

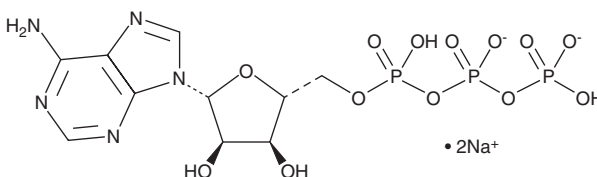
PRODUCT INFORMATION



Adenosine 5'-triphosphate (sodium salt) (solution)

Item No. 40182

CAS Registry No.: 987-65-5
Formal Name: adenosine 5'-(tetrahydrogen triphosphate), disodium salt
Synonyms: 5'-ATP, ATP, NSC 20268
MF: C₁₀H₁₄N₅O₁₃P₃ • 2Na
FW: 551.1
Purity: ≥95%
Supplied as: An aqueous solution titrated with sodium hydroxide to a pH 7.3-7.5



Storage: -20°C
Stability: ≥1 year

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Adenosine 5'-triphosphate (ATP) (sodium salt) (solution) is supplied as an aqueous solution titrated with sodium hydroxide to a pH 7.3-7.5. ATP (sodium salt) (solution) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the aqueous solution of ATP (sodium salt) (solution) should be diluted with the aqueous buffer of choice. The solubility of ATP (sodium salt) (solution) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

ATP is a central component of energy storage and metabolism *in vivo*, providing the metabolic energy to drive metabolic pumps and serving as a coenzyme in a wide array of enzymatic reactions.¹ ATP is a substrate for kinases involved in cell signaling and of adenylate cyclases that produce the second messenger cAMP.¹ It is utilized in various cellular processes including, respiration, biosynthetic reactions, motility, and cell division.

Reference

1. Knowles, J.R. Enzyme-catalyzed phosphoryl transfer reactions. *Annu. Rev. Biochem.* **49**, 877-919 (1980).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/13/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM