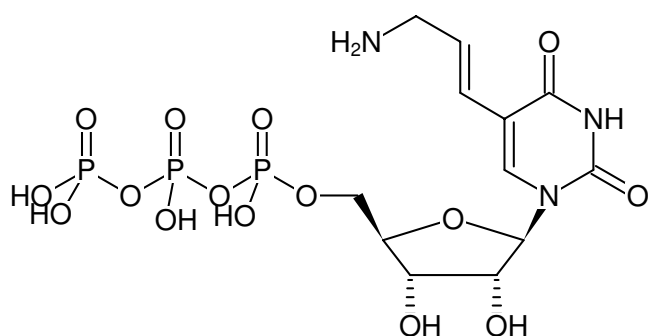




Aminoallyl-UTP - Solution

5-(3-Aminoallyl)-uridine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-821S	50 µl (50 mM)
NU-821L	5 x 50 µl (50 mM)



Structural formula of Aminoallyl-UTP - Solution

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: C₁₂H₂₀N₃O₁₅P₃ (free acid)

Molecular Weight: 539.22 g/mol (free acid)

Exact Mass: 539.01 g/mol (free acid)

CAS#: 112131-73-4

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 50 mM - 55 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{max} 249/289 nm, ε 10.7/7.1 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Applications:

Nucleic Acid Sequence Based Amplification^[1]

Microarray based hybridization detection^[1]

RNA-labeling^[2]

Selected References:

[1] Scheler *et al.* (2009) Fluorescent labeling of NASBA amplified tmRNA molecules for microarray applications. *BMC Biotechnol.* **9**:45.

[2] Hoen *et al.* (2003) Fluorescent labelling of cRNA for microarray applications. *Nucleic Acids Res.* **31** (5):e20.

Labesse *et al.* (2011) Structural and functional characterization of the Mycobacterium tuberculosis uridine monophosphate kinase: insights into the allosteric regulation. *Nucleic Acids Res.* **39** (8):3458.