

# PRODUCT INFORMATION

## 5-(3-Aminoallyl)uridine-5'-O-triphosphate (sodium salt)

Item No. 38423

CAS Registry No.: 85280-66-6

Formal Name: 5-(3-amino-1-propenyl)-uridine  
5'-(tetrahydrogen triphosphate),  
tetrasodium salt

Synonyms: AA-UTP, aminoallyl-UTP

MF:  $C_{12}H_{16}N_3O_{15}P_3 \cdot 4Na$

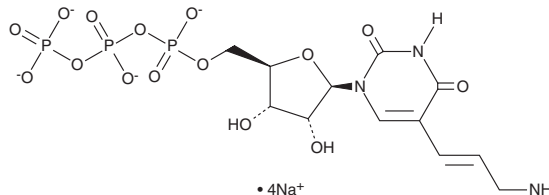
FW: 627.2

Purity:  $\geq 95\%$

Supplied as: A solution in water

Storage:  $-80^{\circ}C$

Stability:  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Description

5-(3-Aminoallyl)uridine-5'-O-triphosphate (aminoallyl-UTP) is an amine-modified nucleotide that has been used in the synthesis of affinity-tagged UTP.<sup>1</sup> Incorporation of aminoallyl-UTP into complementary RNA (cRNA) and bacterial transfer-messenger RNA (tmRNA) has been used in the generation of fluorescently labeled probes for microarrays.<sup>2,3</sup>

### References

1. Langer, P.R., Waldrop, A.A., and Ward, D.C. Enzymatic synthesis of biotin-labeled polynucleotides: Novel nucleic acid affinity probes. *Proc. Natl. Acad. Sci. USA* **78**, 6633-6637 (1981).
2. t Hoen, P.A., de Kort, F., ban Ommen, G.J.B., *et al.* Fluorescent labelling of cRNA for microarray applications. *Nucleic Acids Res.* **31(5)**, e20 (2003).
3. Scheler, O., Glynn, B., Parkel, S., *et al.* Fluorescent labeling of NASBA amplified tmRNA molecules for microarray applications. *BMC Biotechnol.* **9**, 45 (2009).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 04/14/2026

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM