Jemielity *et al.* (2003) Novel anti-reverse cap analogs with superior translational properties. *RNA* **9**:1108.

Peng *et al.* (2002) Synthesis and application of a chain-terminating dinucleotide mRNA cap analog. *Org. Lett.* **4** (2):161.

Stepinski *et al.* (2001) Synthesis and properties of mRNAs containing the novel anti-reverse cap analogs 7-methyl (3'-O-methyl)GpppG and 7-methyl (3'-deoxy)GpppG. *RNA* **7**:1486.



Structural formula of $m_2^{7,3'-O}GP_3G$ (ARCA Cap Analog) - Solution

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at $-20\text{ }^\circ\text{C}$

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: $C_{22}H_{31}N_{10}O_{18}P_3$ (free acid)

Molecular Weight: 816.46 g/mol (free acid)

Exact Mass: 816.10 g/mol (free acid)

CAS#: 400806-46-4 (acid), 400806-60-2 (Trisodium salt)

Purity: $\geq 98\%$ (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 100 mM - 110 mM

pH: 7.25 ± 0.25

Spectroscopic Properties: λ_{max} 255 nm, ϵ 22.6 L $mmol^{-1} cm^{-1}$ (Tris-HCl pH 7.5)