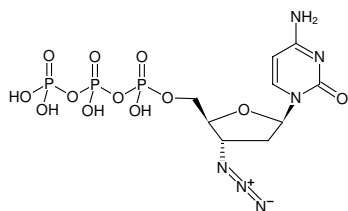


**3'-Azido-2',3'-ddCTP**

3'-Azido-2',3'-dideoxycytosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-246S	10 µl (100 mM)
NU-246L	5 x 10 µl (100 mM)



Structural formula of 3'-Azido-2',3'-ddCTP

**For research use only!****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>12</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 492.17 g/mol (free acid)**Exact Mass:** 492.00 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 100 mM - 110 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)