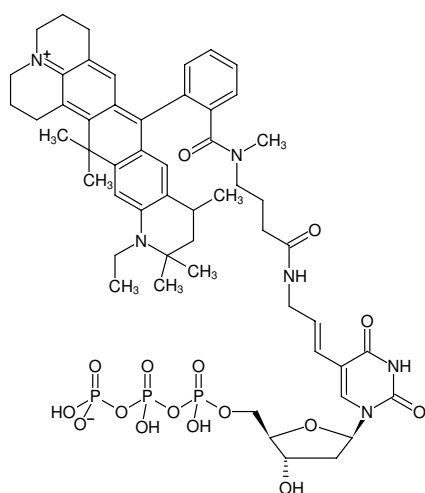




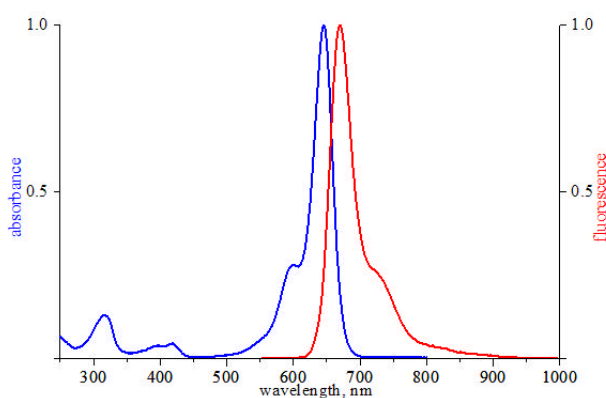
## Aminoallyl-dUTP-ATTO-647N

5-(3-Aminoallyl)-2'-deoxyuridine-5'-triphosphate, labeled with ATTO 647N, Triethylammonium salt

Cat. No.	Amount
NU-803-647N-S	10 $\mu$ l (1 mM)
NU-803-647N-L	5 x 10 $\mu$ l (1 mM)



Structural formula of Aminoallyl-dUTP-ATTO-647N



excitation and emission spectrum of ATTO 647N

**For general laboratory use.**

**Please centrifuge briefly before opening (volume  $\leq$  2 ml).**

**Shipping:** shipped on gel packs

**Storage Conditions:** store at  $-20\text{ }^{\circ}\text{C}$

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

**Shelf Life:** 12 months after date of delivery

**Molecular Formula:**  $\text{C}_{54}\text{H}_{69}\text{N}_6\text{O}_{16}\text{P}_3$  (free acid)

**Molecular Weight:** 1151.08 g/mol (free acid)

**Exact Mass:** 1150.40 g/mol (free acid)

**Purity:**  $\geq 95\%$  (HPLC)

**Form:** solution in 10 mM Tris-HCl

**Color:** blue

**Concentration:** 1.0 mM - 1.1 mM

**pH:**  $7.5 \pm 0.5$

**Spectroscopic Properties:**  $\lambda_{\text{exc}}$  646 nm,  $\lambda_{\text{em}}$  664 nm,  $\epsilon$  150.0 L  $\text{mmol}^{-1}\text{ cm}^{-1}$  (Tris-HCl pH 7.5)

### Applications:

Incorporation into DNA/cDNA by  
- Nick Translation with DNase I/ DNA Polymerase I <sup>in-house data</sup>

### Description:

Aminoallyl-dUTP-ATTO647N is recommended for direct enzymatic labeling of DNA/cDNA by 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-ATTO647N/dTTP ratio for Nick Translation: 30-50% Aminoallyl-dUTP-ATTO647N/ 50% dTTP

*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.*