

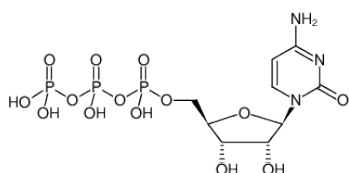
**CTP - Solid**

Cytidine 5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1050-100	100 mg
NU-1050-1G	1 g
NU-1050-10G	10 g
NU-1050-100G	100 g

**Quality Control Specifications:**

in vitro transcription (T7 RNA polymerase): visible RNA fragments after 5 min incubation, Proteases: not detectable



Structural formula of CTP - Solid

**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**Molecular Formula:** C<sub>9</sub>H<sub>16</sub>N<sub>3</sub>O<sub>14</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 483.16 g/mol (free acid)**CAS#:** 36051-68-0**EC number:** 252-849-3**Purity:** ≥ 95 % (HPLC)**Form:** Powder**Color:** White to slightly yellow**Solubility:** Soluble in water. Clear solution.**Spectroscopic Properties:** λ<sub>max</sub> 271 nm, ε 8.9 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.0)