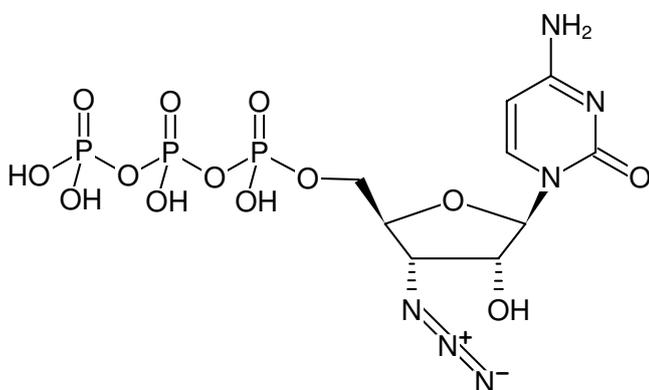


**3'-Azido-3'-dCTP**

3'-Azido-3'-deoxycytidine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-993	100 µl (10 mM)



Structural formula of 3'-Azido-3'-dCTP

**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>9</sub>H<sub>15</sub>N<sub>6</sub>O<sub>13</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 508.17 g/mol (free acid)**Exact Mass:** 507.99 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> 271 nm, ε 8.9 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)