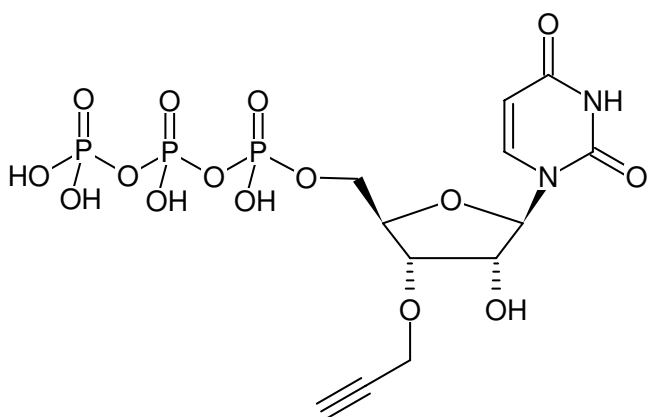


**3'-(O-Propargyl)-UTP**

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

Cat. No.	Amount
NU-948	1 $\mu\text{mol}$



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

**For research use only!****Shipping:** shipped on blue ice**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:**  $\text{C}_{12}\text{H}_{17}\text{N}_2\text{O}_{15}\text{P}_3$  (free acid)**Molecular Weight:** 522.19 g/mol**Exact Mass:** 521.98 g/mol**Purity:**  $\geq 95\%$  (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:**  $\lambda_{\text{max}}$  262 nm,  $\epsilon$  9.8 L  $\text{mmol}^{-1}$   $\text{cm}^{-1}$  (Tris-HCl pH 7.5)