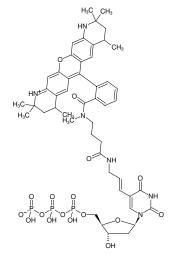




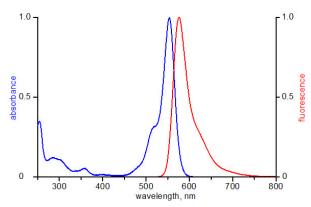
## Aminoallyl-dUTP-ATTO-550

5-(3-Aminoallyl)-2'-deoxyuridine-5'-triphosphate, labeled with ATTO 550, Triethylammonium salt

Cat. No.	Amount
NU-803-550-S	10 μl (1 mM)
NU-803-550-L	5 x 10 μl (1 mM)



Structural formula of Aminoallyl-dUTP-ATTO-550



excitation and emission spectrum of ATTO 550

## For general laboratory use.

**Shipping:** shipped on gel packs **Storage Conditions:** store at -20 °C

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

Shelf Life: 12 months after date of delivery
Molecular Formula: C<sub>49</sub>H<sub>61</sub>N<sub>6</sub>O<sub>17</sub>P<sub>3</sub> (free acid)
Molecular Weight: 1098.97 g/mol (free acid)
Exact Mass: 1098.33 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

Form: filtered solution (30 kDa) in 10 mM Tris-HCl

Color: red-violet

Concentration: 1.0 mM - 1.1 mM

**pH:** 7.5 ±0.5

Spectroscopic Properties:  $\lambda_{exc}$  554 nm,  $\lambda_{em}$  576 nm,

 $\epsilon$  120.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

## **Applications:**

Incorporation into DNA/cDNA by

- Nick Translation with DNAse I DNA Polymerase I in-house data

## **Description:**

Aminoallyl-dUTP-ATTO550 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-ATTO550/dTTP ratio Nick Translation: 30-50% Aminoallyl-dUTP-ATTO550/ 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 very depending on the application and assay conditions. For optimal produdct yields and high incorporation rates an individual optimization of the Dye-labeled-dUTP/dTTP ratio is recommended.