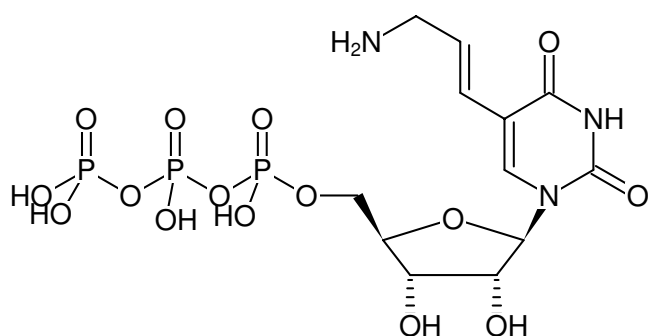




## Aminoallyl-UTP - Solid

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

Cat. No.	Amount
NU-821-1	1 mg
NU-821-5	5 mg



Structural formula of Aminoallyl-UTP - Solid

**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:** C<sub>12</sub>H<sub>17</sub>N<sub>3</sub>O<sub>15</sub>P<sub>3</sub> (Anion)

**Molecular Weight:** 539.22 g/mol (free acid)

**Exact Mass:** 539.01 g/mol (free acid)

**CAS#:** 112131-73-4

**Purity:** ≥ 95 % (HPLC)

**Form:** solid

**Color:** white to off-white

**Spectroscopic Properties:** λ<sub>max</sub> 249/289 nm, ε 10.7/7.1 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

### Applications:

Nucleic Acid Sequence Based Amplification<sup>[1]</sup>

Microarray based hybridization detection<sup>[1]</sup>

RNA-labeling<sup>[2]</sup>

### 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.

Gilles 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.