



Fluorescein-12-ddATP

Fluorescein-X-7-Deaza-7-propargylamino-ddATP

5/6-Fluorescein-X-7-deaza-7-propargylamino-2',3'-dideoxyadenosine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-1612-FAMX	30 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: C₄₁H₄₁N₆O₁₈P₃ (free acid)

Molecular Weight: 998.72 g/mol (free acid)

Exact Mass: 998.17 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

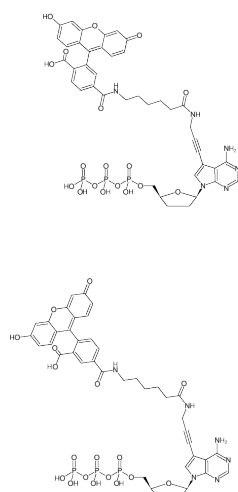
Form: solution in water

Color: yellow-orange

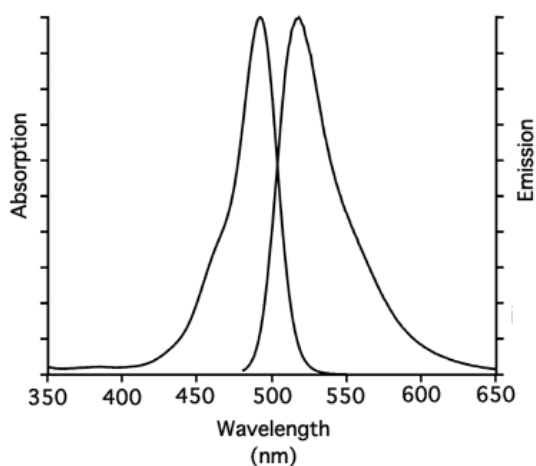
Concentration: 1.0 mM - 1.1 mM

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)



Structural formula of Fluorescein-12-ddATP



excitation and emission spectrum of 5/6-FAM