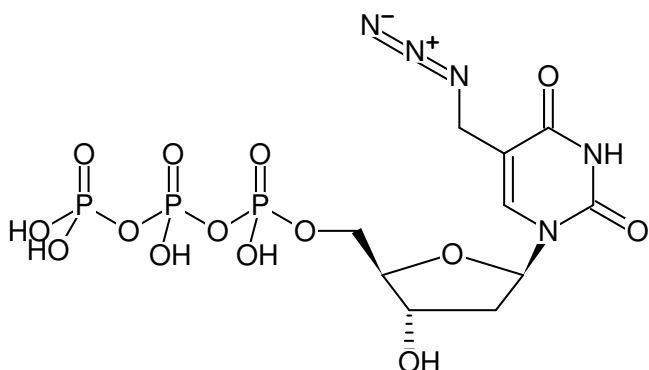


**5-Azidomethyl-dUTP**

5-AmdUTP

5-Azidomethyl-2'-deoxyuridine-5'-triphosphate, Sodium salt

Cat. No.	Amount
CLK-084S	20 µl (10 mM)
CLK-084L	5 x 20 µl (10 mM)



Structural formula of 5-Azidomethyl-dUTP

**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>10</sub>H<sub>16</sub>N<sub>5</sub>O<sub>14</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 523.18 g/mol (free acid)**Exact Mass:** 522.99 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ<sub>max</sub> 267 nm, ε 9.65 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)