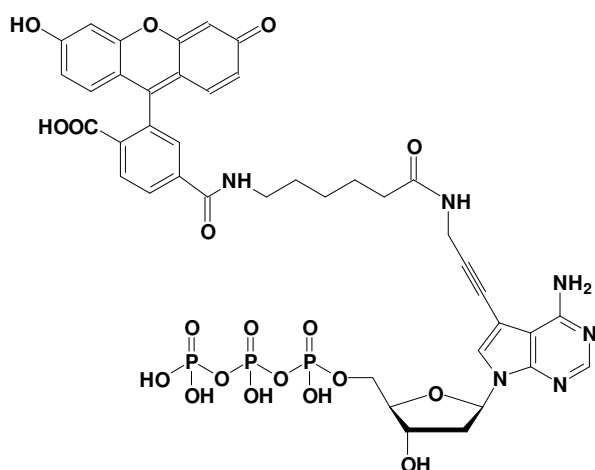


Fluorescein-12-dATP

Fluorescein-X-7-Deaza-7-propargylamino-2'-deoxy-adenosine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-1611-FAMX	25 μ l / 1 mM



structural formula of Fluorescein-12-dATP

Cat. No.: NU-1611-FAMX

Molecular Formula: C₄₁H₄₁N₆O₁₉P₃ (free acid)

Molecular Weight: 1014.72 (free acid)

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

Spectroscopic properties:

λ_{exc} 492 nm; λ_{em} 517 nm; ϵ 83,000 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!