

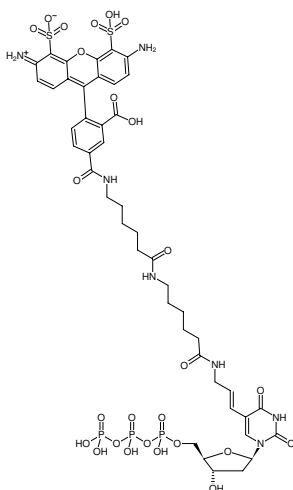


## Aminoallyl-dUTP-XX-AF488

also known as Alexa Fluor 488®-dUTP

5-(3-Aminoallyl)-2'-deoxyuridine-5'-triphosphate, labeled with AF488, Triethylammonium salt

Cat. No.	Amount
NU-803-XX-AF488-S	10 µl (1 mM)
NU-803-XX-AF488-L	5 x 10 µl (1 mM)



Structural formula of Aminoallyl-dUTP-XX-AF488

### 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>45</sub>H<sub>54</sub>N<sub>7</sub>O<sub>26</sub>P<sub>3</sub>S<sub>2</sub> (free acid)

**Molecular Weight:** 1265.99 g/mol (free acid)

**Exact Mass:** 1265.18 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

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

**Color:** yellow

**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 494 nm, λ<sub>em</sub> 515 nm, ε 73.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

### Applications:

Incorporation into DNA/cDNA by

- PCR with *Taq* polymerase <sup>in-house data</sup>

- Nick Translation with DNase I/ DNA Polymerase I <sup>in-house data</sup>

### Description:

Aminoallyl-dUTP-XX-AF488 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. AF488 (also known as Alexa Fluor 488®) is a hydrophilic dye with excellent photostability compared to fluorescein.

Recommended Aminoallyl-dUTP-XX-AF488/dTTP ratio for PCR and Nick Translation: 30-50% Aminoallyl-dUTP-XX-AF488/ 70-50% 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.*

### Related Products:

HighFidelity AF488 PCR Labeling Kit, #APP-101-AF488

HighFidelity GREEN PCR Labeling Testkit, #APP-101-GREEN

AF488 NT Labeling Kit, #PP-305-AF488