

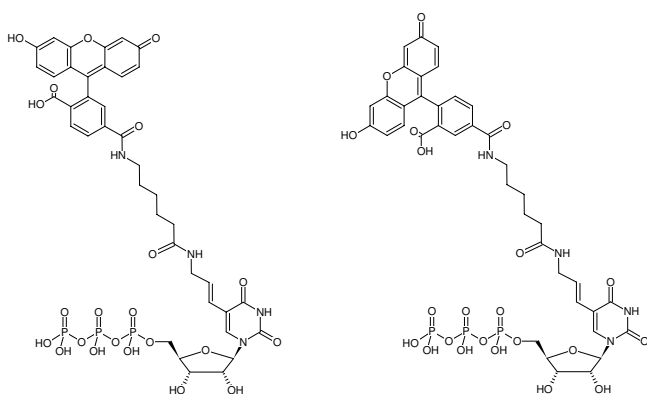


## Fluorescein-12-UTP

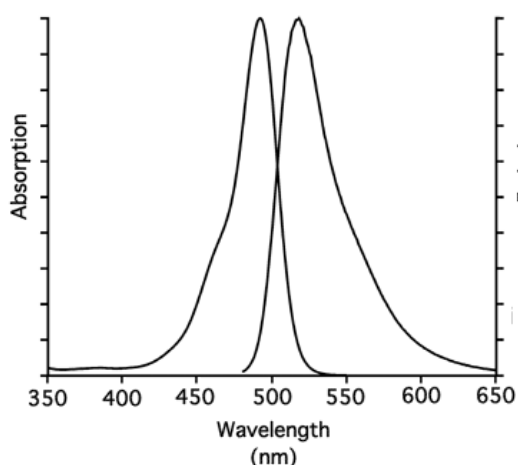
Fluorescein-X-(5-aminoallyl)-UTP

5/6-Fluorescein-X-(5-aminoallyl)-uridine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-821-FAMX-S	10 µl (5 mM)
NU-821-FAMX-L	5 x 10 µl (5 mM)



Structural formula of Fluorescein-12-UTP



excitation and emission spectrum of 5/6-FAM

### For research use only!

**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>39</sub>H<sub>41</sub>N<sub>4</sub>O<sub>22</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 1010.68 g/mol (free acid)

**Exact Mass:** 1010.14 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

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

**Color:** yellow-orange

**Concentration:** 5.0 mM - 5.5 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 492 nm, λ<sub>em</sub> 517 nm, ε 83.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

### Applications:

Genome wide expression profiling<sup>[1]</sup>

Identification of light and temperature entrained oscillating transcripts<sup>[1]</sup>

Analysis of gene regulation<sup>[2]</sup>

Estimation of evolutionary homologies<sup>[2]</sup>

### Related Products:

HighYield T7 RNA Synthesis Kit, #RNT-101

HighYield T7 Fluorescein RNA Labeling Kit, #RNT-101-FAMX

### Selected References:

[1] Collins *et al.* (2010) Genome-wide analyses reveal a role for peptide hormones in planarian germline development. *PLoS Biol.* **8** (10):e1000509.

[2] Dunn *et al.* (2007) Molecular paleoecology: using gene regulatory analysis to address the origins of complex life cycles in the late Precambrian. *Evol Dev.* **9** (1):10.