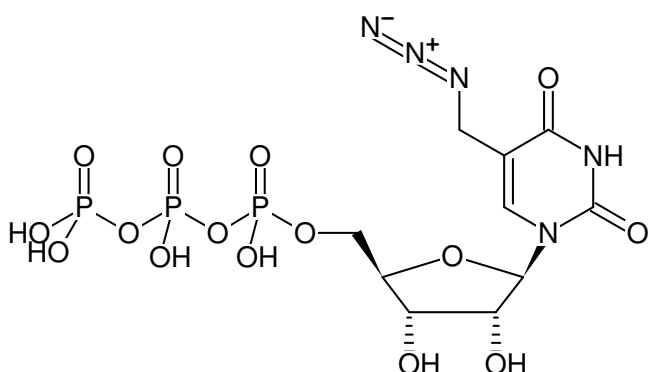


**5-Azidomethyl-UTP**

5-AmUTP

5-Azidomethyl-uridine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
CLK-083S	20 µl (10 mM)
CLK-083L	5 x 20 µl (10 mM)



Structural formula of 5-Azidomethyl-UTP

**For research use only!****Shipping:** shipped on blue ice**Storage Conditions:** store at -20 °C**Additional Storage Conditions:** avoid freeze/thaw cycles

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

**Shelf Life:** 12 months after date of delivery**Molecular Formula:** C<sub>10</sub>H<sub>16</sub>N<sub>5</sub>O<sub>15</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 539.18 g/mol (free acid)**Purity:** ≥ 90 % (HPLC)**Form:** sterile solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ<sub>max</sub> 260 nm, ε 8.8 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)