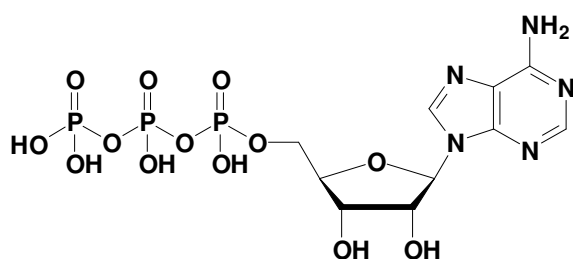


ATP, solution

Adenosine 5'-triphosphate, sodium salt

100 mM

Cat. No.	Amount
NU-1010	1 ml (100 μ mol)



For *in vitro* use only

Quality guaranteed for 12 months

Store at -20°C, short term (up to one week) exposure to ambient temperature possible

Concentration

100 mM +/-2%

Form

clear aqueous solution, pH 8.0 +/-0.2 (4°C)

Purity

>99%

Molecular Formula

C₁₀H₁₃N₅O₁₃P₃ (Anion)

Molecular Weight

504.16 (Anion)

Absorbance

absorbance max: 259 nm (pH 7)

ϵ at absorbance max: 15.1 mmol⁻¹ cm⁻¹

Quality Control Specifications

in vitro transcription:

suitable

contamination with bacterial and human DNA:

not detectable

activity of DNase, Protease or Phosphatase:

not detectable

Description

Ultrapure ATP supplied as clear aqueous solution (pH 8.0).

Applications:

ATP-sensitive calcium channels^[1]

V-ATPases (cellular proton pumps)^[2]

ATP-coupled chromatin remodelling^[3]

ATP-binding cassette transporters^[4]

ATP-grasp enzymes^[5]

Selected References:

[1] Wang *et al.* (2011) The biological effect of endogenous sulfur dioxide in the cardiovascular system. *Eur. J. Pharmacol.* **670(1)**:1.

[2] Scott *et al.* (2011) Duelling functions of the V-ATPase. *EMBO J.* **30(20)**:4113.

[3] Erdel *et al.* (2011) Chromatin remodelling in mammalian cells by ISWI-type complexes—where, when and why? *FEBS J.* **278(19)**:3608.

[4] Gatti *et al.* (2011) Novel insights into targeting ATP-binding cassette transporters for antitumor therapy. *Curr. Med. Chem.* **18(27)**:4237.

[5] Fawaz *et al.* (2011) The ATP-grasp enzymes. *Bioorg. Chem.* **39(5-6)**:185.

Erlich *et al.* (1988) Primer-directed enzymatic amplification of DNA with a thermostable DNA polymerase. *Science* **29 (239)**:487.

Gelfand *et al.* (1991) Detection of specific polymerase chain reaction product by utilizing the 5'-3' exonuclease activity of *Thermus aquaticus* DNA polymerase. *Proc. Natl. Acad. Sci. USA* **88 (16)**:7276.

Sanger *et al.* (1977) DNA sequencing with chain-terminating inhibitors. *Proc. Natl. Acad. Sci. USA* **74**:5463.