

Aminoallyl-dUTP – Solid (1 mg)

Ordering Information

[Order Online »](#)

| | |
|--------------|-----|
| JBS-NU-803-1 | 1mg |
|--------------|-----|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 5-(3-Aminoallyl)-2'-deoxyuridine-5'-triphosphate, Sodium salt

Appearance white to off-white solid

Application Notes Incorporation into DNA/cDNA by PCR with Taq polymerase
Nick Translation with DNase I/ DNA Polymerase I
Primer Extension with Klenow exo-
3'-End Labeling with Terminal deoxynucleotidyl Transferase (TdT)
Reverse Transcription with MMLV Reverse Transcriptase

CAS 116840-18-7

Couple Type Modified nucleotides

Formula $C_{12}H_{20}N_3O_{14}P_3$ (free acid)

MW 523.22 (free acid)

Purity ≥ 95 % (HPLC)

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



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)