



NTP Mix - 25 mM Solution

Premix of 25 mM ATP, CTP, GTP and UTP

Adenosine-5'-triphosphate, Sodium salt; Cytidine-5'-triphosphate, Sodium salt; Guanosine-5'-triphosphate, Sodium salt; Uridine-5'-triphosphate, Sodium salt

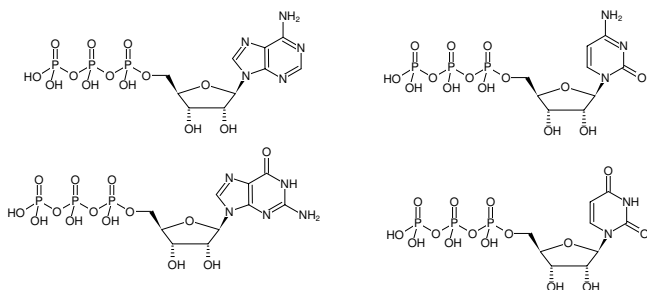
Cat. No.	Amount
NU-1024S	1 ml
NU-1024L	5x 1 ml

Description:

NTP Mix / 25 mM is an equimolar mixture of 25 mM ATP, CTP, GTP and UTP supplied as ultrapure aqueous solution and suitable for all molecular biology applications including *in vitro* transcription and RNA labeling.

Quality Control Specifications:

in vitro transcription: suitable
contamination with bacterial and human DNA: not detectable
activity of DNase, Protease or Phosphatase: not detectable



Structural formula of NTP Mix - 25 mM Solution

For *in vitro* 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. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

Shelf Life: 12 months

Molecular Formula:

ATP: C₁₀H₁₆N₅O₁₃P₃ (free acid)

CTP: C₉H₁₆N₃O₁₄P₃ (free acid)

GTP: C₁₀H₁₆N₅O₁₄P₃ (free acid)

UTP: C₉H₁₅N₂O₁₅P₃ (free acid)

Molecular Weight:

ATP: 507.18 g/mol (free acid)

CTP: 483.16 g/mol (free acid)

GTP: 523.18 g/mol (free acid)

UTP: 484.14 g/mol (free acid)

Purity: ≥ 99 % (HPLC)

Form: clear aqueous solution

Concentration: 25 mM each nucleotide

pH: 8.0 ± 0.2 (22 °C)