

# Aminoallyl-dUTP-XX – ATTO-425, L pack

Aminoallyl-dUTP-XX-ATTO425 is recommended for direct enzymatic labeling of DNA/cDNA e.g. by PCR and Nick Translation. It is incorporated as substitute for its natural counterpart dTTP. The resulting Dye-labeled DNA/cDNA probes are ideally suited for fluorescence hybridization applications such as FISH or microarray-based gene expression profiling. Optimal substrate properties and thus labeling efficiency is ensured by an optimized linker attached to the C5 position of uridine. Recommended Aminoallyl-dUTP-XX-ATTO425/dTTP ratio for PCR and Nick Translation: 20-30% Aminoallyl-dUTP-XX-ATTO425/ 80-70% dTTP (PCR), 30-50% Aminoallyl-dUTP-XX-ATTO425/ 70-50% dTTP (Nick Translation) Please note: Protect the Dye-labeled dUTP from exposure to light and carry out experimental procedures in low light conditions. The optimal final concentration of the Dye-labeled dUTP may vary depending on the application and assay conditions. For optimal product yields and high incorporation rates an individual optimization of the Dye-labeled-dUTP/dTTP ratio is recommended.

## Ordering Information

[Order Online »](#)

JBS-NU-803-XX-425-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

**Alternative Name** 5-(3-Aminoallyl)-2'-deoxyuridine-5'-triphosphate, labeled with ATTO 425, Triethylammonium salt

**Appearance** green-yellow solution

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

**Concentration** 1 mM – 1.1 mM

**Couple Type** Modified nucleotides

**Emission Maximum** 485 nm

**Excitation Maximum** 439 nm

**Formula**  $C_{46}H_{67}N_6O_{21}P_3$  (free acid)

**Formulation** sterile solution in 10 mM Tris-HCl

**MW** 1132.97 (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-fr@enzolifesciences.com](mailto:info-fr@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)