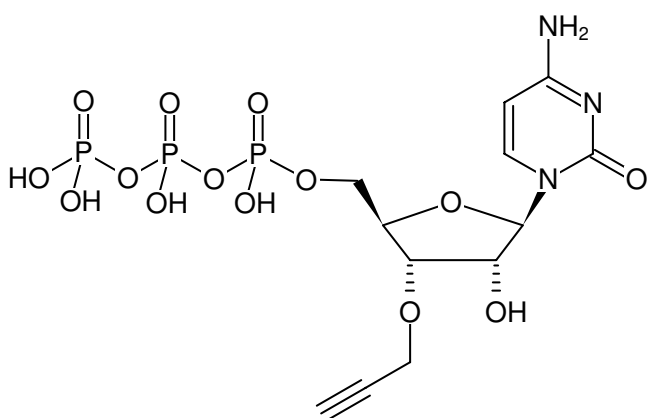




### 3'-(O-Propargyl)-CTP

3'-(O-Propargyl)-cytosine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-947	1 $\mu$ mol



Structural formula of 3'-(O-Propargyl)-CTP

#### For general laboratory use.

**Shipping:** shipped on gel packs

**Storage Conditions:** store at -20 °C

**Additional Storage Conditions:** Short term exposure (up to 1 week cumulative) to ambient temperature possible.

**Shelf Life:** 12 months after date of delivery

**Molecular Formula:** C<sub>12</sub>H<sub>18</sub>N<sub>3</sub>O<sub>14</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 521.20 g/mol

**Exact Mass:** 521.00 g/mol

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

**Form:** clear sticky semisolid

**Color:** white to off-white

**Spectroscopic Properties:**  $\lambda_{\text{max}}$  271 nm,  $\epsilon$  8.9 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)