

Allylamine-dUTP

Allylamine-2'-deoxyuridine-5'-triphosphate can replace TTP in reactions where it serves as a substrate for *E. coli* DNA polymerase 1 (holoenzyme and Klenow fragment), terminal deoxynucleotide transferase, T4 and Taq DNA polymerases, and reverse transcriptases (from AMV and M-MuLV). Allylamine-dUTP is incorporated into DNA by a variety of labeling reactions including nick translation, random primed DNA synthesis and 3' end labeling reactions. Allylamine-labeled DNA can be efficiently conjugated to the active ester form of signal generating moieties such as fluorescent dyes, biotin and other haptens to produce labeled probes for hybridization/detection assays. The efficient incorporation of Allylamine-dUTP, in contrast to many dye-labeled nucleotides, provides for greater incorporation of signaling moieties and stronger signals.

Citations: 1

[View Online »](#)

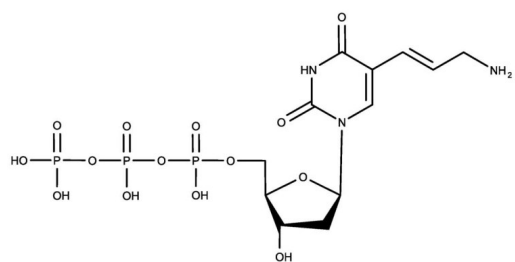
Ordering Information

[Order Online »](#)

ENZ-42861	2.5μmol
-----------	---------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Allylamino-2'-deoxyuridine-5'-triphosphate
Appearance	Clear, colorless solution, free of solid particulates.
Concentration	50mM
Extinction Coefficient	8,100 M ⁻¹ cm ⁻¹ (290 nm, water)
Formula	C ₁₂ H ₂₀ N ₃ O ₁₄ P ₃
Formulation	Liquid. Solution in water.
MW	523.2 (free acid)
Purity	≥93% (HPLC)
Purity Detail	Purified by ion-exchange chromatography.
Technical Info / Product Notes	Several of Enzo's products and product applications are covered by US and foreign patents and patents pending.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)