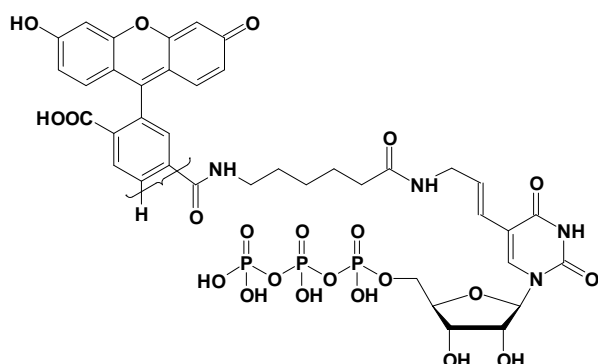


Fluorescein-12-UTP

5/6-Fluorescein-X-5-Aminoallyl-uridine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-821-FAMX	30 μ l / 1 mM



Cat. No.: NU-821-FAMX

Molecular Formula: C₃₉H₄₁N₄O₂₂P₃ (free acid)

Molecular Weight: 1010.68 (free acid)

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

Spectroscopic properties:

λ_{exc} 495 nm; λ_{em} 520 nm; ϵ 70,000 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:

Genome wide expression profiling^[1]

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

Analysis of gene regulation^[2]

Estimation of evolutionary homologies^[2]

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.