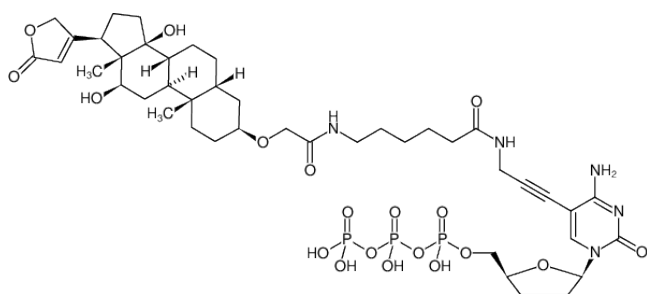


**DIG-11-ddCTP**

Digoxigenin-3-O-methylcarbonyl- $\epsilon$ -aminocaproyl-[5-(3-aminopropargyl)-2',3'-dideoxycytidine-5'-triphosphate], Triethylammoniumsalt

Cat. No.	Amount
NU-850-DIGX	25 $\mu$ l (1 mM)



Structural formula of DIG-11-ddCTP

**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>43</sub>H<sub>64</sub>N<sub>5</sub>O<sub>19</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 1047.92 g/mol (free acid)

**Exact Mass:** 1047.34 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:**  $\lambda_{\max}$  294 nm,  $\epsilon$  9.3 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)