

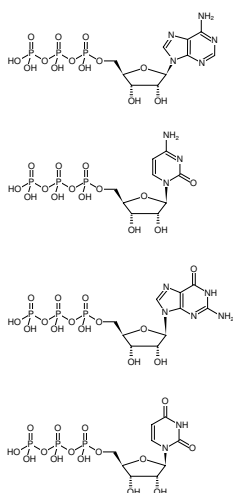


NTP Bundle

4 x 100 mM (ATP, CTP, GTP, 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-1014S	4 x 200 µl (4 x 20 µmol)
NU-1014L	4 x 1 ml (4 x 100 µmol)



Structural formula of NTP Bundle

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.

Shelf Life: 12 months after date of delivery

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: 100 mM ±2 %

pH: 8.0 ±0.2 (22 °C)

Spectroscopic Properties: ATP: λ_{max} 259 nm, ε 15.1 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.0)

CTP: λ_{max} 271 nm, ε 8.9 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.0)

GTP: λ_{max} 252 nm, ε 14.2 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.0)

UTP: λ_{max} 262 nm, ε 9.8 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.0)



NTP Bundle

4 x 100 mM (ATP, CTP, GTP, UTP)

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

Description:

NTP Bundle contains four separate solutions of ultrapure ATP, CTP, GTP and UTP supplied as clear aqueous solutions (pH 8.0).

ATP (#NU-1010)

CTP (#NU-1011)

GTP (#NU-1012)

UTP (#NU-1013)

Quality Control Specifications:

in vitro transcription: suitable

contamination with bacterial and human DNA: not detectable

activity of DNase, Protease or Phosphatase: not detectable