

7-Propargylamino-7-deaza-dATP - DY 751

7-Deaza-7-propargylamino-2'-deoxy-adenosine-5'-triphosphate, labeled with DY 751,
 Triethylammonium salt

Cat. No.	Amount
NU-1611-751	40 μ l (1 mM)

Cat. No.: NU-1611-751

Molecular Formula: C₅₇H₇₁N₆O₂₀P₃S₂ (free acid)

Molecular Weight: 1317.26 (free acid)

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

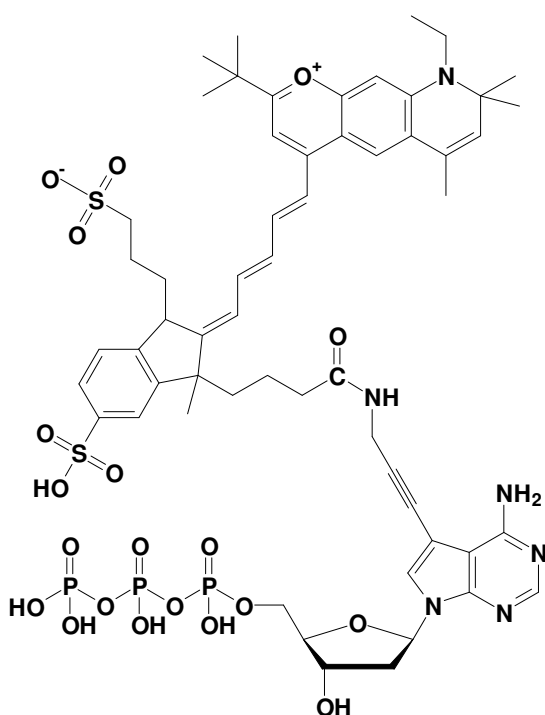
Spectroscopic properties:

λ_{exc} 751 nm; λ_{em} 779 nm; ϵ 220,000 cm⁻¹ M⁻¹

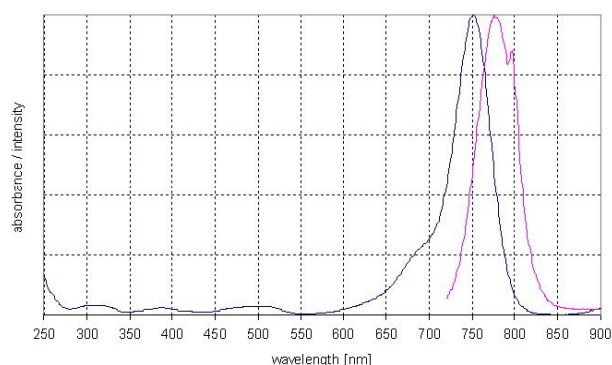
Storage conditions:

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!



structural formula of 7-Propargylamino-7-deaza-dATP - DY 751



excitation and emission spectrum of DY 751