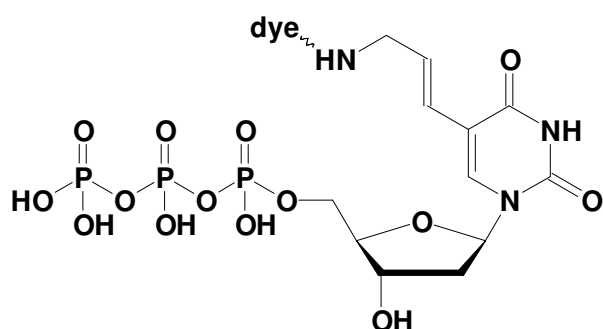


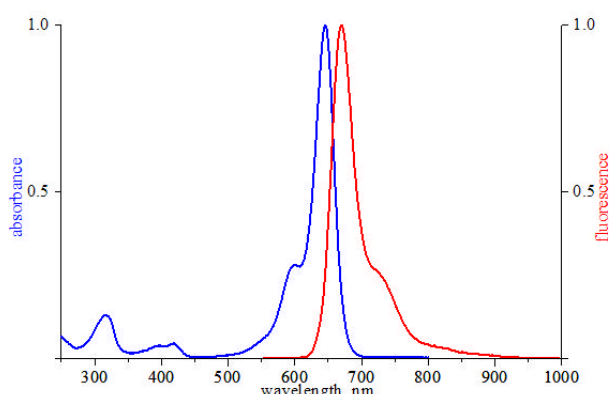
Aminoallyl-dUTP - ATTO 647N

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

Cat. No.	Amount
NU-803-647N	10 μ l / 1 mM



structural formula of Aminoallyl-dUTP



excitation and emission spectrum of ATTO 647N

Cat. No.: NU-803-647N

Molecular Formula: C₁₂H₂₀N₃O₁₄P₃ - ATTO 647N
 (free acid)

Molecular Weight: 1151.22 (free acid)

Purity: > 95%, clear aqueous solution, pH 7.5

Spectroscopic properties:

λ_{exc} 644 nm; λ_{em} 669 nm; ϵ 150,000 cm⁻¹ M⁻¹

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

For research use only!

Applications:

FISH^[1-4]

Microarray based hybridization detection^[3, 5]

Differential display^[6]

DNA-labeling^[7, 8]

Nick translation^[1, 9]

Selected References:

[1] Kocanova *et al.* (2010) Activation of estrogen-responsive genes does not require their nuclear co-localization. *PLoS Genet.* **6** (4):e1000922.

[2] Cox *et al.* (2004) Fluorescent DNA hybridization probe preparation using amine modification and reactive dye coupling. *BioTechniques* **36**:114.

[3] Williams *et al.* (2006) Genomic DNA as a general cohybridization standard for ratiometric microarrays. *Methods in Enzymology* **410**:237.

[4] Gao *et al.* (2007) Identification of the differential expressive tumor associated genes in rectal cancers by cDNA microarray. *Gastroenterology* **13**:341.

Aminoallyl-dUTP - ATTO 647N

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

[5] Wen *et al.* (2004) A visual DNA chip for simultaneous detection of hepatitis B virus, hepatitis C virus and human immunodeficiency virus type-1. *Biosen. Bioelek.* **19**:685.

[6] Cheviron *et al.* (2008) Transcriptomic variation and plasticity in rufous-collared sparrows (*Zonotrichia capensis*) along an altitudinal gradient. *Mol. Ecology* **17**:4556.

[7] Dirsch *et al.* (2007) Probe production for in situ hybridization by PCR and subsequent covalent labelling with fluorescent dyes. *Appl. Immunochem.* **15**:332.

[8] Cherkasov *et al.* (2010) New Nucleotide Analogues with Enhanced Signal Properties. *Bioconjugate Chem.* **21**:122.

[9] Nimmakayalu *et al.* (2000) Simple method for preparation of fluor/hapten-labeled dUTP. *Biotechniques* **28**:518.

Mayr JA *et al.* (2011) Heterozygous mutation in the X chromosomal NDUFA1 gene in a girl with complex I deficiency. *Mol Genet Metab.* **103 (4)**:358.