

# Aminoallyl-dUTP-XX- ATTO-643 (L pack)

## Ordering Information

[Order Online »](#)

JBS-NU-803-XX-643-L	5x10µl
---------------------	--------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Appearance blue solution

Application Notes Incorporation into DNA/cDNA by  
– PCR with Taq polymerase  
– Nick Translation with DNase I/ DNA Polymerase I

Concentration 1.0 mM – 1.1 mM

Couple Type Modified nucleotides

Emission Maximum 665 nm

Excitation Maximum 643 nm

Formula  $C_{24}H_{41}N_5O_{16}P_3$  – ATTO 643 (free acid)

Formulation sterile solution in 10 mM Tris-HCl

MW 1567.90 (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)