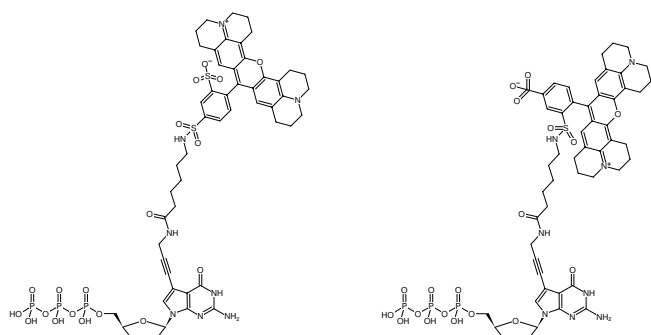




7-Propargylamino-7-deaza-ddGTP-Texas Red

7-Deaza-7-propargylamino-2',3'-dideoxyguanosine-5'-triphosphate, labeled with Texas Red, Triethylammonium salt

Cat. No.	Amount
NU-1618-TXR	240 µl (0,5 mM)



Structural formula of 7-Propargylamino-7-deaza-ddGTP-Texas Red

For research use only!

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: C₅₁H₅₉N₈O₁₉P₃S₂ (free acid)

Molecular Weight: 1245.11 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: red-violet

Concentration: 0.50 mM - 0.55 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{exc} 588 nm, λ_{em} 609 nm, ε 80.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)