

**tC⁰TP**tC⁰ (1,3-diaza-2-oxophenoxazine) ribonucleoside triphosphate, Ammonium salt

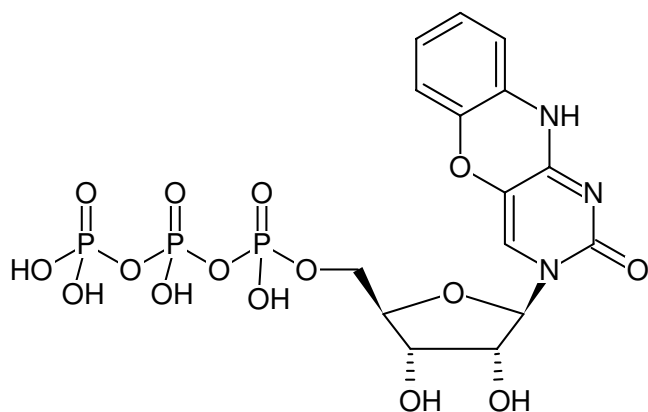
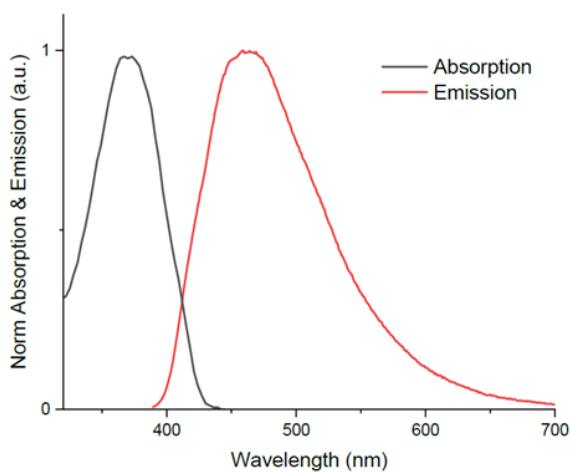
Cat. No.	Amount
NU-315	10 µl (10 mM)

For general laboratory use.**Please centrifuge briefly before opening (volume ≤2 ml).****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:** C₁₅H₁₈N₃O₁₅P₃ (free acid)**Molecular Weight:** 573.24 g/mol (free acid)**Exact Mass:** 573.00 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in 10 mM Tris-HCl**Color:** yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{exc} 360 nm, λ_{em} 460 nm, ε 10.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Applications:**

Incorporation into RNA/cRNA

- Random labeling via in vitro Transcription with T7 and SP6 RNA Polymerase^[1]- 3'-End Labeling of RNA oligonucleotides with Terminal deoxynucleotidyl Transferase (TdT)^[1]**Selected References:**[1] Baladi *et al.* (2021) Stealth Fluorescence Labeling for Live Microscopy Imaging of mRNA Delivery. *J Am Chem Soc.* **143(14)**:5413-5424.Structural formula of tC⁰TPexcitation and emission spectrum of tC⁰TP