

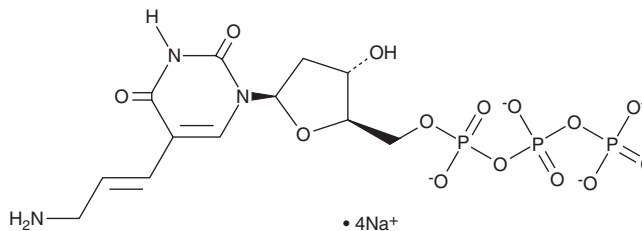
PRODUCT INFORMATION



5-(3-Aminoallyl)-2'-deoxyuridine-5'-O-triphosphate (sodium salt)

Item No. 38361

CAS Registry No.: 85280-65-5
Formal Name: 5-(3-amino-1-propen-1-yl)-2'-deoxy-uridine 5'-(tetrahydrogen triphosphate), tetrasodium salt
Synonyms: AA-dUTP, aminoallyl-dUTP
MF: C₁₂H₁₆N₃O₁₄P₃ • 4Na
FW: 611.2
Purity: ≥98%
Supplied as: A solution in water
Storage: -80°C
Stability: ≥2 years



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

Description

5-(3-Aminoallyl)-2'-deoxyuridine-5'-O-triphosphate (aminoallyl-dUTP) is an amine-modified nucleotide that has been used in the synthesis of fluorescently labeled dUTP.¹ Incorporation of aminoallyl-dUTP into primers has been used in the generation of fluorescently labeled probes for DNA microarrays and FRET-based assays to measure the activity of reverse transcriptases.^{2,3}

References

1. Kuznetsova, V.E., Spitsyn, M.A., Shershov, V.E., *et al.* Novel fluorescently labeled nucleotides: Synthesis, spectral properties and application in polymerase chain reaction. *Mendeleev Commun.* **26(2)**, 95-98 (2016).
2. Xiang, C.C., Kozhich, O.A., Chen, M., *et al.* Amine-modified random primers to label probes for DNA microarrays. *Nat. Biotechnol.* **20(7)**, 738-742 (2002).
3. Krebs, J.F. and Kore, A.R. Novel FRET-based assay to detect reverse transcriptase activity using modified dUTP analogues. *Bioconjug. Chem.* **19(1)**, 185-191 (2008).

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, 05/01/2023

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