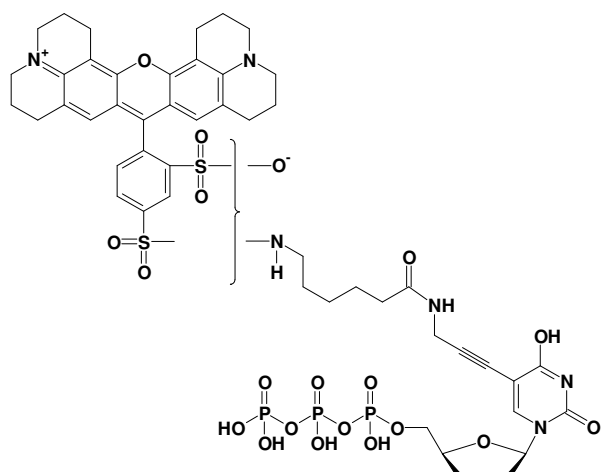


5-Propargylamino-ddUTP - Texas Red

5-Propargylamino- 2', 3'-Dideoxy-uridine-5'-triphosphate, labeled with Texas Red, Triethylammonium salt

Cat. No.	Amount
NU-1619-TXR	240 μ l (0,5 mM)



structural formula of 5-Propargylamino-ddUTP - Texas Red

Cat. No.: NU-1619-TXR

Molecular Formula: C₄₉H₅₇N₆O₂₀P₃S₂ (free acid)

Molecular Weight: 1207,06 (free acid)

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

Spectroscopic properties:

λ_{exc} 588 nm; λ_{em} 609 nm; ϵ 80.000 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!