

PRODUCT INFORMATION



2-Aminoadenosine-5'-O-triphosphate (sodium salt)

Item No. 38424

Formal Name: 2-amino-adenosine 5'-(tetrahydrogen triphosphate), tetrasodium salt

Synonyms: 2-amino ATP, 2,6-Diaminopurine Riboside-5'-O-triphosphate

MF: $C_{10}H_{13}N_6O_{13}P_3 \cdot 4Na$

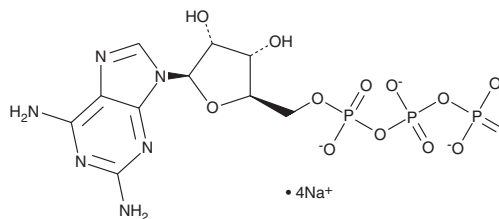
FW: 610.1

Purity: $\geq 95\%$

Supplied as: A 10 mM solution in water

Storage: $-80^{\circ}C$

Stability: ≥ 4 years



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

Description

2-Aminoadenosine-5'-O-triphosphate (2-amino ATP) is a modified nucleotide.¹ It promotes microtubule assembly *in vitro*.² 2-amino ATP has been used as an alternative substrate for T4 DNA ligase in ligase-catalyzed oligonucleotide polymerizations with 96.6% fidelity.³ Incorporation of 2-amino ATP into complementary RNA (cRNA) increases signal detection in microarrays.¹

References

1. Nguyen, A., Zhao, C., Dorris, D., *et al.* Quantitative assessment of the use of modified nucleoside triphosphates in expression profiling: Differential effects on signal intensities and impacts on expression ratios. *BMC Biotechnol.* **2**, 14 (2002).
2. Muraoka, M. and Sakai, H. Effects of purinenucleotide analogues on microtubule assembly. *Cell Struct. Funct.* **24(5)**, 305-312 (1999).
3. Lei, Y. and Hili, R. Structure-activity relationships of the ATP cofactor in ligase-catalysed oligonucleotide polymerisations. *Org. Biomol. Chem.* **15(11)**, 2349-2352 (2017).

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, 04/14/2026

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