

PRODUCT INFORMATION



3'-Amino-3'-deoxyadenosine-5'-O-diphosphate (sodium salt)

Item No. 38363

Formal Name: 3'-amino-3'-deoxy-adenosine
5'-(trihydrogen diphosphate),
trisodium salt

MF: $C_{10}H_{13}N_6O_9P_2 \cdot 3Na$

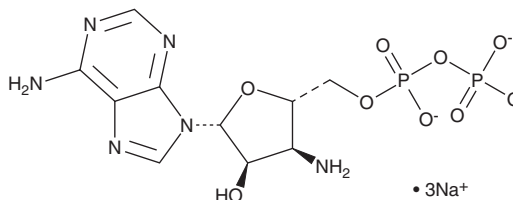
FW: 492.2

Purity: $\geq 98\%$

Supplied as: A 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

3'-Amino-3'-deoxyadenosine-5'-O-diphosphate is an amine-modified nucleotide that has been used in the synthesis of fluorescently labeled ADP.¹ It also has been used in the synthesis of phosphorylated probes for GTP pyrophosphokinase domains.²

References

1. Webb, M.R., Reid, G.P., Munasinghe, V.R.N., *et al.* A series of related nucleotide analogues that aids optimization of fluorescence signals in probing the mechanism of P-loop ATPases, such as actomyosin. *Biochemistry* **43**(45), 14463-14471 (2004).
2. Mojir, V., Roghanian, M., Tamman, H., *et al.* Nonhydrolysable analogues of (p)ppGpp and (p)ppApp alarmone nucleotides as novel molecular tools. *ACS Chem. Biol.* **16**(9), 1680-1691 (2021).

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, 04/13/2026

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