

**Ant-m⁷GTP**

2'/3'-O-Anthraniloyl-7-methylguanosine-5'-triphosphate, Triethylammonium salt

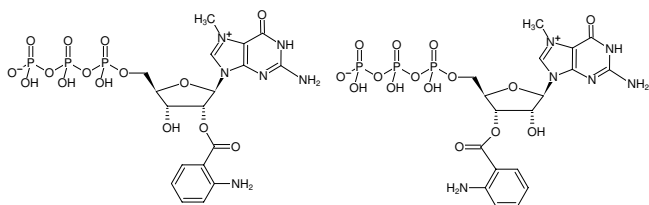
Cat. No.	Amount
NU-227S	100 µl (10 mM)
NU-227L	5 x 100 µl (10 mM)

Selected References:

Gunawardana *et al.* (2015) Efficient preparation and properties of mRNAs containing a fluorescent cap analog: Anthraniloyl-m⁷GpppG *Translation (Austin)*. **3 (1)**:e988538.

Wei *et al.* (1998) Wheat germ poly (A) binding protein enhances the binding affinity of eukaryotic initiation factor 4F and (iso)4F for cap analogues. *Biochemistry* **37 (7)**:1910.

Ren *et al.* (1996) Synthesis of a fluorescent 7-methylguanosine analog and a fluorescence spectroscopic study of its reaction with wheatgerm cap binding proteins. *Nucleic Acids Res.* **24 (18)**:3629.

Structural formula of Ant-m⁷GTP**For research use only!****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:** C₁₈H₂₄N₆O₁₅P₃ (free acid)**Molecular Weight:** 656.33 g/mol (free acid)**Exact Mass:** 656.04 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} 258/280/332 nm, ε 9.8/8.0/4.6 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5), λ_{exc} 332 nm, λ_{em} 423 nm