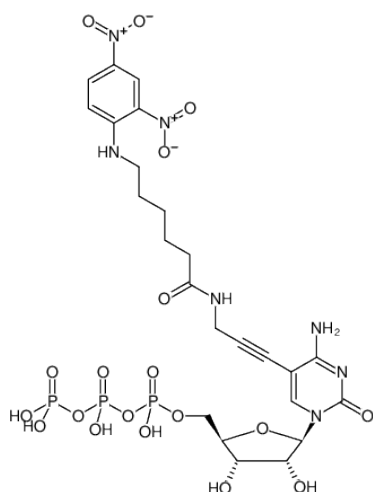


**5-Propargylamino-CTP-DNP**

DNP-11-CTP

5-Propargylamino-cytidine-5'-triphosphate, labeled with Dinitrophenol, Triethylammonium salt

Cat. No.	Amount
NU-831-DNP	30 µl (10 mM)



Structural formula of 5-Propargylamino-CTP-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:** C<sub>24</sub>H<sub>32</sub>N<sub>7</sub>O<sub>19</sub>P<sub>3</sub>**Molecular Weight:** 815.47 g/mol**Exact Mass:** 815.10 g/mol**Purity:** ≥ 95 % (HPLC)**Form:** filtered solution (30 kDa) in 10 mM Tris-HCl**Color:** yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ<sub>max</sub> 361 nm, ε 17.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)