



## NTP Bundle

ATP, CTP, GTP and UTP 100 mM solutions

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

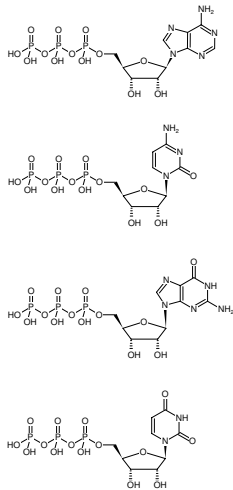
Cat. No.	Amount
NU-1014S	4 x 200 µl (4 x 20 µmol)
NU-1014L	4 x 1 ml (4 x 100 µmol)

(Tris-HCl pH 7.0)

CTP:  $\lambda_{\max}$  271 nm,  $\epsilon$  8.9 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.0)

GTP:  $\lambda_{\max}$  252 nm,  $\epsilon$  14.2 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.0)

UTP:  $\lambda_{\max}$  262 nm,  $\epsilon$  9.8 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.0)



Structural formula of NTP Bundle

**For general laboratory use.**

**Shipping:** shipped on gel packs

**Storage Conditions:** store at -20 °C

**Shelf Life:** 12 months after date of delivery

**Molecular Formula:**

ATP: C<sub>10</sub>H<sub>16</sub>N<sub>5</sub>O<sub>13</sub>P<sub>3</sub> (free acid)

CTP: C<sub>9</sub>H<sub>16</sub>N<sub>3</sub>O<sub>14</sub>P<sub>3</sub> (free acid)

GTP: C<sub>10</sub>H<sub>16</sub>N<sub>5</sub>O<sub>14</sub>P<sub>3</sub> (free acid)

UTP: C<sub>9</sub>H<sub>15</sub>N<sub>2</sub>O<sub>15</sub>P<sub>3</sub> (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:  $\lambda_{\max}$  259 nm,  $\epsilon$  15.1 L mmol<sup>-1</sup> cm<sup>-1</sup>



## NTP Bundle

ATP, CTP, GTP and UTP 100 mM solutions

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