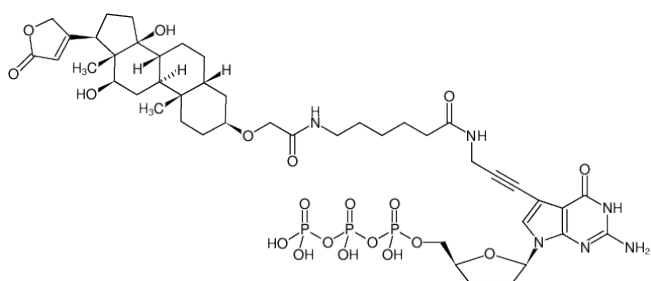


**DIG-11-ddGTP**

Digoxigenin-  
3-O-methylcarbonyl- $\epsilon$ -aminocaproyl-[7-deaza-7-propargylamino-2',3'-dideoxyguanosine-5'-triphosphate],  
Triethylammonium salt

Cat. No.	Amount
NU-1618-DIGX	25 $\mu$ l (1 mM)



Structural formula of DIG-11-ddGTP

**For general laboratory use.**

**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:** C<sub>45</sub>H<sub>65</sub>N<sub>6</sub>O<sub>19</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 1086.96 g/mol (free acid)

**Exact Mass:** 1086.35 g/mol (free acid)

**Purity:**  $\geq$  95 % (HPLC)

**Form:** solution in water

**Color:** colorless to slightly yellow

**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5  $\pm$  0.5

**Spectroscopic Properties:**  $\lambda_{\text{max}}$  272 nm,  $\epsilon$  11.9 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)