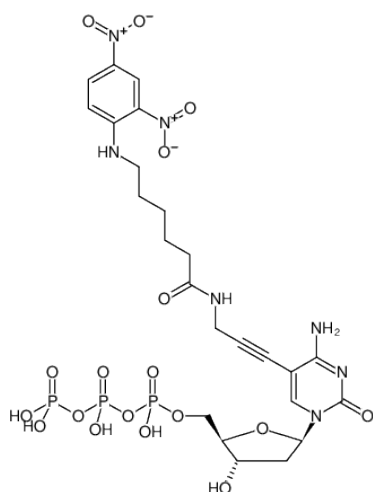


**5-Propargylamino-dCTP-DNP**

DNP-11-dCTP

5-Propargylamino-2'-deoxycytidine-5'-triphosphate, labeled with Dinitrophenol, Triethylammonium salt

Cat. No.	Amount
NU-809-DNP	25 µl (1 mM)



Structural formula of 5-Propargylamino-dCTP-DNP

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:** C₂₄H₃₂N₇O₁₈P₃**Molecular Weight:** 799.47 g/mol**Exact Mass:** 799.10 g/mol**Purity:** ≥ 95 % (HPLC)**Form:** filtered solution (30 kDa) in 10 mM Tris-HCl**Color:** yellow**Concentration:** 1.0 mM - 1.1 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 361 nm, ε 17.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)