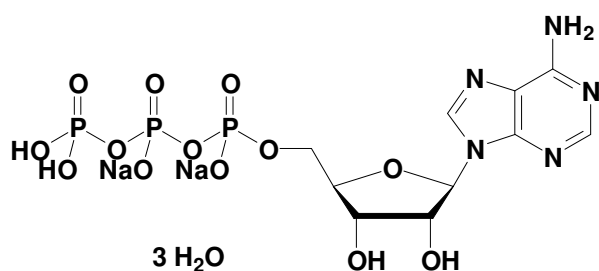


ATP, lyophilized Adenosine 5'-triphosphate, sodium salt

Cat. No.	Amount
NU-1010-10G	10 g, lyophilized
NU-1010-100G	100 g, lyophilized



Molecular Formula: C₁₀H₁₄N₅Na₂O₁₃P₃ × 3H₂O

Molecular Weight: 605.18

Absorbance

absorbance max: 259 nm (pH 7)

ε at absorbance max: 15.1 mmol⁻¹ cm⁻¹

Purity: >95%

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C.

For research use only!

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.