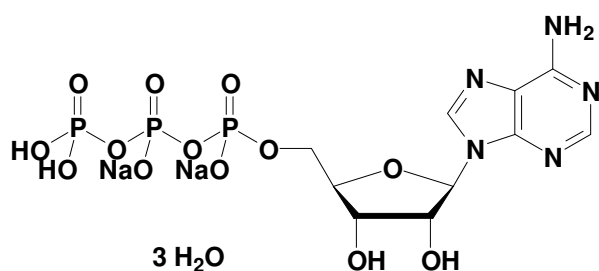


ATP, lyophilized

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

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!