

**5'-Cy3-ApU**

Cy3-labeled ApU
Triethylammonium salt

Cat. No.	Amount
NU-917-CY3	10 µl (5 mM)

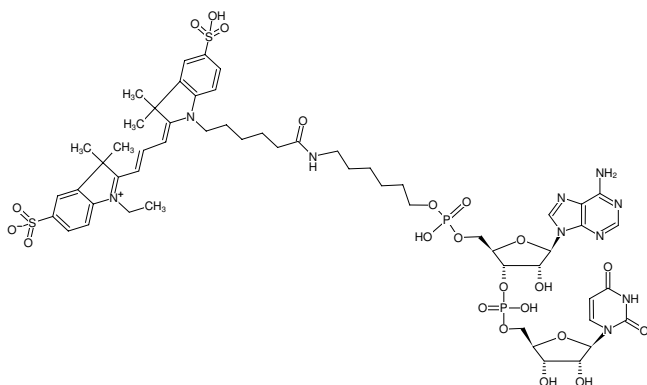
Applications:

Initiator dinucleotide for the 5'-labeling of RNA by *in vitro* transcription

Selected References:

Kang *et al.* (2020) Transcription reinitiation by recycling RNA polymerase that diffuses on DNA after releasing terminated RNA. *Nat Commun.* **11(1)**:450.

Kang *et al.* (2021) Hopping and Flipping of RNA Polymerase on DNA during Recycling for Reinitiation after Intrinsic Termination in Bacterial Transcription. *Int J Mol Sci.* **22(5)**:2398.



Structural formula of 5'-Cy3-ApU

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: C₅₆H₇₄N₁₀O₂₂P₂S₂ (free acid)

Molecular Weight: 1365.32 g/mol (free acid)

Exact Mass: 1364.39 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: filtered solution (30 kDa) in water

Color: pink

Concentration: 5 mM - 5.5 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{exc} 550 nm, λ_{em} 570 nm, ε 150.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)