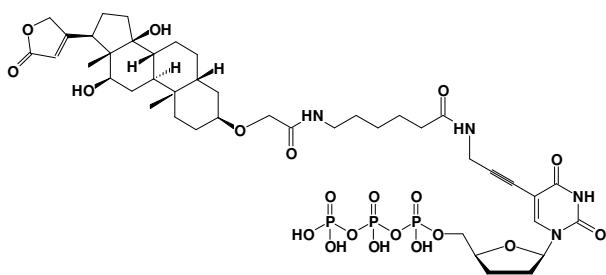


Digoxigenin-11-ddUTP

Digoxigenin-3-O-methylcarbonyl- ϵ -aminocaproyl-[5-(3-aminopropargyl)-2',3'-dideoxy-uridine-5'-triphosphate], Triethylammoniumsalt

Cat. No.	Amount
NU-1619-DIGX	25 μ l / 1mM



Cat. No.: NU-1619-DIGX

Molecular Formula: C₄₃H₆₃N₄O₂₀P₃ (free acid)

Molecular Weight: 1048.90 (free acid)

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

Spectroscopic properties: λ_{\max} 290 nm;
 ϵ 13000 M⁻¹cm⁻¹ Tris HCl pH 7.5

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!