

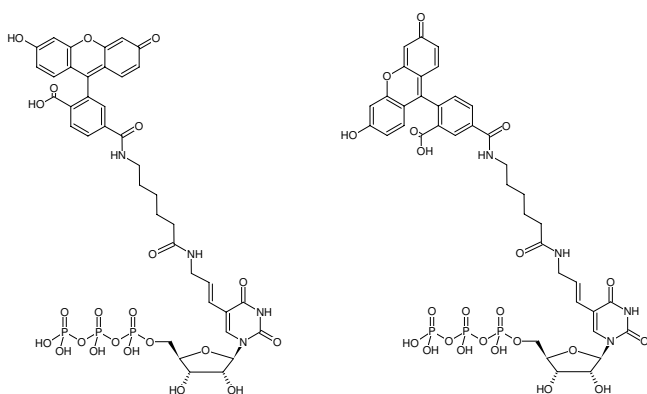


Fluorescein-12-UTP

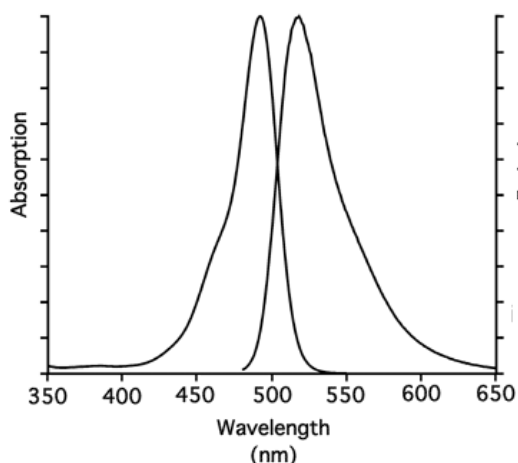
Fluorescein-X-(5-aminoallyl)-UTP

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 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: C₃₉H₄₁N₄O₂₂P₃ (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_{exc} 492 nm, λ_{em} 517 nm, ε 83.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Applications:

Genome wide expression profiling^[1]

Identification of light and temperature entrained oscillating transcripts^[1]

Analysis of gene regulation^[2]

Estimation of evolutionary homologies^[2]

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.