

# PRODUCT INFORMATION



## 3'-Amino-3'-deoxyadenosine-5'-O-triphosphate (sodium salt)

Item No. 38364

**Formal Name:** 3'-amino-3'-deoxy-adenosine 5'-(tetrahydrogen triphosphate), tetrasodium salt

**Synonyms:** 3'-Deoxy-3'-amino ATP, 3'-NH<sub>2</sub>-dATP

**MF:** C<sub>10</sub>H<sub>13</sub>N<sub>6</sub>O<sub>12</sub>P<sub>3</sub> • 4Na

**FW:** 594.1

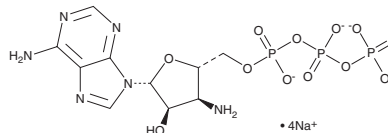
**Purity:** ≥98%

**Supplied as:** A solution in water

**Storage:** -80°C

**Stability:** ≥2 years

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



### Description

3'-Amino-3'-deoxyadenosine-5'-O-triphosphate is an amine-modified nucleotide that has been used as a probe for quantifying ATP substrate specificity of kinases and ligases.<sup>1</sup> It terminates RNA strand synthesis catalyzed by RNA polymerase in *E. coli*.<sup>2</sup>

### References

1. Kinoshita, Y. and Nishigaki, K. Unexpectedly general replaceability of ATP in ATP-requiring enzymes. *J. Biochem.* **122**(1), 205-211 (1997).
2. Kutateladze, T., Beabealashvili, R., Azhayev, A., *et al.* 3'-deoxy-3'-aminonucleoside 5'-triphosphates — terminators of RNA synthesis, catalyzed by DNA-dependent RNA polymerase from *Escherichia coli*. *FEBS Lett.* **153**(2), 420-426 (1983).

#### 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, 03/23/2023

#### 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