



Fluorescein-12-dATP

Fluorescein-X-7-Deaza-7-propargylamino-dATP

5/6-Fluorescein-X-7-deaza-7-propargylamino-2'-deoxyadenosine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-1611-FAMX	100 µl (1 mM)

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: C₄₁H₄₁N₆O₁₉P₃ (free acid)

Molecular Weight: 1014.72 g/mol (free acid)

Exact Mass: 1014.16 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: yellow-orange

Concentration: 1.0 mM - 1.1 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:

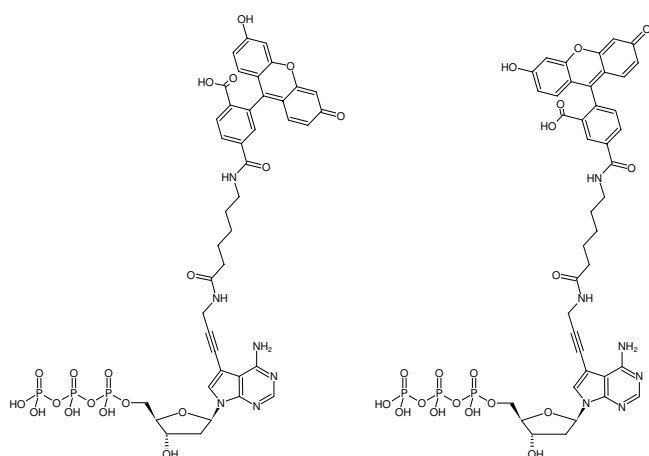
Incorporation by primer extension^[2]

Incorporation by PCR^[1]

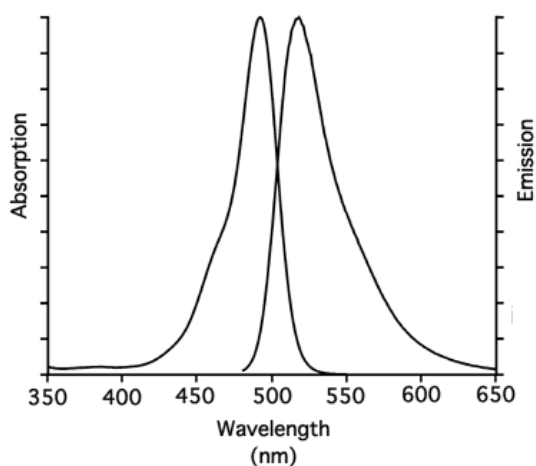
Selected References:

[1] Vrána *et al.* (2000) Flow sorting of mitotic chromosomes in common wheat (*Triticum aestivum* L.). *Genetics* **156** (4):2033.

[2] Bulyk *et al.* (1999) Quantifying DNA-protein interactions by double-stranded DNA arrays. *Nat. Biotechnol.* **17** (6):573.



Structural formula of Fluorescein-12-dATP



excitation and emission spectrum of 5/6-FAM