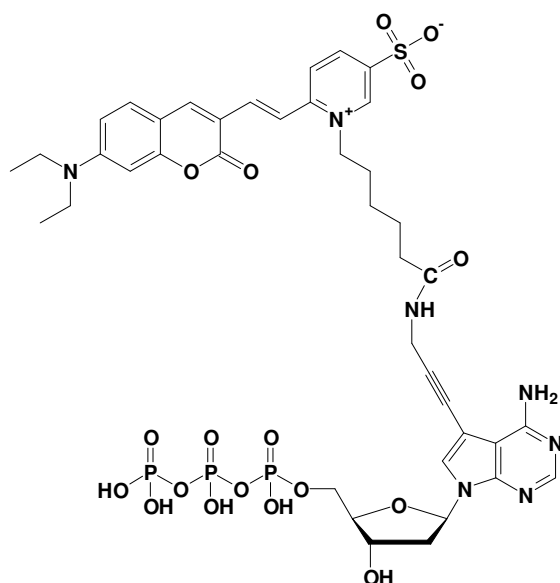


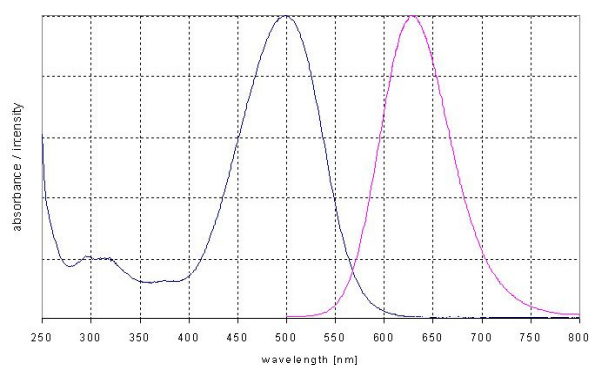
7-Propargylamino-7-deaza-dATP - DY 480XL

7-Deaza-7-propargylamino-2'-deoxy-adenosine-5'-triphosphate, labeled with DY 480XL, Triethylammonium salt

Cat. No.	Amount
NU-1611-480XL	10 μ l / (1 mM)



structural formula of 7-Propargylamino-7-deaza-dATP - DY 480XL



excitation and emission spectrum of DY 480XL

Cat. No.: NU-1611-480XL

Molecular Formula: C₄₀H₄₈N₇O₁₈P₃S (free acid)

Molecular Weight: 1039,83 (free acid)

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

Spectroscopic properties:

λ_{exc} 500 nm; λ_{em} 630 nm; ϵ 50,000 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!